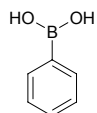




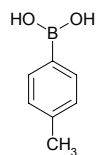
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	P20009	121.93



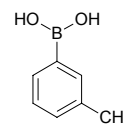
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	393622	135.96



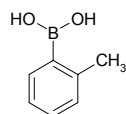
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	393614	135.96



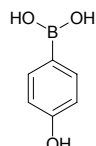
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	393606	135.96



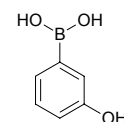
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	523976	137.93



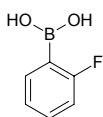
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	523968	137.93



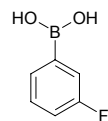
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	445223	139.92



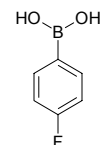
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	441643	139.92



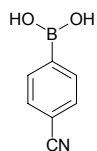
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	417556	139.92



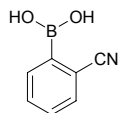
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	521418	146.94



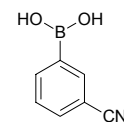
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	521396	146.94



**Sigma - Aldrich
SAR Boronic Acids**

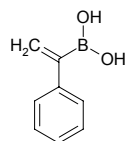
Brand	Catalog Number	MW
ALDRICH	513016	146.94





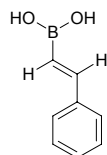
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	571350	147.97



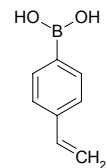
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	473790	147.97



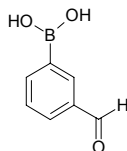
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	417580	147.97



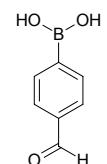
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	441651	149.94



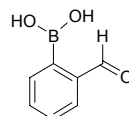
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	431966	149.94



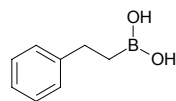
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	431958	149.94



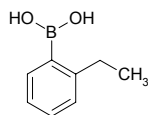
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	588423	149.99



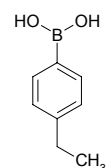
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	521523	149.99



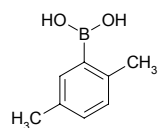
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	499536	149.99



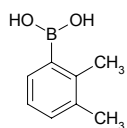
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	483516	149.99



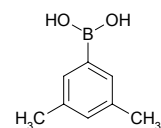
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	483508	149.99



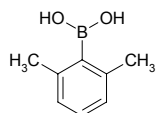
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	480088	149.99

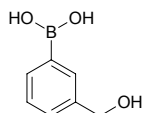


**Sigma - Aldrich
SAR Boronic Acids**

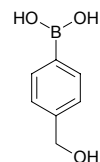
Brand	Catalog Number	MW
ALDRICH	480061	149.99

**Sigma - Aldrich
SAR Boronic Acids**

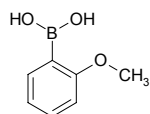
Brand	Catalog Number	MW
ALDRICH	512834	151.96

**Sigma - Aldrich
SAR Boronic Acids**

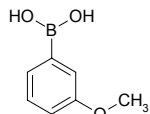
Brand	Catalog Number	MW
ALDRICH	512338	151.96

**Sigma - Aldrich
SAR Boronic Acids**

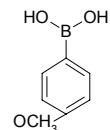
Brand	Catalog Number	MW
ALDRICH	445231	151.96

**Sigma - Aldrich
SAR Boronic Acids**

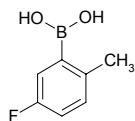
Brand	Catalog Number	MW
ALDRICH	441686	151.96

**Sigma - Aldrich
SAR Boronic Acids**

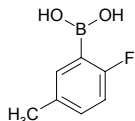
Brand	Catalog Number	MW
ALDRICH	417599	151.96

**Sigma - Aldrich
SAR Boronic Acids**

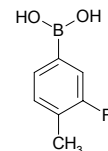
Brand	Catalog Number	MW
ALDRICH	567442	153.95

**Sigma - Aldrich
SAR Boronic Acids**

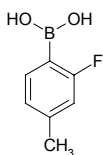
Brand	Catalog Number	MW
ALDRICH	567434	153.95

**Sigma - Aldrich
SAR Boronic Acids**

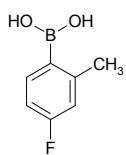
Brand	Catalog Number	MW
ALDRICH	567426	153.95

**Sigma - Aldrich
SAR Boronic Acids**

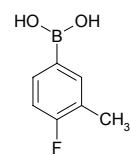
Brand	Catalog Number	MW
ALDRICH	567418	153.95

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	565652	153.95

**Sigma - Aldrich
SAR Boronic Acids**

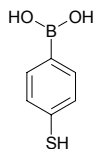
Brand	Catalog Number	MW
ALDRICH	483567	153.95





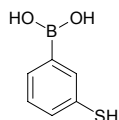
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	524018	154.00



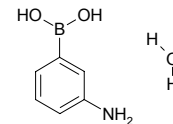
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	523992	154.00



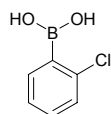
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	287512	154.96



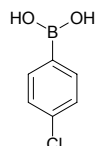
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	445215	156.38



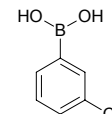
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	417548	156.38



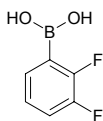
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	417521	156.38



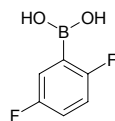
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	514039	157.91



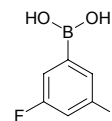
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	514020	157.91



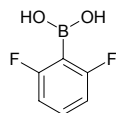
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	471925	157.91



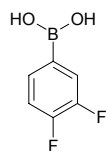
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	470791	157.91



