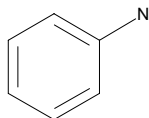




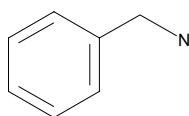
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	242284	93.13



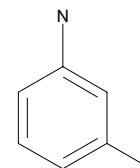
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	B16305	107.16



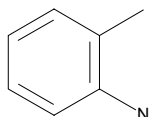
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	511218	107.16



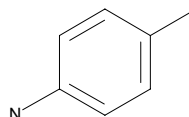
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	185426	107.16



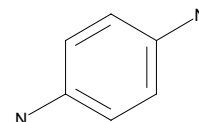
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	236314	107.16



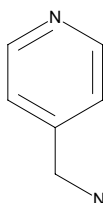
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	275158	108.14



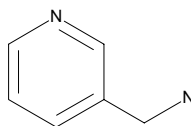
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	A65603	108.14



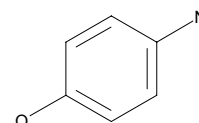
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	A65409	108.14



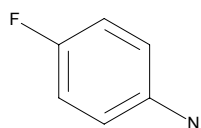
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	245992	109.13



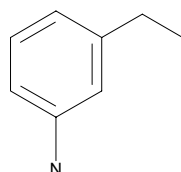
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	F3800	111.12



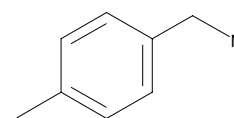
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	175498	121.18



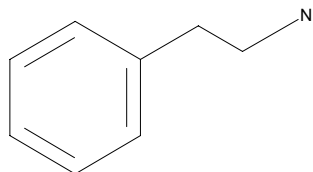
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	M31201	121.18

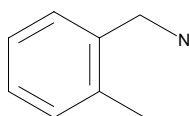


**Sigma - Aldrich
SAR Amines**

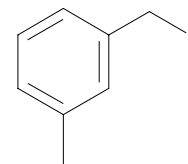
Brand	Catalog Number	MW
Aldrich	407267	121.18

**Sigma - Aldrich
SAR Amines**

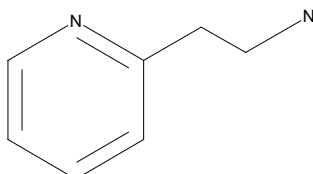
Brand	Catalog Number	MW
Aldrich	127132	121.18

**Sigma - Aldrich
SAR Amines**

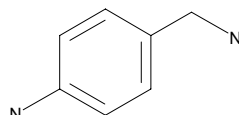
Brand	Catalog Number	MW
Aldrich	126829	121.18

**Sigma - Aldrich
SAR Amines**

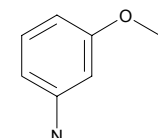
Brand	Catalog Number	MW
Aldrich	A55306	122.17

**Sigma - Aldrich
SAR Amines**

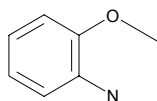
Brand	Catalog Number	MW
Aldrich	368466	122.17

**Sigma - Aldrich
SAR Amines**

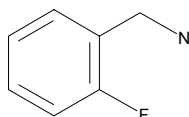
Brand	Catalog Number	MW
Aldrich	A88204	123.16

**Sigma - Aldrich
SAR Amines**

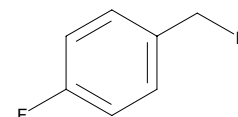
Brand	Catalog Number	MW
Aldrich	A88182	123.16

**Sigma - Aldrich
SAR Amines**

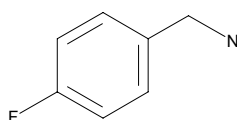
Brand	Catalog Number	MW
Aldrich	162485	125.15

**Sigma - Aldrich
SAR Amines**

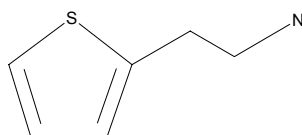
Brand	Catalog Number	MW
Aldrich	162493	125.15

**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	162493	125.15

**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	423270	127.21

**Sigma - Aldrich
SAR Amines**

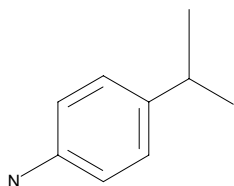
Brand	Catalog Number	MW
Aldrich	459917	127.57





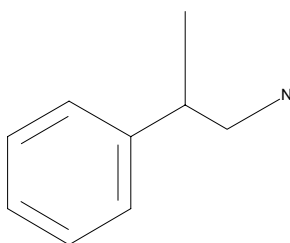
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	175439	135.21



