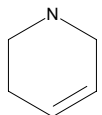




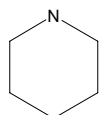
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	134759	83.13



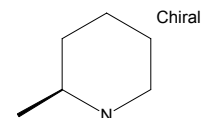
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	411027	85.15



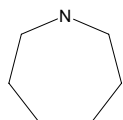
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	522902	99.18



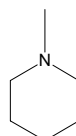
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	H10401	99.18



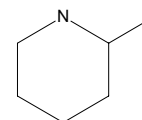
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	M72609	99.18



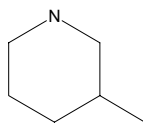
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	M72803	99.18



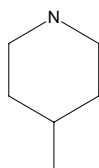
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	M73001	99.18



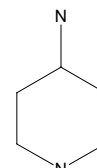
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	M73206	99.18



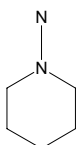
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	561479	100.16



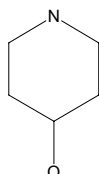
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	A75900	100.16



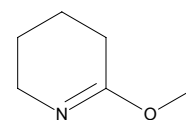
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	128775	101.15



**Sigma - Aldrich Building Blocks**

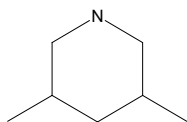
Brand	Catalog Number	MW
Aldrich	539198	113.16





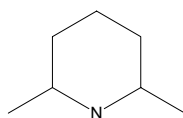
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	186104	113.20



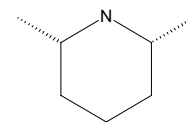
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	540811	113.20



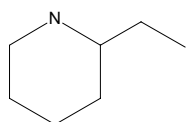
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	D180300	113.20



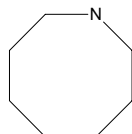
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	E45805	113.20



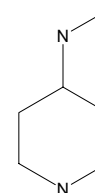
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	H2155	113.20



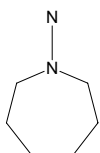
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	590495	114.19



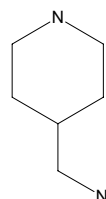
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	A56450	114.19



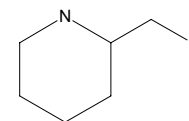
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	A65158	114.19



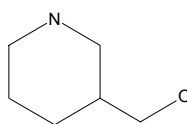
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	155225	115.18



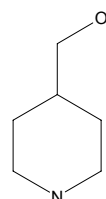
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	155233	115.18



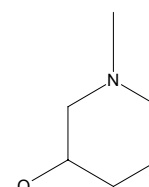
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	497312	115.18



**Sigma - Aldrich Building Blocks**

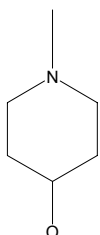
Brand	Catalog Number	MW
Aldrich	H42001	115.18





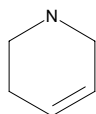
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	H42206	115.18



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	439266	119.60

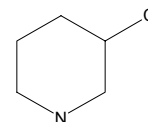


Cl



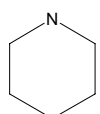
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	595365	119.60



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	P46105	121.61

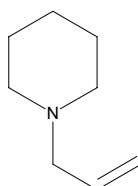


Cl



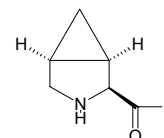
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	586463	125.22



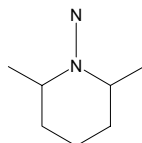
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	391190	127.14



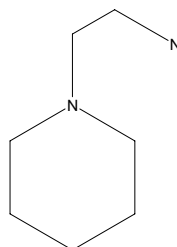
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	133450	128.22



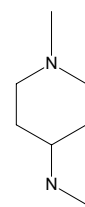
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	141666	128.22



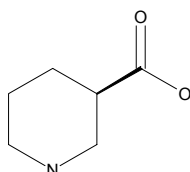
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	221406	128.22



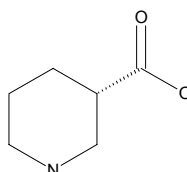
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	535664	129.16



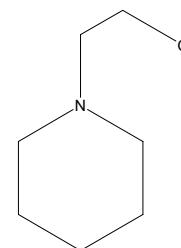
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	535672	129.16



