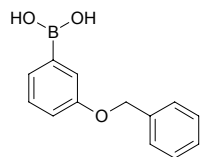




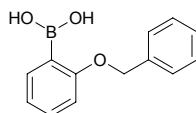
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	526339	228.06



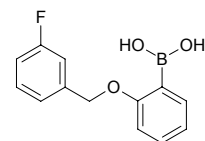
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	521337	228.06



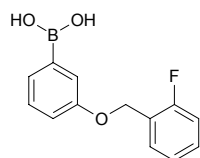
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	657492	246.05



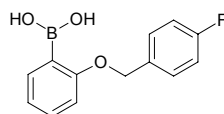
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	652113	246.05



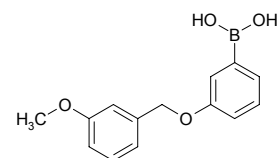
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	652105	246.05



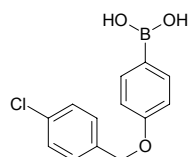
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	657395	258.08



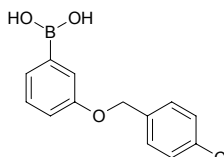
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	651044	262.50



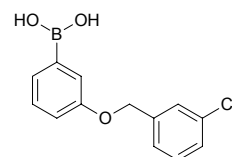
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	651036	262.50



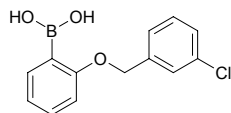
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	645265	262.50



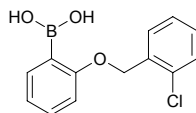
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	645249	262.50



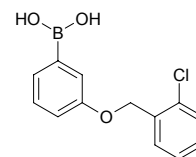
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	639443	262.50



**Sigma - Aldrich
SAR Boronic Acids**

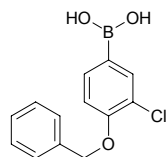
Brand	Catalog Number	MW
ALDRICH	635731	262.50





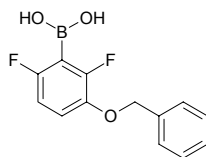
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	595616	262.50



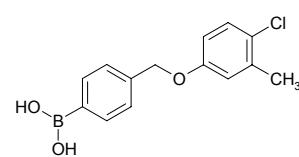
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	635707	264.04



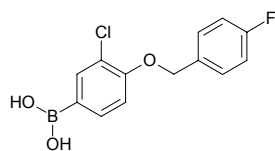
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	645931	276.53



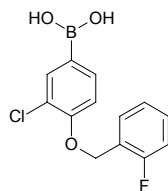
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	645397	280.49



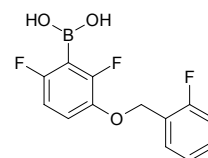
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	639575	280.49



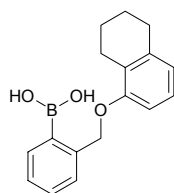
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	651346	282.03



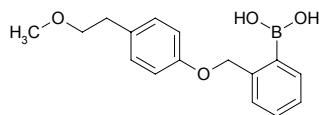
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	652229	282.15



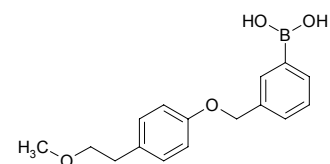
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	658065	286.14



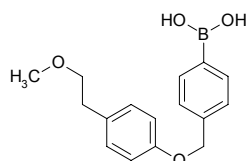
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	657506	286.14



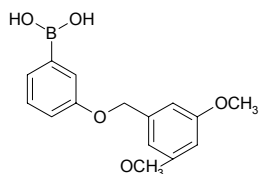
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	652083	286.14



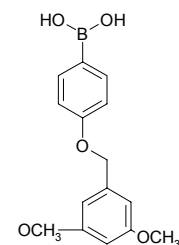
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	635766	288.11



**Sigma - Aldrich
SAR Boronic Acids**

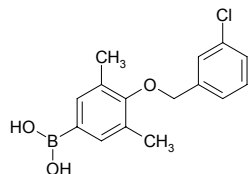
Brand	Catalog Number	MW
ALDRICH	635758	288.11