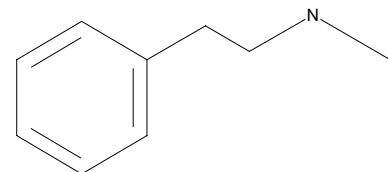
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	180076	135.21



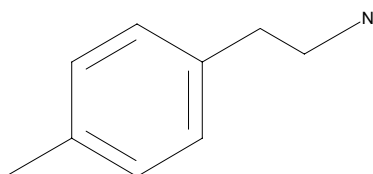
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	M68423	135.21



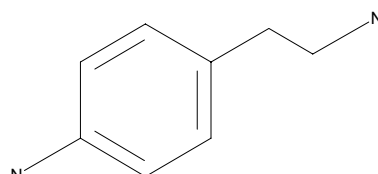
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	132020	135.21



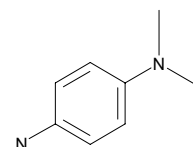
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	123056	136.20



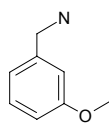
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	193992	136.20



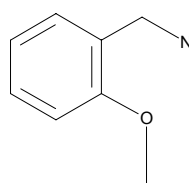
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	159891	137.18



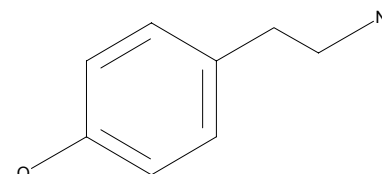
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	159883	137.18



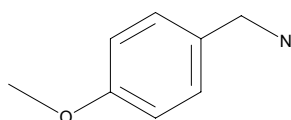
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	T90344	137.18



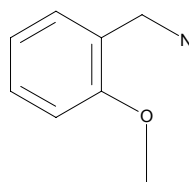
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	M11103	137.18



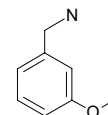
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	159883	137.18



**Sigma - Aldrich
SAR Amines**

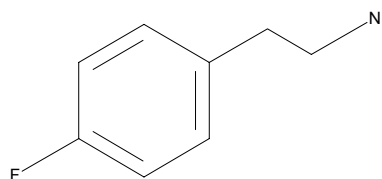
Brand	Catalog Number	MW
Aldrich	159891	137.18





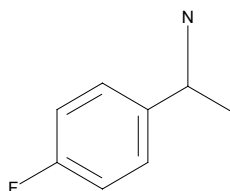
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	361828	139.17



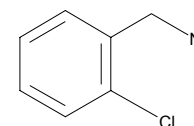
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	154040	139.17



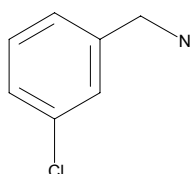
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	C27204	141.60



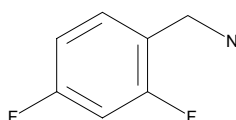
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	127167	141.60



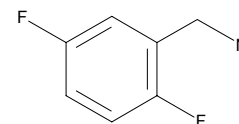
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	264377	143.14



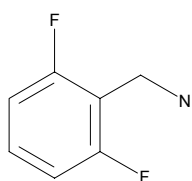
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	264385	143.14



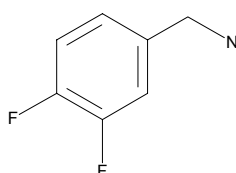
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	264393	143.14



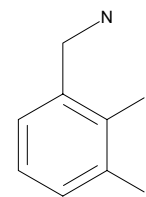
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	264407	143.14



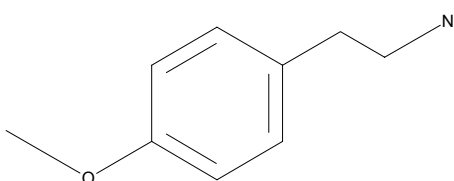
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	538612	143.14



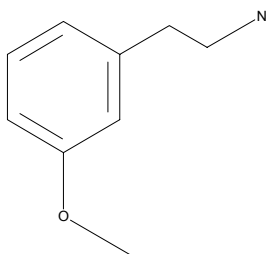
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	187305	151.21



**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	270229	151.21



**Sigma - Aldrich
SAR Amines**

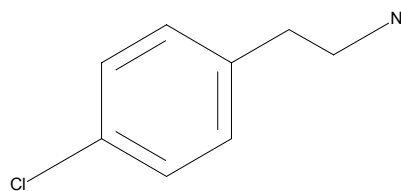
Brand	Catalog Number	MW
Aldrich	476021	151.21