**Sigma - Aldrich Building Blocks**

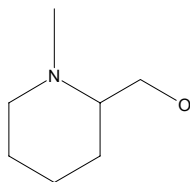
Brand	Catalog Number	MW
Aldrich	116068	129.20





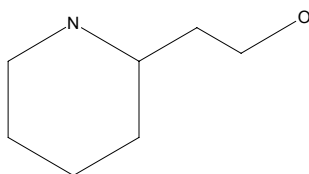
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	155241	129.20



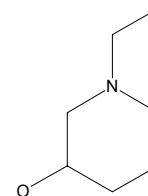
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	433594	129.20



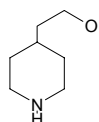
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	E32207	129.20



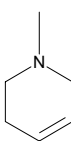
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	P46156	129.20



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	332380	133.62

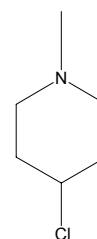


Cl



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	C56301	133.62



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	293369	136.63

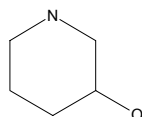


Cl



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	174416	137.61

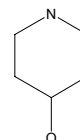


Cl



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	391468	137.61

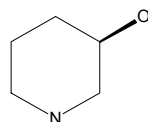


Cl



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	410462	137.61

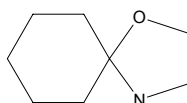


Cl



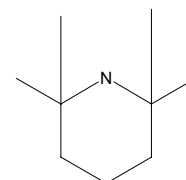
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	555800	141.21



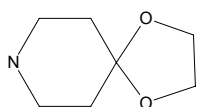
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	115754	141.26

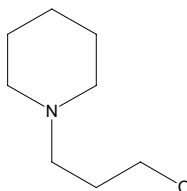


**Sigma - Aldrich Building Blocks**

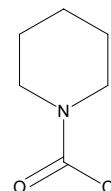
Brand	Catalog Number	MW
Aldrich	178365	143.19

**Sigma - Aldrich Building Blocks**

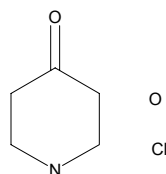
Brand	Catalog Number	MW
Aldrich	152935	143.23

**Sigma - Aldrich Building Blocks**

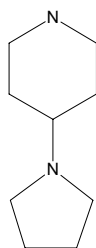
Brand	Catalog Number	MW
Aldrich	518522	147.61

**Sigma - Aldrich Building Blocks**

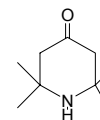
Brand	Catalog Number	MW
Aldrich	151769	153.61

**Sigma - Aldrich Building Blocks**

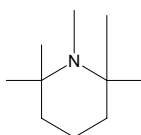
Brand	Catalog Number	MW
Aldrich	437352	154.26

**Sigma - Aldrich Building Blocks**

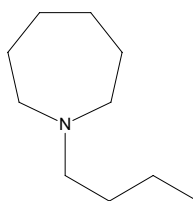
Brand	Catalog Number	MW
Aldrich	459119	155.24

**Sigma - Aldrich Building Blocks**

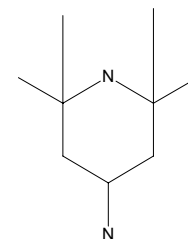
Brand	Catalog Number	MW
Aldrich	274631	155.29

**Sigma - Aldrich Building Blocks**

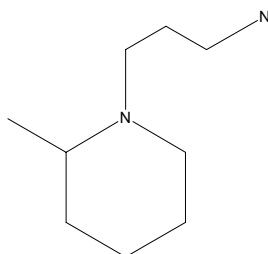
Brand	Catalog Number	MW
Aldrich	571954	155.29

**Sigma - Aldrich Building Blocks**

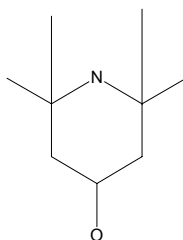
Brand	Catalog Number	MW
Aldrich	115738	156.27

**Sigma - Aldrich Building Blocks**

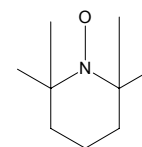
Brand	Catalog Number	MW
Aldrich	186112	156.27