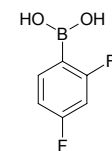
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	465089	157.91



**Sigma - Aldrich
SAR Boronic Acids**

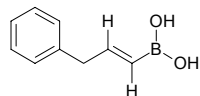
Brand	Catalog Number	MW
ALDRICH	465070	157.91





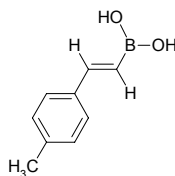
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	642606	162.00



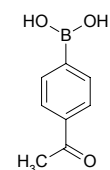
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	568139	162.00



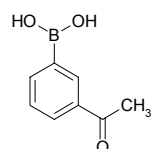
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	470821	163.97



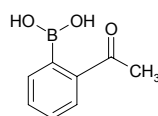
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	470813	163.97



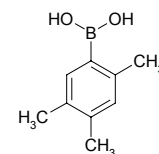
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	470805	163.97



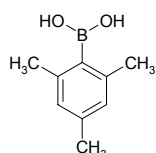
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	542326	164.01



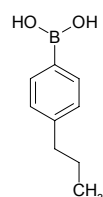
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	542318	164.01



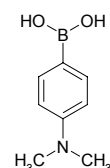
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	521507	164.01



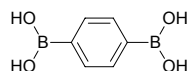
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	483532	165.00



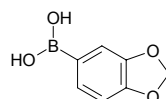
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	417130	165.75



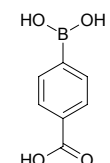
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	499994	165.94



**Sigma - Aldrich
SAR Boronic Acids**

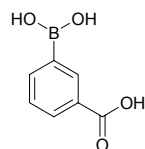
Brand	Catalog Number	MW
ALDRICH	456772	165.94





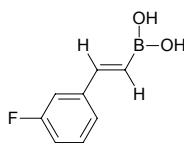
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	456764	165.94



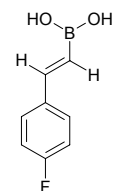
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	637971	165.96



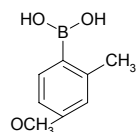
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	518972	165.96



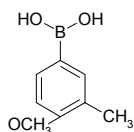
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	639370	165.99



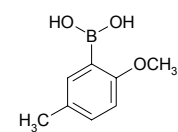
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	595500	165.99



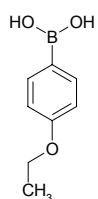
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	567507	165.99



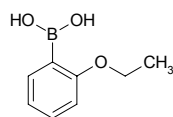
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	455539	165.99



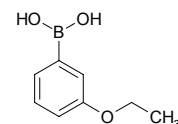
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	455520	165.99



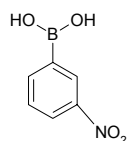
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	441635	165.99



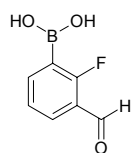
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	325104	166.93



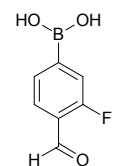
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	645788	167.93



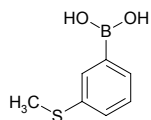
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	542296	167.93

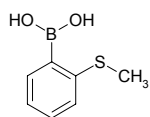


**Sigma - Aldrich
SAR Boronic Acids**

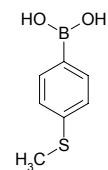
Brand	Catalog Number	MW
ALDRICH	526002	168.02

**Sigma - Aldrich
SAR Boronic Acids**

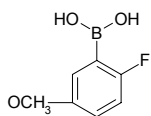
Brand	Catalog Number	MW
ALDRICH	521248	168.02

**Sigma - Aldrich
SAR Boronic Acids**

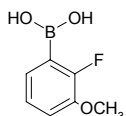
Brand	Catalog Number	MW
ALDRICH	456802	168.02

**Sigma - Aldrich
SAR Boronic Acids**

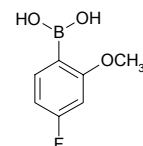
Brand	Catalog Number	MW
ALDRICH	597112	169.95

**Sigma - Aldrich
SAR Boronic Acids**

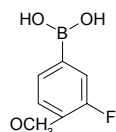
Brand	Catalog Number	MW
ALDRICH	594253	169.95

**Sigma - Aldrich
SAR Boronic Acids**

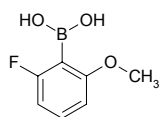
Brand	Catalog Number	MW
ALDRICH	564494	169.95

**Sigma - Aldrich
SAR Boronic Acids**

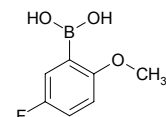
Brand	Catalog Number	MW
ALDRICH	564036	169.95

**Sigma - Aldrich
SAR Boronic Acids**

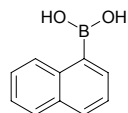
Brand	Catalog Number	MW
ALDRICH	512761	169.95

**Sigma - Aldrich
SAR Boronic Acids**

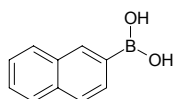
Brand	Catalog Number	MW
ALDRICH	483540	169.95

**Sigma - Aldrich
SAR Boronic Acids**

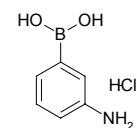
Brand	Catalog Number	MW
ALDRICH	N257	171.99

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	480134	171.99

**Sigma - Aldrich
SAR Boronic Acids**

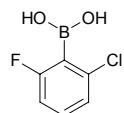
Brand	Catalog Number	MW
ALDRICH	410705	173.41





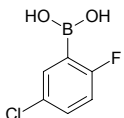
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	566071	174.37



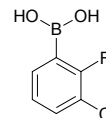
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	557234	174.37