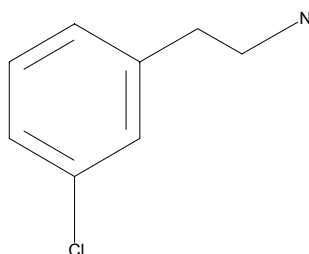
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	C65408	155.63



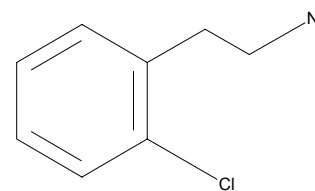
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	340847	155.63



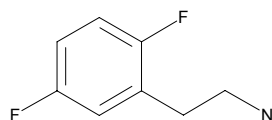
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	243736	155.63



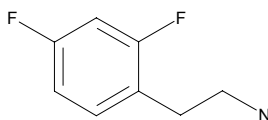
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
SALOR	L300276	157.16



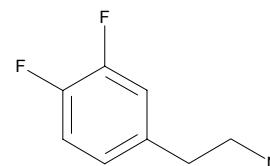
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
SALOR	L300241	157.16



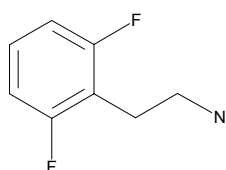
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
SALOR	L300233	157.16



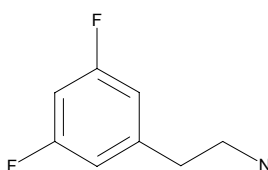
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
SALOR	L300187	157.16



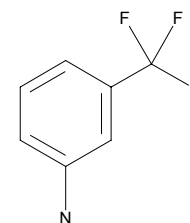
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
SALOR	L300179	157.16



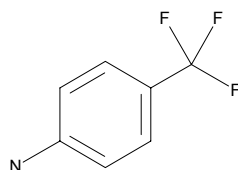
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	A41801	161.13



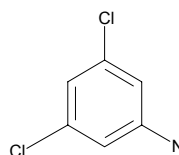
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	224936	161.13



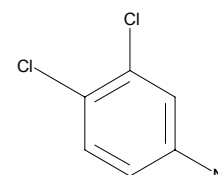
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	D55792	162.02



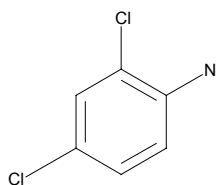
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	437778	162.02

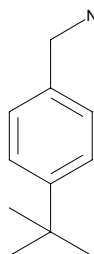


**Sigma - Aldrich SAR Amines**

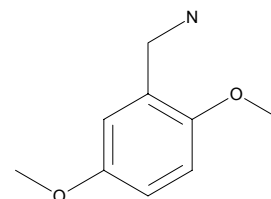
Brand	Catalog Number	MW
Aldrich	112151	162.02

**Sigma - Aldrich SAR Amines**

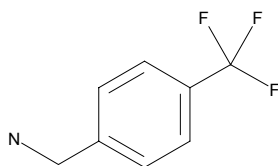
Brand	Catalog Number	MW
Aldrich	631280	163.26

**Sigma - Aldrich SAR Amines**

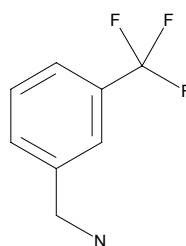
Brand	Catalog Number	MW
Aldrich	540315	167.21

**Sigma - Aldrich SAR Amines**

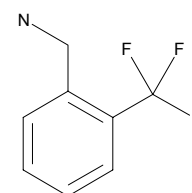
Brand	Catalog Number	MW
Aldrich	263508	175.15

**Sigma - Aldrich SAR Amines**

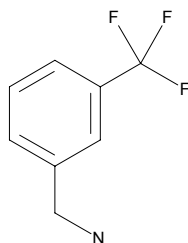
Brand	Catalog Number	MW
Aldrich	263494	175.15

**Sigma - Aldrich SAR Amines**

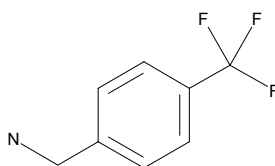
Brand	Catalog Number	MW
Aldrich	263486	175.15

**Sigma - Aldrich SAR Amines**

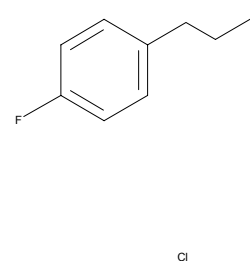
Brand	Catalog Number	MW
Aldrich	263494	175.15

