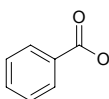




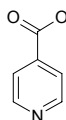
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	242381	122.12



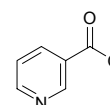
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	I17508	123.11



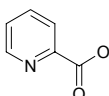
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	N7850	123.11



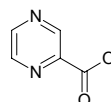
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	P42800	123.11



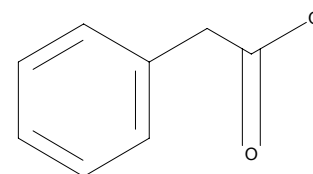
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	P56100	124.10



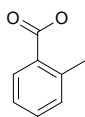
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	P16621	136.15



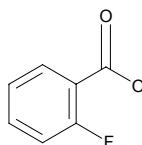
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	T36404	136.15



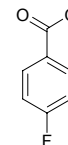
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	412244	140.12



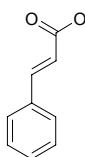
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	418846	140.12



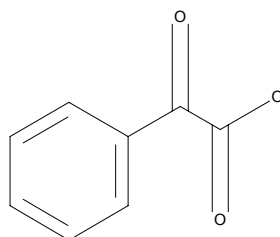
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	133760	148.16



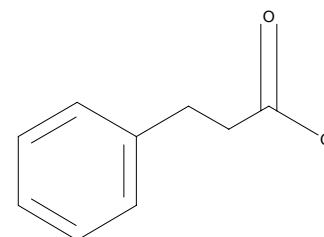
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	B13055	150.14



**Sigma - Aldrich
SAR COOH**

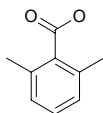
Brand	Catalog Number	MW
ALDRICH	135232	150.18





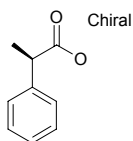
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	156906	150.18



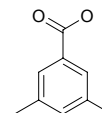
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	279897	150.18



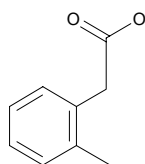
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	D149608	150.18



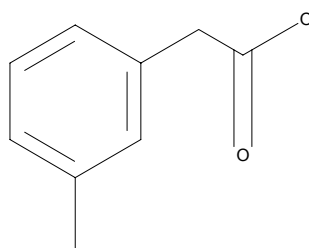
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	T38083	150.18



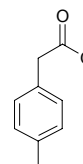
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	T38091	150.18



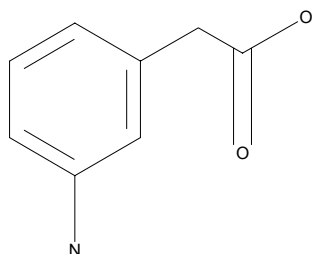
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	T38105	150.18



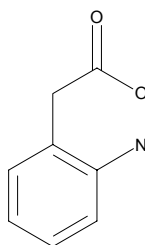
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	335444	151.17



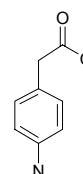
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	552836	151.17



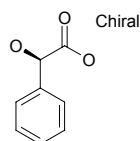
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	A71352	151.17



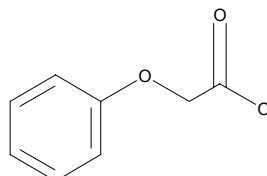
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	154210	152.15



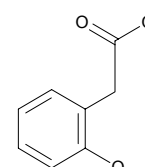
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	158518	152.15



**Sigma - Aldrich
SAR COOH**

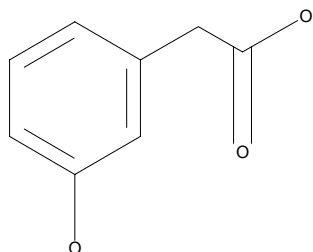
Brand	Catalog Number	MW
ALDRICH	H49804	152.15





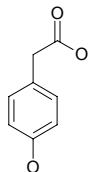
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	H49901	152.15



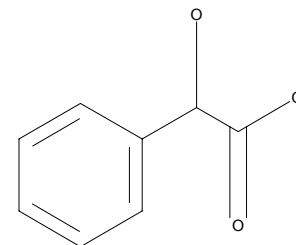
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	H50004	152.15