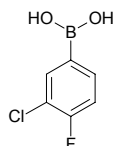
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	557226	174.37



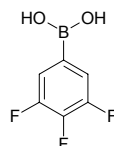
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	512230	174.37



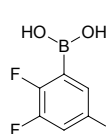
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	524700	175.90



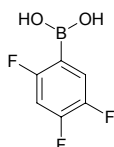
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	524697	175.90



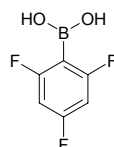
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	524689	175.90



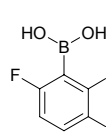
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	524107	175.90



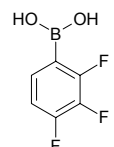
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	524093	175.90



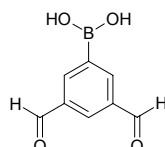
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	524085	175.90



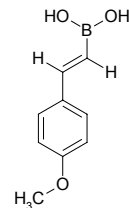
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	567485	177.95



**Sigma - Aldrich
SAR Boronic Acids**

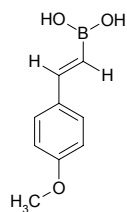
Brand	Catalog Number	MW
ALDRICH	518980	178.00





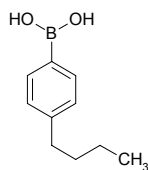
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	518980	178.00



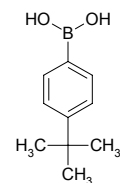
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	521493	178.04



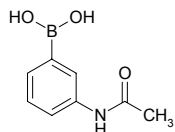
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	480053	178.04



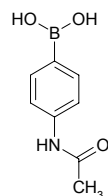
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	566012	178.98



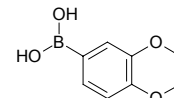
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	565806	178.98



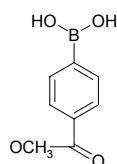
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	635995	179.97



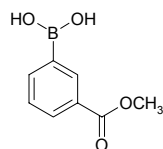
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	594539	179.97



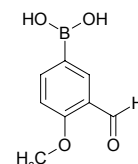
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	591130	179.97



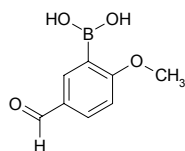
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	512869	179.97



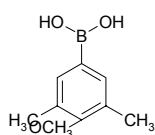
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	512257	179.97



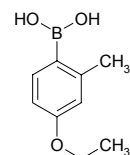
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	598062	180.01



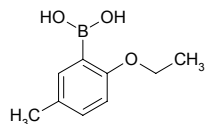
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	594830	180.01

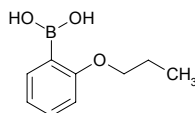


**Sigma - Aldrich
SAR Boronic Acids**

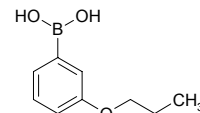
Brand	Catalog Number	MW
ALDRICH	567493	180.01

**Sigma - Aldrich
SAR Boronic Acids**

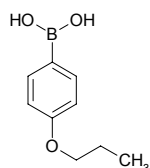
Brand	Catalog Number	MW
ALDRICH	557722	180.01

**Sigma - Aldrich
SAR Boronic Acids**

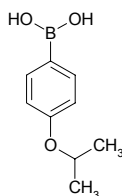
Brand	Catalog Number	MW
ALDRICH	557714	180.01

**Sigma - Aldrich
SAR Boronic Acids**

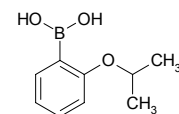
Brand	Catalog Number	MW
ALDRICH	557706	180.01

**Sigma - Aldrich
SAR Boronic Acids**

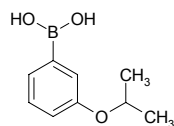
Brand	Catalog Number	MW
ALDRICH	542474	180.01

**Sigma - Aldrich
SAR Boronic Acids**

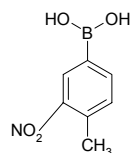
Brand	Catalog Number	MW
ALDRICH	542466	180.01

**Sigma - Aldrich
SAR Boronic Acids**

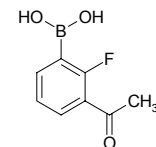
Brand	Catalog Number	MW
ALDRICH	542458	180.01

**Sigma - Aldrich
SAR Boronic Acids**

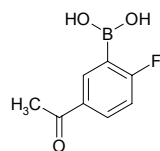
Brand	Catalog Number	MW
ALDRICH	521477	180.96

**Sigma - Aldrich
SAR Boronic Acids**

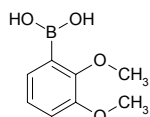
Brand	Catalog Number	MW
ALDRICH	651249	181.96

**Sigma - Aldrich
SAR Boronic Acids**

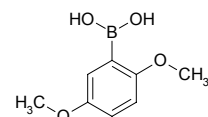
Brand	Catalog Number	MW
ALDRICH	639583	181.96

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	557730	181.99

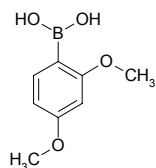
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	483494	181.99

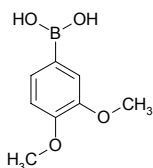


**Sigma - Aldrich
SAR Boronic Acids**

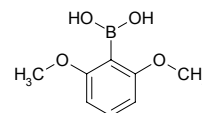
Brand	Catalog Number	MW
ALDRICH	483486	181.99

**Sigma - Aldrich
SAR Boronic Acids**

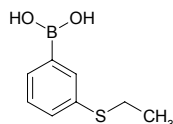
Brand	Catalog Number	MW
ALDRICH	480118	181.99