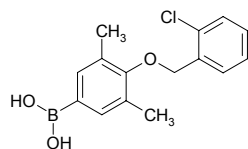
Sigma - Aldrich SAR Boronic Acids

Brand	Catalog Number	MW
ALDRICH	645443	290.56



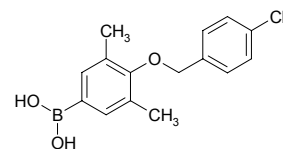
Sigma - Aldrich SAR Boronic Acids

Brand	Catalog Number	MW
ALDRICH	645435	290.56



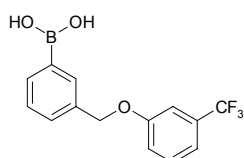
Sigma - Aldrich SAR Boronic Acids

Brand	Catalog Number	MW
ALDRICH	645354	290.56



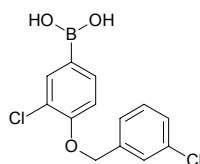
Sigma - Aldrich SAR Boronic Acids

Brand	Catalog Number	MW
ALDRICH	651311	296.06



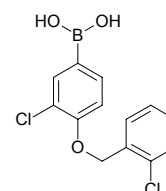
Sigma - Aldrich SAR Boronic Acids

Brand	Catalog Number	MW
ALDRICH	639540	296.95



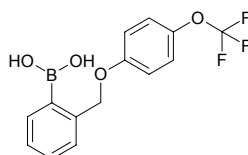
Sigma - Aldrich SAR Boronic Acids

Brand	Catalog Number	MW
ALDRICH	639532	296.95



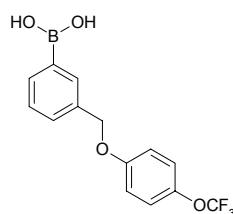
Sigma - Aldrich SAR Boronic Acids

Brand	Catalog Number	MW
ALDRICH	652156	312.06



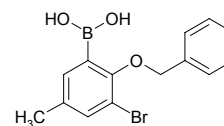
Sigma - Aldrich SAR Boronic Acids

Brand	Catalog Number	MW
ALDRICH	652075	312.06



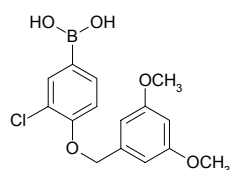
Sigma - Aldrich SAR Boronic Acids

Brand	Catalog Number	MW
ALDRICH	639435	320.98



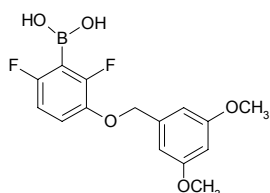
Sigma - Aldrich SAR Boronic Acids

Brand	Catalog Number	MW
ALDRICH	639516	322.56



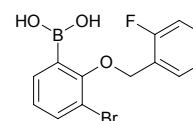
Sigma - Aldrich SAR Boronic Acids

Brand	Catalog Number	MW
ALDRICH	651109	324.09



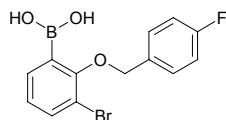
Sigma - Aldrich SAR Boronic Acids

Brand	Catalog Number	MW
ALDRICH	645834	324.94

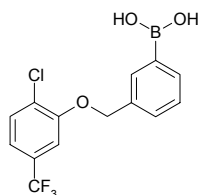


**Sigma - Aldrich
SAR Boronic Acids**

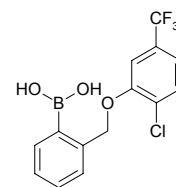
Brand	Catalog Number	MW
ALDRICH	645761	324.94

**Sigma - Aldrich
SAR Boronic Acids**

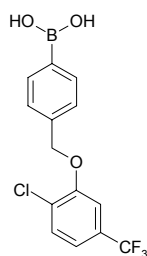
Brand	Catalog Number	MW
ALDRICH	657530	330.50

**Sigma - Aldrich
SAR Boronic Acids**

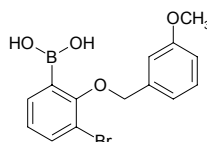
Brand	Catalog Number	MW
ALDRICH	652245	330.50

**Sigma - Aldrich
SAR Boronic Acids**

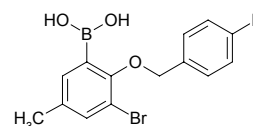
Brand	Catalog Number	MW
ALDRICH	652237	330.50

**Sigma - Aldrich
SAR Boronic Acids**

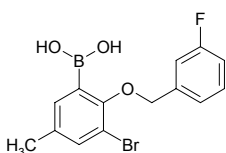
Brand	Catalog Number	MW
ALDRICH	645850	336.98

**Sigma - Aldrich
SAR Boronic Acids**

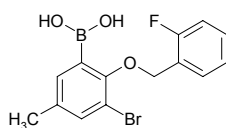
Brand	Catalog Number	MW
ALDRICH	645427	338.97

**Sigma - Aldrich
SAR Boronic Acids**

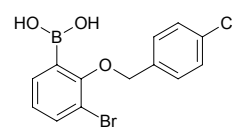
Brand	Catalog Number	MW
ALDRICH	645419	338.97

**Sigma - Aldrich
SAR Boronic Acids**

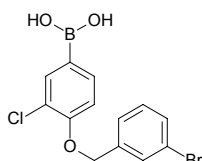
Brand	Catalog Number	MW
ALDRICH	645400	338.97

**Sigma - Aldrich
SAR Boronic Acids**

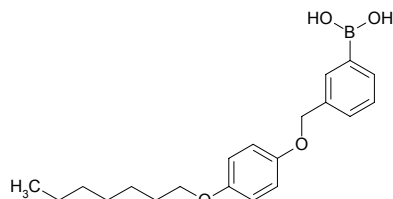
Brand	Catalog Number	MW
ALDRICH	645842	341.40

**Sigma - Aldrich
SAR Boronic Acids**

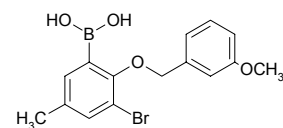
Brand	Catalog Number	MW
ALDRICH	639524	341.40

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	651214	342.25

**Sigma - Aldrich
SAR Boronic Acids**

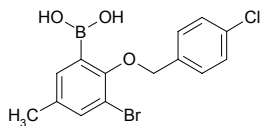
Brand	Catalog Number	MW
ALDRICH	639478	351.01





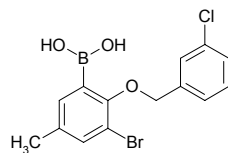
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	645508	355.43



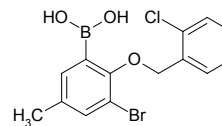
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	645494	355.43



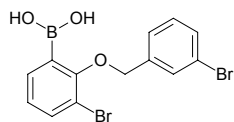
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	645486	355.43



**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	639567	385.85



**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	645478	399.88