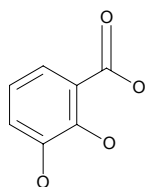
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	M2101	152.15



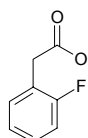
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	126209	154.12



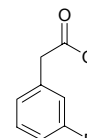
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	208949	154.14



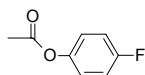
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	248045	154.14



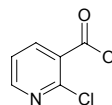
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	469238	154.14



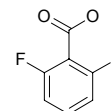
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	150339	157.56



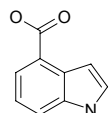
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	190039	158.11



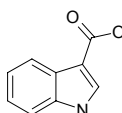
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	246263	161.16



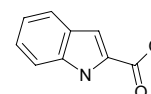
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	284734	161.16



**Sigma - Aldrich
SAR COOH**

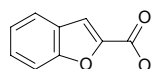
Brand	Catalog Number	MW
ALDRICH	I5109	161.16





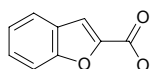
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	307270	162.15



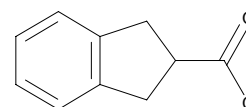
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	307270	162.15



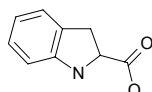
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	595551	162.19



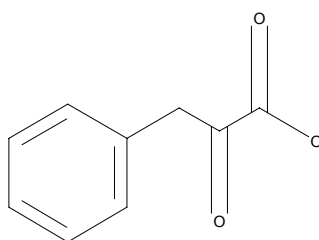
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	302244	163.18



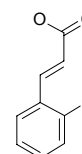
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	286958	164.16



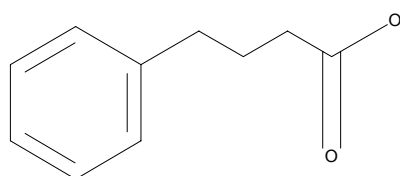
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	H22809	164.16



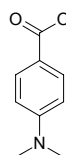
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	P21005	164.21



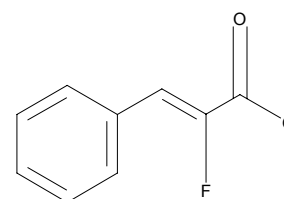
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	D139459	165.19



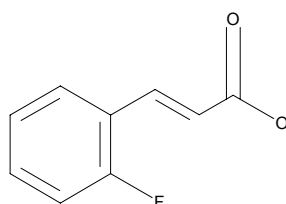
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	163848	166.15



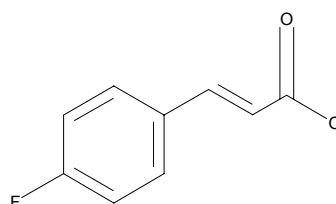
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	222712	166.15



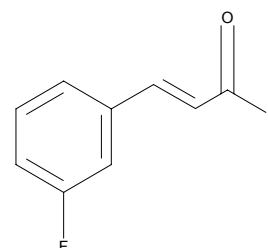
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	222720	166.15



**Sigma - Aldrich
SAR COOH**

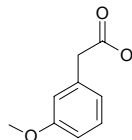
Brand	Catalog Number	MW
ALDRICH	290483	166.15





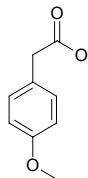
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	M19007	166.18



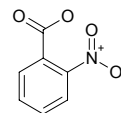
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	M19201	166.18



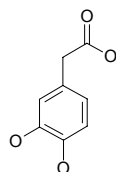
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	127698	167.12



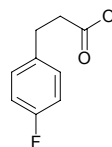
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	850217	168.15



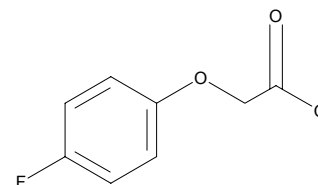
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	560502	168.17



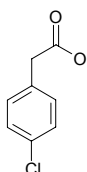
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	850276	170.14



