



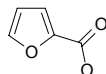
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	C116602	86.09



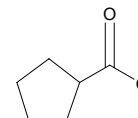
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	F20505	112.09



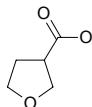
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	C112003	114.15



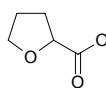
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	339954	116.12



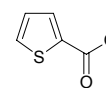
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	341517	116.12



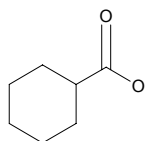
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	T32603	128.15



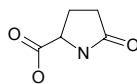
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	101834	128.17



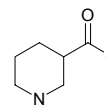
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	292915	129.12



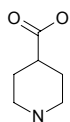
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	211672	129.16



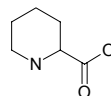
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	I18008	129.16



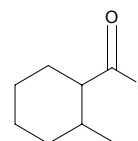
**Sigma - Aldrich
SAR COOH**

Brand	Catalog Number	MW
ALDRICH	P45850	129.16



**Sigma - Aldrich
SAR COOH**

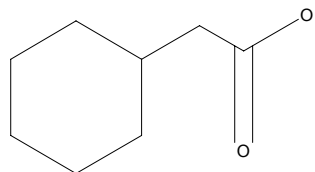
Brand	Catalog Number	MW
ALDRICH	330604	142.20





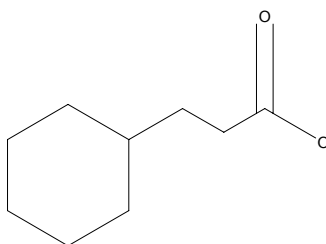
Sigma - Aldrich
SAR COOH

Brand	Catalog Number	MW
ALDRICH	C104507	142.20



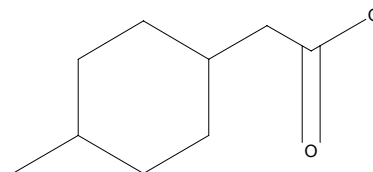
Sigma - Aldrich
SAR COOH

Brand	Catalog Number	MW
ALDRICH	161470	156.23



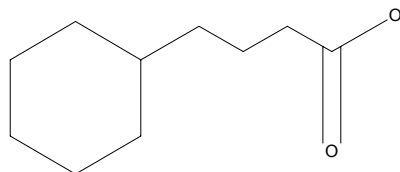
Sigma - Aldrich
SAR COOH

Brand	Catalog Number	MW
ALDRICH	332852	156.23



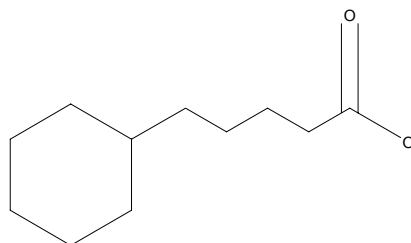
Sigma - Aldrich
SAR COOH

Brand	Catalog Number	MW
ALDRICH	127612	170.25



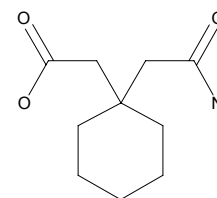
Sigma - Aldrich
SAR COOH

Brand	Catalog Number	MW
ALDRICH	331562	184.28



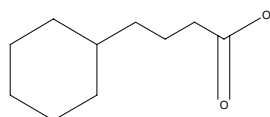
Sigma - Aldrich
SAR COOH

Brand	Catalog Number	MW
ALDRICH	553735	199.25



Sigma - Aldrich
SAR COOH

Brand	Catalog Number	MW
ALDRICH	228311	210.36



K



Sigma - Aldrich
SAR COOH

Brand	Catalog Number	MW
ALDRICH	333840	224.35