**Sigma - Aldrich
SAR Boronic Acids**

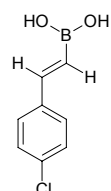
Brand	Catalog Number	MW
ALDRICH	480096	181.99

**Sigma - Aldrich
SAR Boronic Acids**

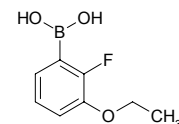
Brand	Catalog Number	MW
ALDRICH	597295	182.05

**Sigma - Aldrich
SAR Boronic Acids**

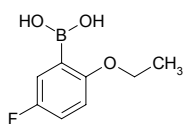
Brand	Catalog Number	MW
ALDRICH	523569	182.42

**Sigma - Aldrich
SAR Boronic Acids**

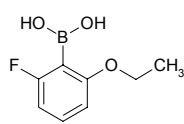
Brand	Catalog Number	MW
ALDRICH	639389	183.98

**Sigma - Aldrich
SAR Boronic Acids**

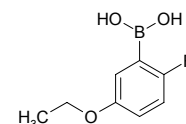
Brand	Catalog Number	MW
ALDRICH	594954	183.98

**Sigma - Aldrich
SAR Boronic Acids**

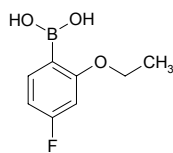
Brand	Catalog Number	MW
ALDRICH	594504	183.98

**Sigma - Aldrich
SAR Boronic Acids**

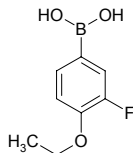
Brand	Catalog Number	MW
ALDRICH	594393	183.98

**Sigma - Aldrich
SAR Boronic Acids**

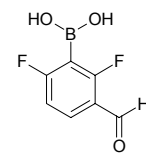
Brand	Catalog Number	MW
ALDRICH	564508	183.98

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	564044	183.98

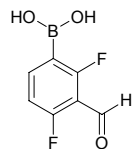
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	652210	185.92

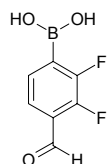


**Sigma - Aldrich
SAR Boronic Acids**

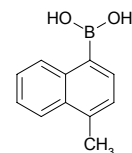
Brand	Catalog Number	MW
ALDRICH	597406	185.92

**Sigma - Aldrich
SAR Boronic Acids**

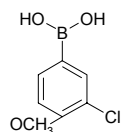
Brand	Catalog Number	MW
ALDRICH	571377	185.92

**Sigma - Aldrich
SAR Boronic Acids**

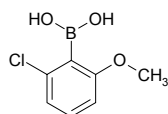
Brand	Catalog Number	MW
ALDRICH	521450	186.02

**Sigma - Aldrich
SAR Boronic Acids**

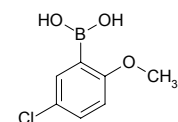
Brand	Catalog Number	MW
ALDRICH	564486	186.40

**Sigma - Aldrich
SAR Boronic Acids**

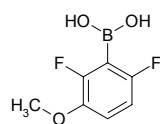
Brand	Catalog Number	MW
ALDRICH	512753	186.40

**Sigma - Aldrich
SAR Boronic Acids**

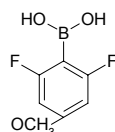
Brand	Catalog Number	MW
ALDRICH	512249	186.40

**Sigma - Aldrich
SAR Boronic Acids**

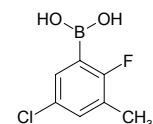
Brand	Catalog Number	MW
ALDRICH	652164	187.94

**Sigma - Aldrich
SAR Boronic Acids**

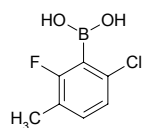
Brand	Catalog Number	MW
ALDRICH	593060	187.94

**Sigma - Aldrich
SAR Boronic Acids**

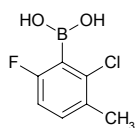
Brand	Catalog Number	MW
ALDRICH	557277	188.40

**Sigma - Aldrich
SAR Boronic Acids**

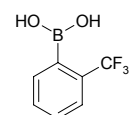
Brand	Catalog Number	MW
ALDRICH	557269	188.40

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	557250	188.40

**Sigma - Aldrich
SAR Boronic Acids**

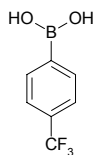
Brand	Catalog Number	MW
ALDRICH	445193	189.93





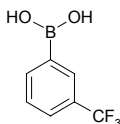
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	439320	189.93



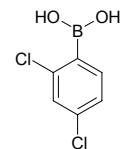
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	432032	189.93



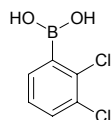
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	521388	190.82



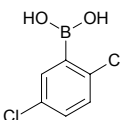
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	514047	190.82



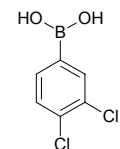
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	512311	190.82



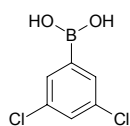
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	471917	190.82



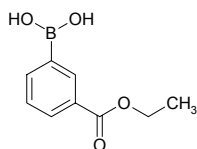
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	445207	190.82



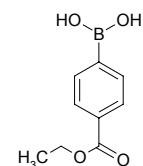
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	574651	194.00



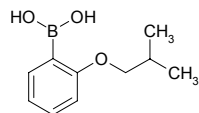
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	574643	194.00



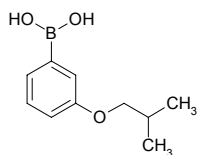
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	645745	194.04



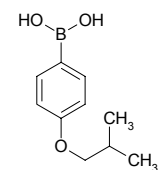
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	645737	194.04



**Sigma - Aldrich
SAR Boronic Acids**

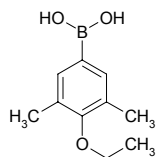
Brand	Catalog Number	MW
ALDRICH	639605	194.04