**Sigma - Aldrich SAR Amines**

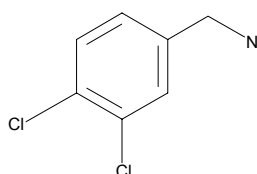
Brand	Catalog Number	MW
Aldrich	263508	175.15

**Sigma - Aldrich SAR Amines**

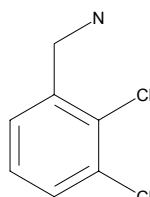
Brand	Catalog Number	MW
Aldrich	180394	175.63

**Sigma - Aldrich SAR Amines**

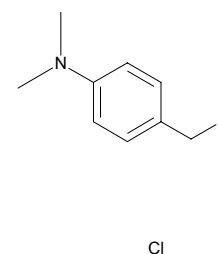
Brand	Catalog Number	MW
Aldrich	D58600	176.05

**Sigma - Aldrich SAR Amines**

Brand	Catalog Number	MW
Aldrich	539341	176.05

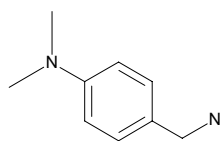
**Sigma - Aldrich SAR Amines**

Brand	Catalog Number	MW
Aldrich	285633	186.69



**Sigma - Aldrich
SAR Amines**

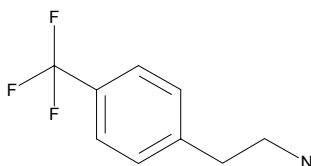
Brand	Catalog Number	MW
Aldrich	285633	186.69



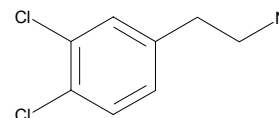
Cl

**Sigma - Aldrich
SAR Amines**

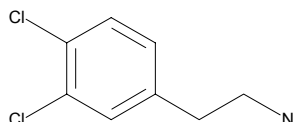
Brand	Catalog Number	MW
SALOR	L300225	189.18

**Sigma - Aldrich
SAR Amines**

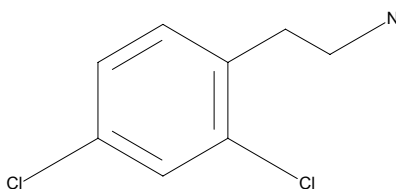
Brand	Catalog Number	MW
SALOR	L300292	190.07

**Sigma - Aldrich
SAR Amines**

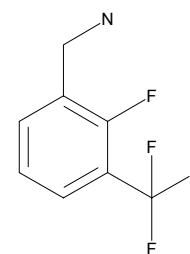
Brand	Catalog Number	MW
Aldrich	560421	190.07

**Sigma - Aldrich
SAR Amines**

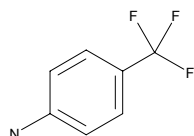
Brand	Catalog Number	MW
Aldrich	357502	190.07

**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	537861	193.15

**Sigma - Aldrich
SAR Amines**

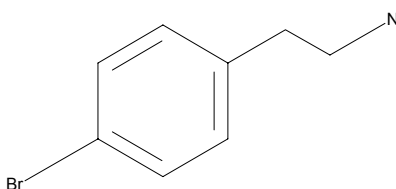
Brand	Catalog Number	MW
Aldrich	172871	197.59



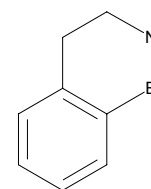
Cl

**Sigma - Aldrich
SAR Amines**

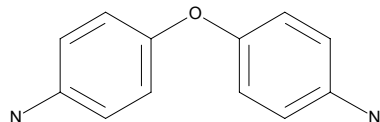
Brand	Catalog Number	MW
Aldrich	189308	200.08

**Sigma - Aldrich
SAR Amines**

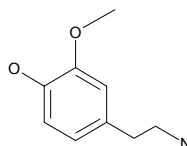
Brand	Catalog Number	MW
Aldrich	534099	200.08

**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	248398	200.24

**Sigma - Aldrich
SAR Amines**

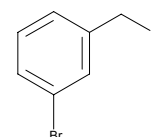
Brand	Catalog Number	MW
Aldrich	164313	203.67



Cl

**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	316288	222.51

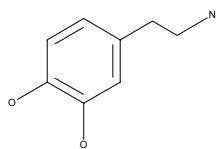


Cl



Sigma - Aldrich
SAR Amines

Brand	Catalog Number	MW
Aldrich	161136	234.09



Br