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	115746	157.26

**Sigma - Aldrich Building Blocks**

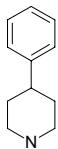
Brand	Catalog Number	MW
Aldrich	426369	157.26





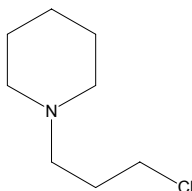
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	639869	161.25



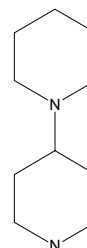
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	590533	161.68



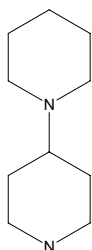
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	150053	168.28



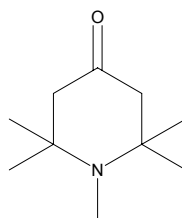
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	534498	168.28



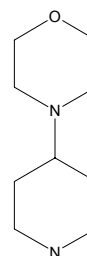
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	536903	169.27



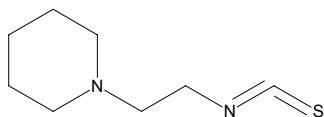
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	578045	170.26



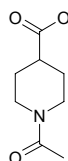
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	560456	170.28



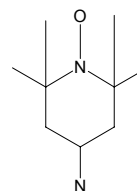
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	647756	171.20



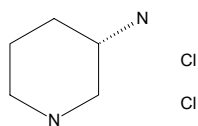
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	163945	172.27



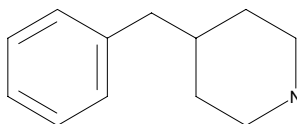
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	536660	173.09



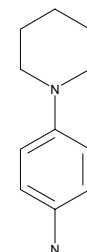
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	142360	175.28



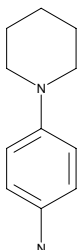
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	556629	176.26

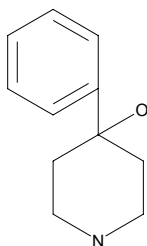


**Sigma - Aldrich Building Blocks**

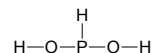
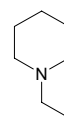
Brand	Catalog Number	MW
Aldrich	556629	176.26

**Sigma - Aldrich Building Blocks**

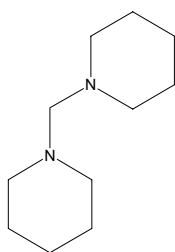
Brand	Catalog Number	MW
Aldrich	H52201	177.25

**Sigma - Aldrich Building Blocks**

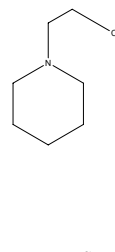
Brand	Catalog Number	MW
Aldrich	436178	179.20

**Sigma - Aldrich Building Blocks**

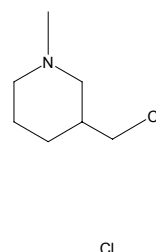
Brand	Catalog Number	MW
Aldrich	170208	182.31

**Sigma - Aldrich Building Blocks**

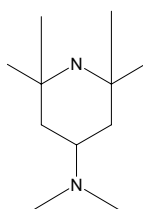
Brand	Catalog Number	MW
Aldrich	C42602	184.11

**Sigma - Aldrich Building Blocks**

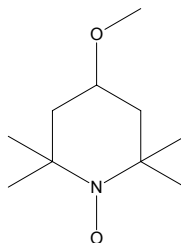
Brand	Catalog Number	MW
Aldrich	C53809	184.11

**Sigma - Aldrich Building Blocks**

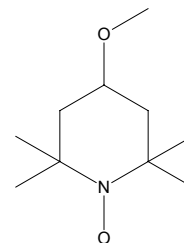
Brand	Catalog Number	MW
Aldrich	298573	184.33

**Sigma - Aldrich Building Blocks**

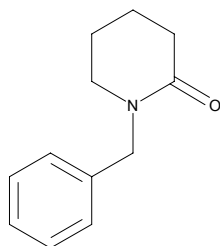
Brand	Catalog Number	MW
Aldrich	514586	187.28

**Sigma - Aldrich Building Blocks**

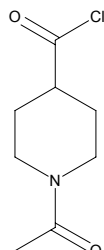
Brand	Catalog Number	MW
Aldrich	514586	187.28