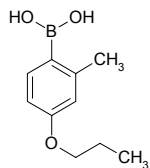
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	597945	194.04



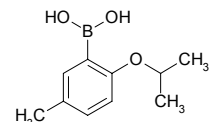
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	594725	194.04



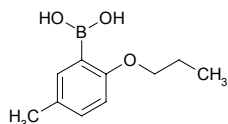
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	565660	194.04



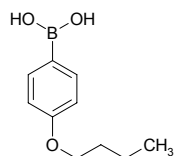
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	565644	194.04



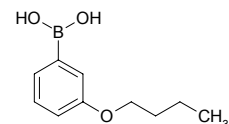
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	542504	194.04



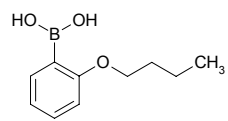
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	542490	194.04



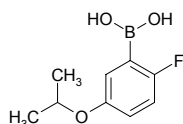
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	542482	194.04



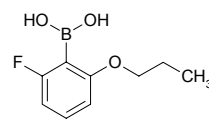
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	645206	198.00



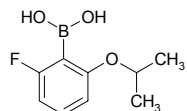
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	639419	198.00



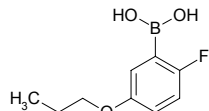
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	639397	198.00



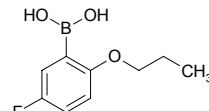
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	594288	198.00



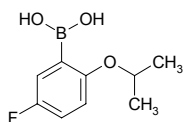
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	565776	198.00

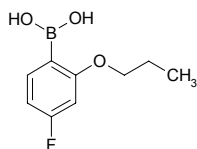


**Sigma - Aldrich
SAR Boronic Acids**

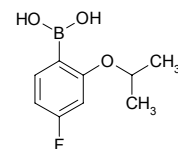
Brand	Catalog Number	MW
ALDRICH	564559	198.00

**Sigma - Aldrich
SAR Boronic Acids**

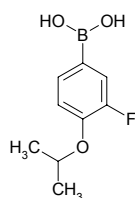
Brand	Catalog Number	MW
ALDRICH	564524	198.00

**Sigma - Aldrich
SAR Boronic Acids**

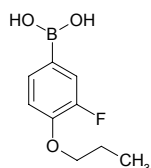
Brand	Catalog Number	MW
ALDRICH	564516	198.00

**Sigma - Aldrich
SAR Boronic Acids**

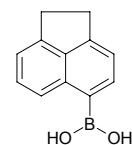
Brand	Catalog Number	MW
ALDRICH	564060	198.00

**Sigma - Aldrich
SAR Boronic Acids**

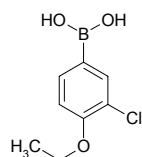
Brand	Catalog Number	MW
ALDRICH	564052	198.00

**Sigma - Aldrich
SAR Boronic Acids**

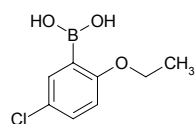
Brand	Catalog Number	MW
ALDRICH	567477	198.03

**Sigma - Aldrich
SAR Boronic Acids**

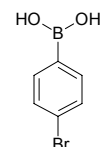
Brand	Catalog Number	MW
ALDRICH	564478	200.43

**Sigma - Aldrich
SAR Boronic Acids**

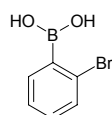
Brand	Catalog Number	MW
ALDRICH	542547	200.43

**Sigma - Aldrich
SAR Boronic Acids**

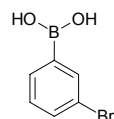
Brand	Catalog Number	MW
ALDRICH	B75956	200.83

**Sigma - Aldrich
SAR Boronic Acids**

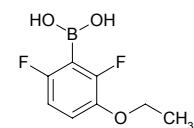
Brand	Catalog Number	MW
ALDRICH	473804	200.83

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	441627	200.83

**Sigma - Aldrich
SAR Boronic Acids**

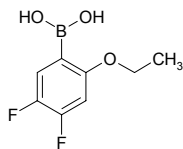
Brand	Catalog Number	MW
ALDRICH	650951	201.97





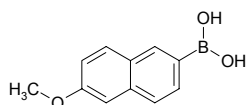
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	650943	201.97



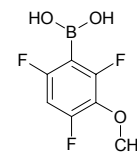
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	521892	202.02



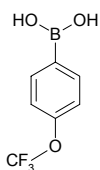
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	652202	205.93



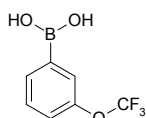
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	510130	205.93



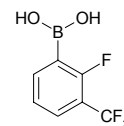
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	510122	205.93



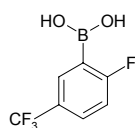
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	558206	207.92



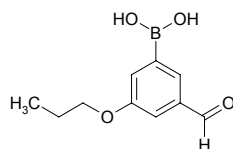
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	558184	207.92



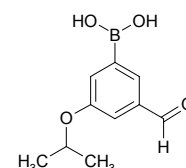
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	657638	208.02



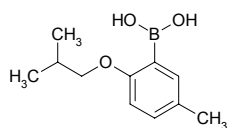
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	657557	208.02



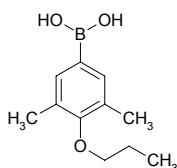
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	651230	208.07



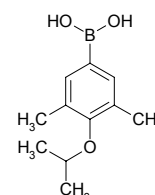
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	597848	208.07



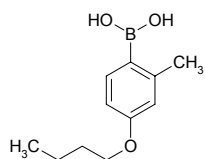
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	597732	208.07