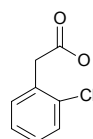
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	139262	170.60



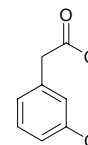
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	190632	170.60



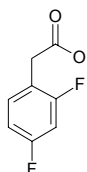
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	C63359	170.60



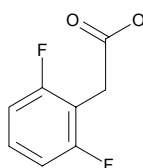
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	264474	172.13



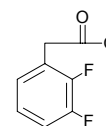
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	264482	172.13



**Sigma - Aldrich
SAR COOH**

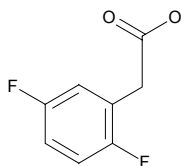
Brand	Catalog Number	MW
ALDRICH	290416	172.13





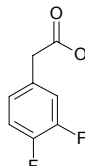
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	290424	172.13



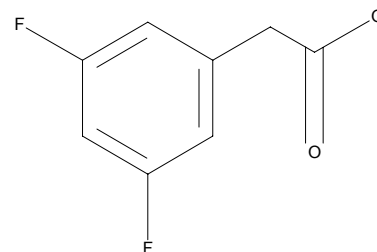
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	290432	172.13



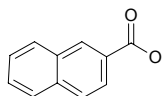
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	290440	172.13



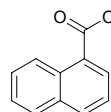
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	180246	172.19



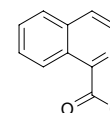
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	N1909	172.19



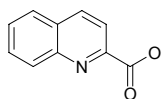
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	150134	173.17



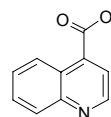
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	160660	173.17



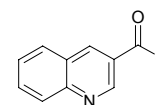
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	174823	173.17



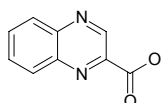
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	177148	173.17



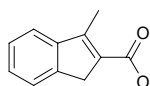
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	293407	174.16



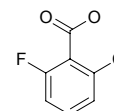
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	188425	174.20



**Sigma - Aldrich
SAR COOH**

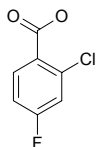
Brand	Catalog Number	MW
ALDRICH	178039	174.56





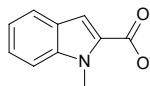
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	297836	174.56



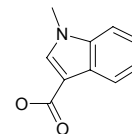
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	134155	175.19



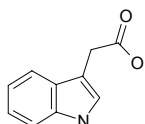
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	465313	175.19



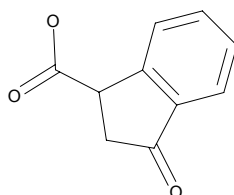
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	13750	175.19



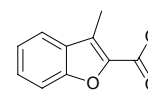
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	410772	176.17



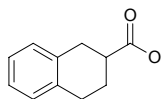
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	530611	176.17



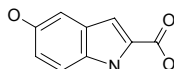
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	178683	176.22



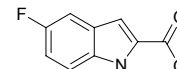
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	143510	177.16



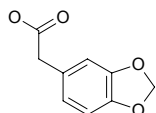
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	265128	179.15



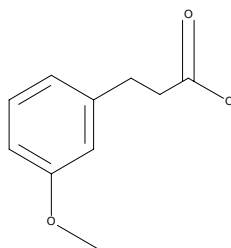
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	329673	180.16



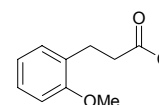
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	349763	180.21



**Sigma - Aldrich
SAR COOH**

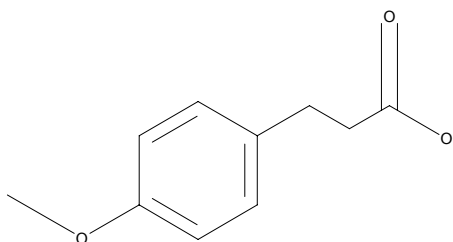
Brand	Catalog Number	MW
ALDRICH	M23500	180.21





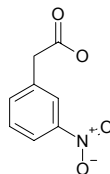
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	M23527	180.21