**Sigma - Aldrich Building Blocks**

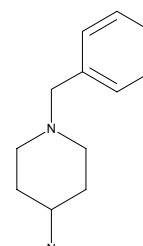
Brand	Catalog Number	MW
Aldrich	496898	189.26

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	632821	189.64

**Sigma - Aldrich Building Blocks**

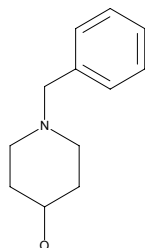
Brand	Catalog Number	MW
Aldrich	195812	190.29





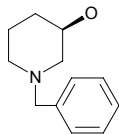
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	152986	191.28



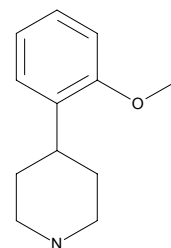
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	455172	191.28



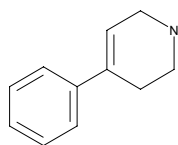
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	556483	191.28



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	134740	195.69

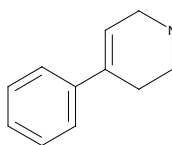


Cl



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	416835	195.69

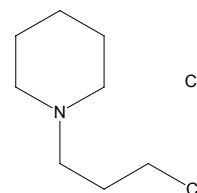


Cl



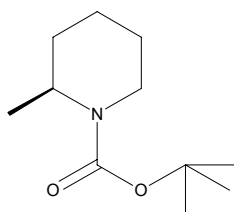
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	596310	198.14



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	522880	199.30

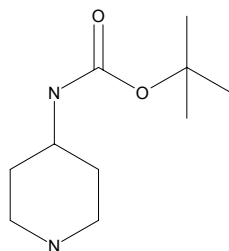


Chiral



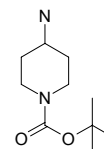
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	540935	200.28



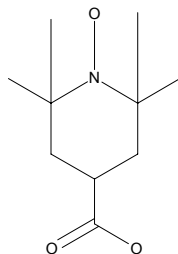
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	640042	200.28



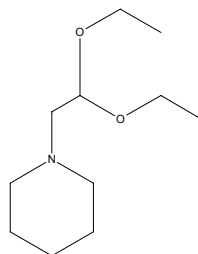
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	382000	201.27



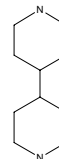
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	P46202	201.31



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	180742	204.75

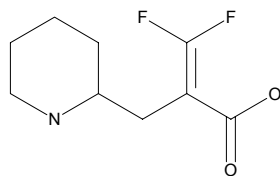


Cl



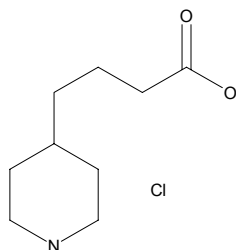
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	630470	205.21



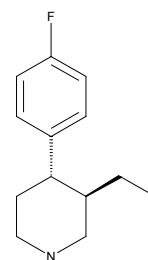
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	540838	207.70



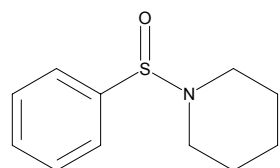
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	556807	209.27



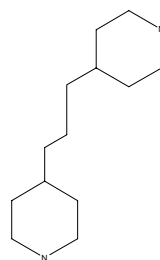
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	630233	209.31



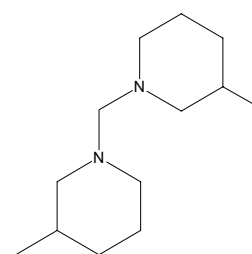
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	121207	210.37



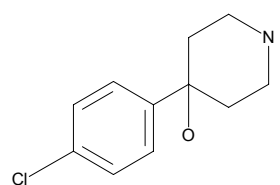
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	192252	210.37



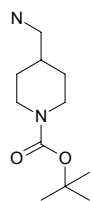
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	C66056	211.69



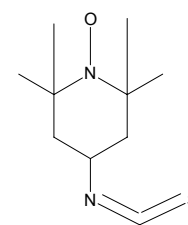
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	641472	214.31