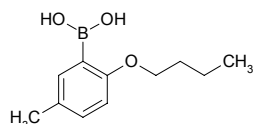


**Sigma - Aldrich
SAR Boronic Acids**

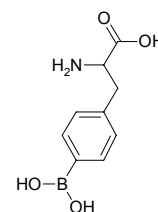
Brand	Catalog Number	MW
ALDRICH	594628	208.07

**Sigma - Aldrich
SAR Boronic Acids**

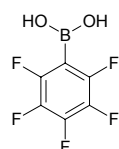
Brand	Catalog Number	MW
ALDRICH	565679	208.07

**Sigma - Aldrich
SAR Boronic Acids**

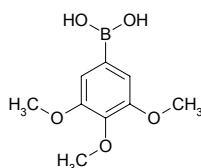
Brand	Catalog Number	MW
ALDRICH	512680	209.01

**Sigma - Aldrich
SAR Boronic Acids**

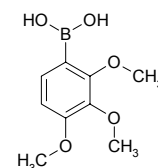
Brand	Catalog Number	MW
ALDRICH	465097	211.88

**Sigma - Aldrich
SAR Boronic Acids**

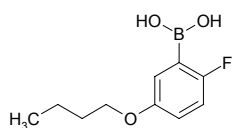
Brand	Catalog Number	MW
ALDRICH	651648	212.01

**Sigma - Aldrich
SAR Boronic Acids**

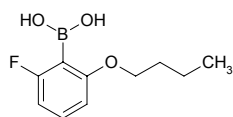
Brand	Catalog Number	MW
ALDRICH	512281	212.01

**Sigma - Aldrich
SAR Boronic Acids**

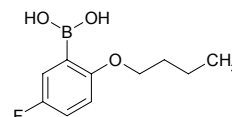
Brand	Catalog Number	MW
ALDRICH	645230	212.03

**Sigma - Aldrich
SAR Boronic Acids**

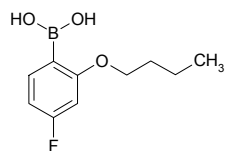
Brand	Catalog Number	MW
ALDRICH	639427	212.03

**Sigma - Aldrich
SAR Boronic Acids**

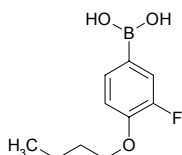
Brand	Catalog Number	MW
ALDRICH	564540	212.03

**Sigma - Aldrich
SAR Boronic Acids**

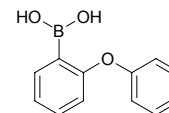
Brand	Catalog Number	MW
ALDRICH	564532	212.03

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	564079	212.03

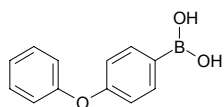
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	521485	214.03

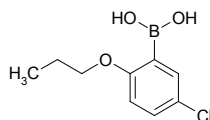


**Sigma - Aldrich
SAR Boronic Acids**

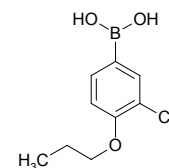
Brand	Catalog Number	MW
ALDRICH	480142	214.03

**Sigma - Aldrich
SAR Boronic Acids**

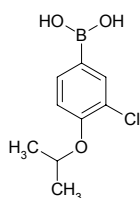
Brand	Catalog Number	MW
ALDRICH	645192	214.46

**Sigma - Aldrich
SAR Boronic Acids**

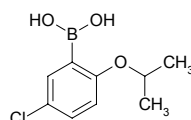
Brand	Catalog Number	MW
ALDRICH	564451	214.46

**Sigma - Aldrich
SAR Boronic Acids**

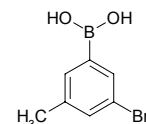
Brand	Catalog Number	MW
ALDRICH	564443	214.46

**Sigma - Aldrich
SAR Boronic Acids**

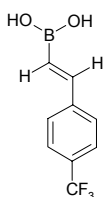
Brand	Catalog Number	MW
ALDRICH	542555	214.46

**Sigma - Aldrich
SAR Boronic Acids**

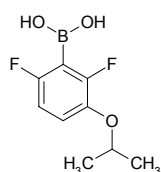
Brand	Catalog Number	MW
ALDRICH	645311	214.86

**Sigma - Aldrich
SAR Boronic Acids**

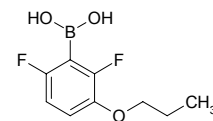
Brand	Catalog Number	MW
ALDRICH	519022	215.97

**Sigma - Aldrich
SAR Boronic Acids**

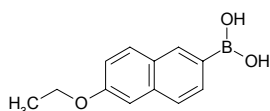
Brand	Catalog Number	MW
ALDRICH	652180	215.99

**Sigma - Aldrich
SAR Boronic Acids**

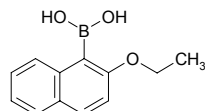
Brand	Catalog Number	MW
ALDRICH	652172	215.99

**Sigma - Aldrich
SAR Boronic Acids**

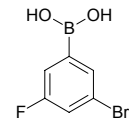
Brand	Catalog Number	MW
ALDRICH	521906	216.05

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	521884	216.05

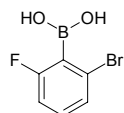
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	645346	218.82

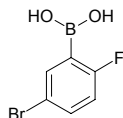


**Sigma - Aldrich
SAR Boronic Acids**

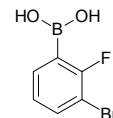
Brand	Catalog Number	MW
ALDRICH	593737	218.82