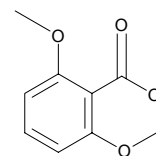
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	103608	181.15



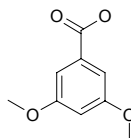
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	D131601	182.18



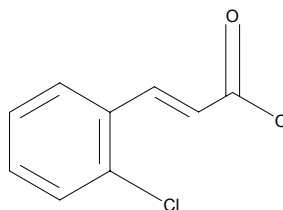
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	D132004	182.18



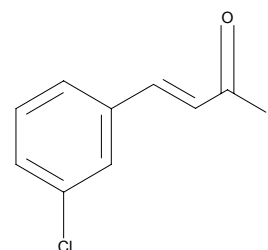
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	C31392	182.61



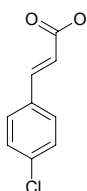
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	C31406	182.61



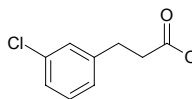
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	C31600	182.61



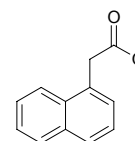
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	631302	184.62



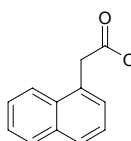
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	317918	186.21



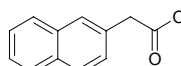
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	N3804	186.21



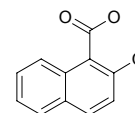
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	N4002	186.21



**Sigma - Aldrich
SAR COOH**

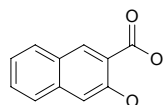
Brand	Catalog Number	MW
ALDRICH	H45809	188.18





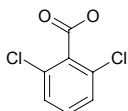
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	H46007	188.18



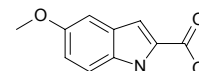
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	D57450	191.01



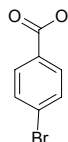
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	M14951	191.19



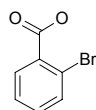
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	108510	201.02



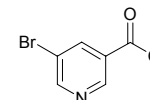
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	137677	201.02



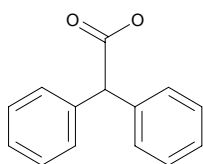
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	228435	202.01



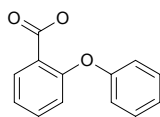
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	D204307	212.25



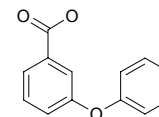
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	153176	214.22



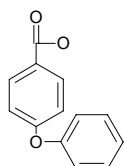
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	190276	214.22



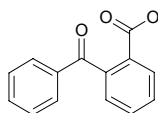
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	246182	214.22



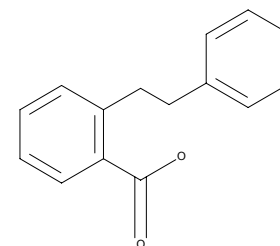
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	B12385	226.23



**Sigma - Aldrich
SAR COOH**

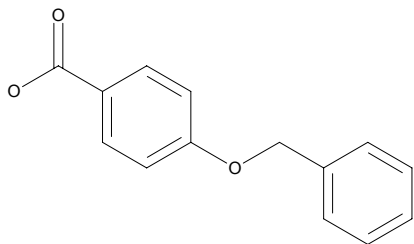
Brand	Catalog Number	MW
ALDRICH	151513	226.28





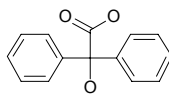
Sigma - Aldrich
SAR COOH

Brand	Catalog Number	MW
ALDRICH	512354	228.25



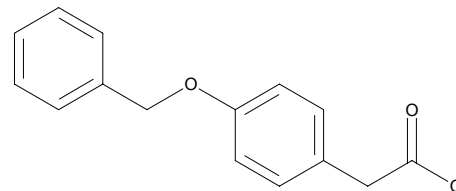
Sigma - Aldrich
SAR COOH

Brand	Catalog Number	MW
ALDRICH	B5194	228.25



Sigma - Aldrich
SAR COOH

Brand	Catalog Number	MW
ALDRICH	555398	242.28



Sigma - Aldrich
SAR COOH

Brand	Catalog Number	MW
ALDRICH	514594	291.15