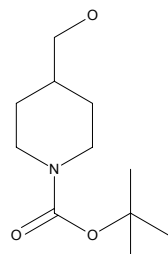
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	253286	214.33



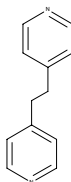
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	556017	215.29



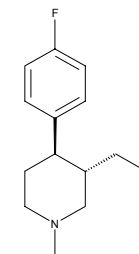
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	214140	220.70



**Sigma - Aldrich Building Blocks**

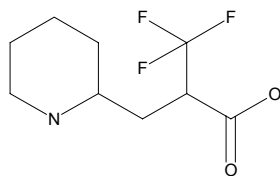
Brand	Catalog Number	MW
Aldrich	568058	223.29





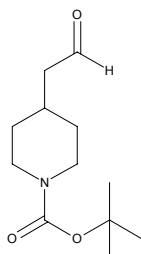
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	559369	225.21



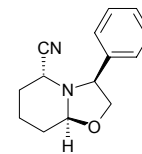
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	562300	227.31



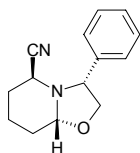
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	642002	228.30



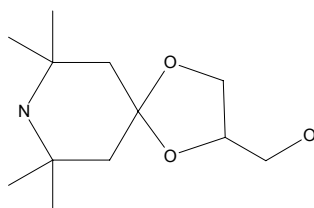
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	642118	228.30



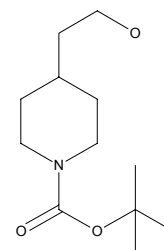
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	370002	229.32



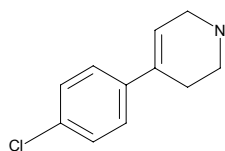
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	547247	229.32



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	C68407	230.14

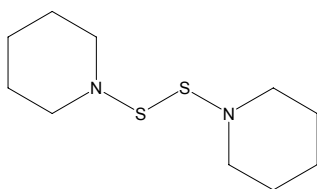


Cl



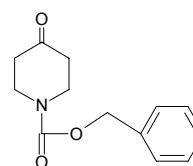
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	511153	232.41



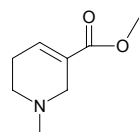
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	464643	233.27



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	444243	236.11

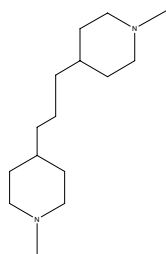


Br



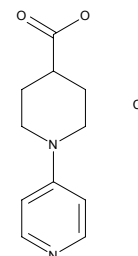
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	192260	238.42



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	572438	242.71

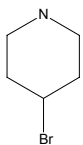


Cl



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	422320	244.96

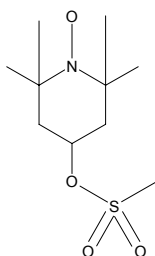


Br



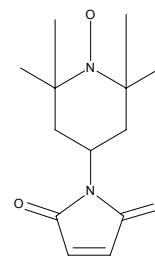
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	412686	251.35



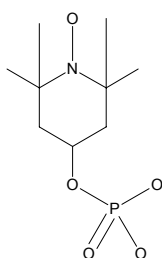
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	253359	252.32



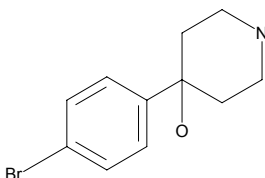
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	253642	253.24



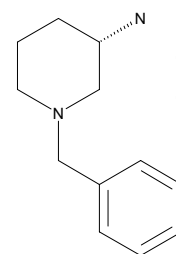
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	407445	256.14



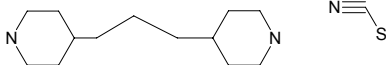
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	536679	263.21



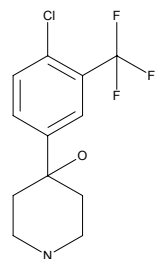
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	344850	269.46



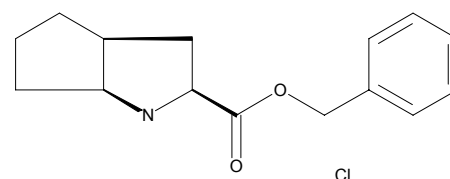
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	556823	279.69