**Sigma - Aldrich
SAR Boronic Acids**

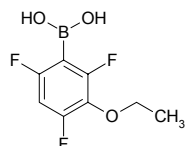
Brand	Catalog Number	MW
ALDRICH	593621	218.82

**Sigma - Aldrich
SAR Boronic Acids**

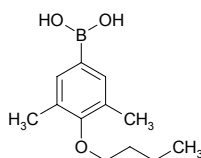
Brand	Catalog Number	MW
ALDRICH	558214	218.82

**Sigma - Aldrich
SAR Boronic Acids**

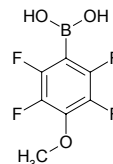
Brand	Catalog Number	MW
ALDRICH	657247	219.96

**Sigma - Aldrich
SAR Boronic Acids**

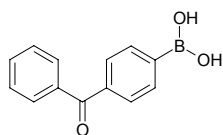
Brand	Catalog Number	MW
ALDRICH	597511	222.09

**Sigma - Aldrich
SAR Boronic Acids**

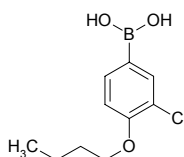
Brand	Catalog Number	MW
ALDRICH	657301	223.92

**Sigma - Aldrich
SAR Boronic Acids**

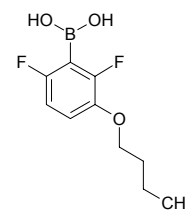
Brand	Catalog Number	MW
ALDRICH	526010	226.04

**Sigma - Aldrich
SAR Boronic Acids**

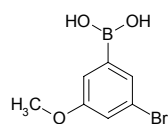
Brand	Catalog Number	MW
ALDRICH	564435	228.49

**Sigma - Aldrich
SAR Boronic Acids**

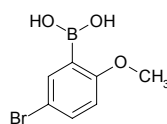
Brand	Catalog Number	MW
ALDRICH	652199	230.02

**Sigma - Aldrich
SAR Boronic Acids**

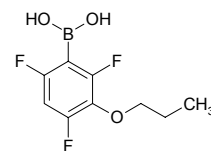
Brand	Catalog Number	MW
ALDRICH	652326	230.86

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	512702	230.86

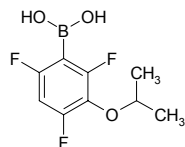
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	657379	233.98

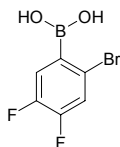


**Sigma - Aldrich
SAR Boronic Acids**

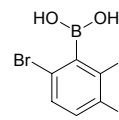
Brand	Catalog Number	MW
ALDRICH	657360	233.98

**Sigma - Aldrich
SAR Boronic Acids**

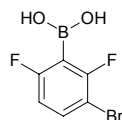
Brand	Catalog Number	MW
ALDRICH	645281	236.81

**Sigma - Aldrich
SAR Boronic Acids**

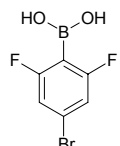
Brand	Catalog Number	MW
ALDRICH	635774	236.81

**Sigma - Aldrich
SAR Boronic Acids**

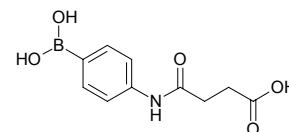
Brand	Catalog Number	MW
ALDRICH	557242	236.81

**Sigma - Aldrich
SAR Boronic Acids**

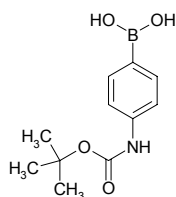
Brand	Catalog Number	MW
ALDRICH	557218	236.81

**Sigma - Aldrich
SAR Boronic Acids**

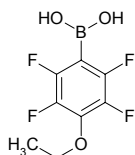
Brand	Catalog Number	MW
ALDRICH	578754	237.02

**Sigma - Aldrich
SAR Boronic Acids**

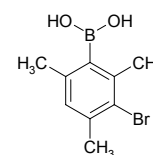
Brand	Catalog Number	MW
ALDRICH	565814	237.07

**Sigma - Aldrich
SAR Boronic Acids**

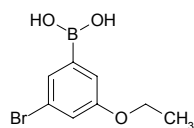
Brand	Catalog Number	MW
ALDRICH	657298	237.95

**Sigma - Aldrich
SAR Boronic Acids**

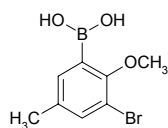
Brand	Catalog Number	MW
ALDRICH	593850	242.91

**Sigma - Aldrich
SAR Boronic Acids**

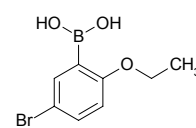
Brand	Catalog Number	MW
ALDRICH	651192	244.88

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	596183	244.88

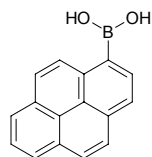
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	512826	244.88

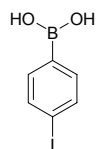


**Sigma - Aldrich
SAR Boronic Acids**

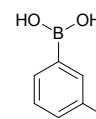
Brand	Catalog Number	MW
ALDRICH	542873	246.08

**Sigma - Aldrich
SAR Boronic Acids**

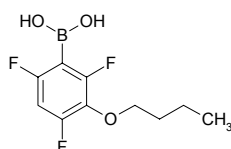
Brand	Catalog Number	MW
ALDRICH	471933	247.83

**Sigma - Aldrich
SAR Boronic Acids**

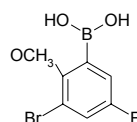
Brand	Catalog Number	MW
ALDRICH	441678	247.83

**Sigma - Aldrich
SAR Boronic Acids**

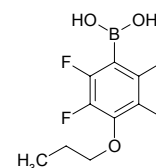
Brand	Catalog Number	MW
ALDRICH	657352	248.01

**Sigma - Aldrich
SAR Boronic Acids**

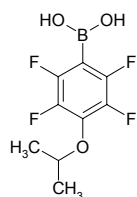
Brand	Catalog Number	MW
ALDRICH	645176	248.85

**Sigma - Aldrich
SAR Boronic Acids**

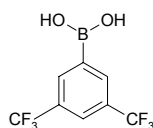
Brand	Catalog Number	MW
ALDRICH	657271	251.97

**Sigma - Aldrich
SAR Boronic Acids**

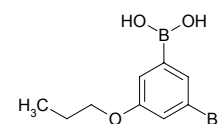
Brand	Catalog Number	MW
ALDRICH	657263	251.97

**Sigma - Aldrich
SAR Boronic Acids**

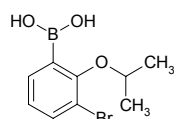
Brand	Catalog Number	MW
ALDRICH	471070	257.93

**Sigma - Aldrich
SAR Boronic Acids**

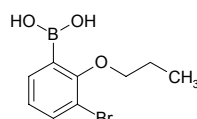
Brand	Catalog Number	MW
ALDRICH	657514	258.91

**Sigma - Aldrich
SAR Boronic Acids**

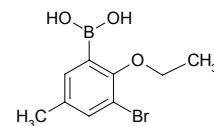
Brand	Catalog Number	MW
ALDRICH	597171	258.91

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	597074	258.91

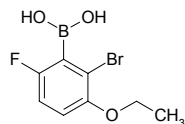
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	596299	258.91

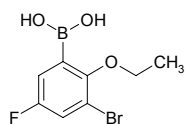


**Sigma - Aldrich
SAR Boronic Acids**

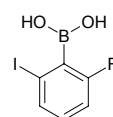
Brand	Catalog Number	MW
ALDRICH	645710	262.87

**Sigma - Aldrich
SAR Boronic Acids**

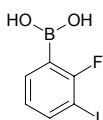
Brand	Catalog Number	MW
ALDRICH	542512	262.87

**Sigma - Aldrich
SAR Boronic Acids**

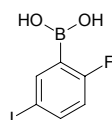
Brand	Catalog Number	MW
ALDRICH	639451	265.82

**Sigma - Aldrich
SAR Boronic Acids**

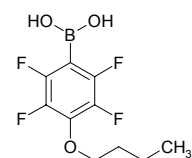
Brand	Catalog Number	MW
ALDRICH	593524	265.82

**Sigma - Aldrich
SAR Boronic Acids**

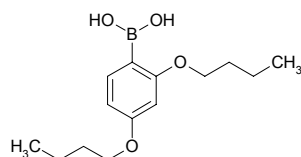
Brand	Catalog Number	MW
ALDRICH	593419	265.82

**Sigma - Aldrich
SAR Boronic Acids**

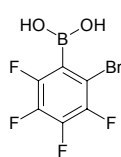
Brand	Catalog Number	MW
ALDRICH	657255	266.00

**Sigma - Aldrich
SAR Boronic Acids**

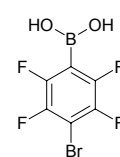
Brand	Catalog Number	MW
ALDRICH	650994	266.15

**Sigma - Aldrich
SAR Boronic Acids**

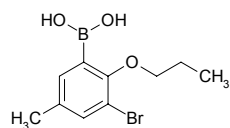
Brand	Catalog Number	MW
ALDRICH	645303	272.79

**Sigma - Aldrich
SAR Boronic Acids**

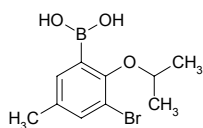
Brand	Catalog Number	MW
ALDRICH	593966	272.79

**Sigma - Aldrich
SAR Boronic Acids**

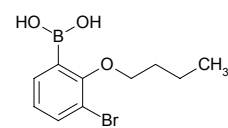
Brand	Catalog Number	MW
ALDRICH	596744	272.94

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	596620	272.94

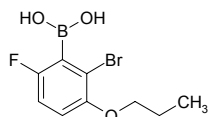
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	592129	272.94

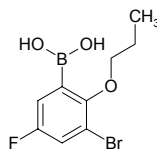


**Sigma - Aldrich
SAR Boronic Acids**

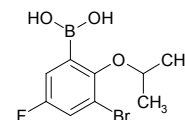
Brand	Catalog Number	MW
ALDRICH	645729	276.90

**Sigma - Aldrich
SAR Boronic Acids**

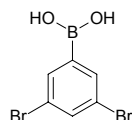
Brand	Catalog Number	MW
ALDRICH	565784	276.90

**Sigma - Aldrich
SAR Boronic Acids**

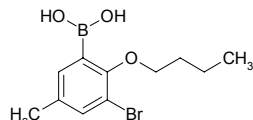
Brand	Catalog Number	MW
ALDRICH	542520	276.90

**Sigma - Aldrich
SAR Boronic Acids**

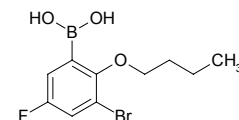
Brand	Catalog Number	MW
ALDRICH	499501	279.72

**Sigma - Aldrich
SAR Boronic Acids**

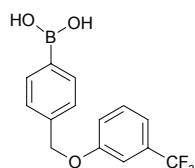
Brand	Catalog Number	MW
ALDRICH	596841	286.96

**Sigma - Aldrich
SAR Boronic Acids**

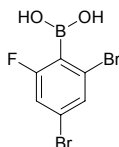
Brand	Catalog Number	MW
ALDRICH	542539	290.93

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	651222	296.06

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	651257	297.71

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	595756	300.95