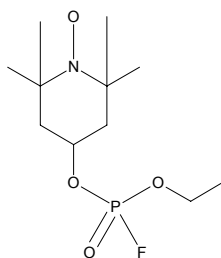
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	570907	281.79



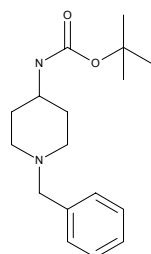
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	253553	283.28



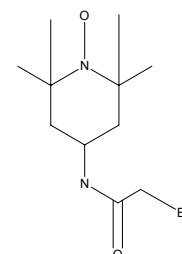
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	540943	290.41



**Sigma - Aldrich Building Blocks**

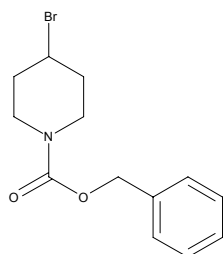
Brand	Catalog Number	MW
Aldrich	253227	293.21





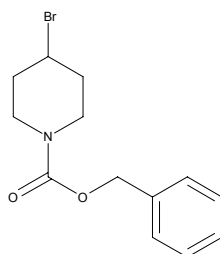
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	569224	298.18



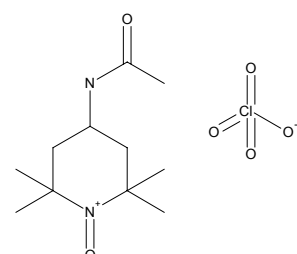
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	569224	298.18



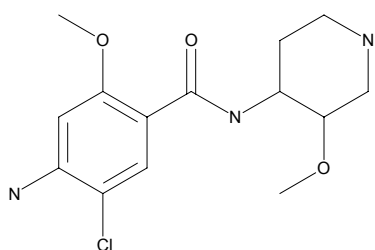
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	517704	312.75



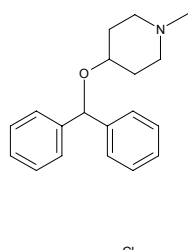
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	552380	313.79



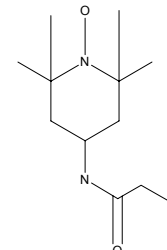
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	280550	317.86



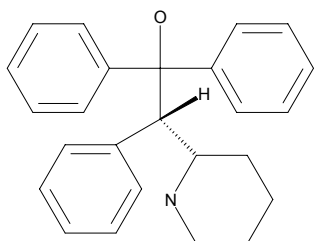
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	253367	340.21



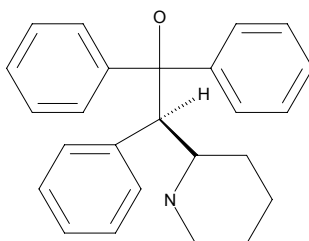
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	544817	357.50



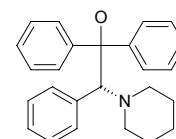
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	544825	357.50



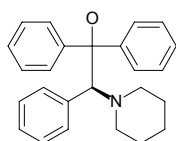
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	647462	357.50



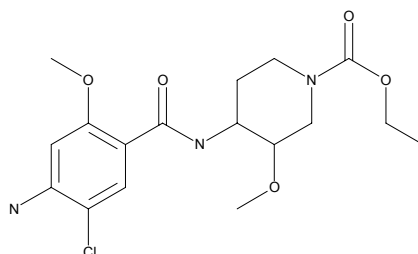
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	647470	357.50



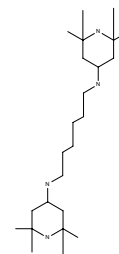
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	552399	385.85



### Sigma - Aldrich Building Blocks

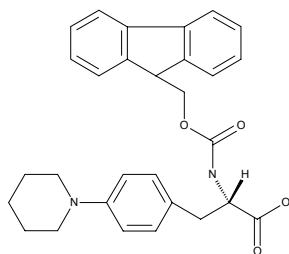
Brand	Catalog Number	MW
Aldrich	389420	394.69





**Sigma - Aldrich  
Building Blocks**

Brand	Catalog Number	MW
Aldrich	468754	470.57



**Sigma - Aldrich  
Building Blocks**

Brand	Catalog Number	MW
Aldrich	510319	484.42

