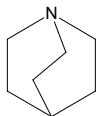




Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	197602	111.19



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	290734	112.18



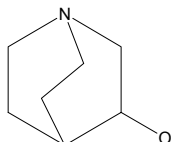
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	378704	125.22



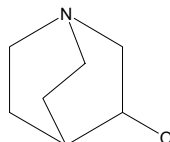
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	253340	127.19



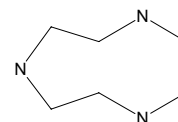
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	Q1875	127.19



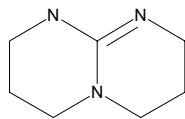
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	311308	129.21



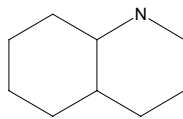
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	345571	139.20



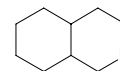
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	125741	139.24



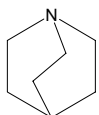
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	262595	139.24



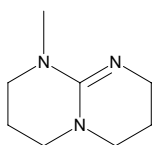
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	135917	147.65



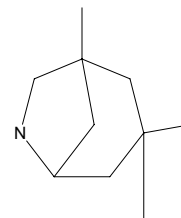
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	359505	153.23



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	194816	153.27

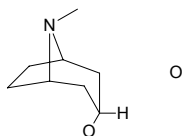


Cl



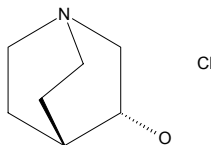
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	T89400	159.23



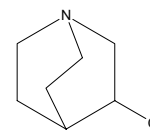
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	532908	163.65



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	Q1883	163.65

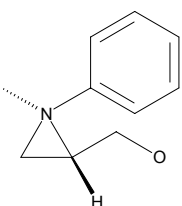


Cl



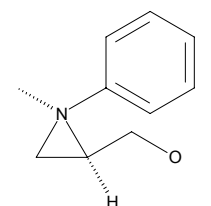
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	570532	165.24



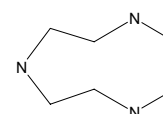
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	570567	165.24



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	329495	165.67

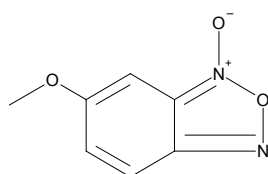


Cl



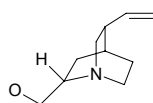
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	391409	166.14



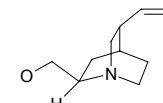
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	472530	167.25



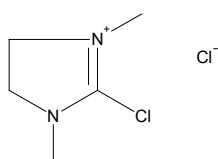
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	472549	167.25



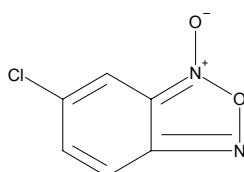
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	529249	169.05



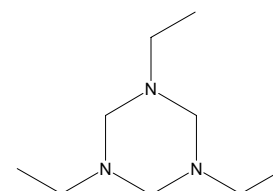
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	391395	170.56



Sigma - Aldrich Building Blocks

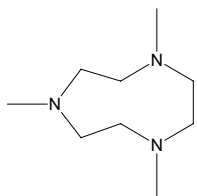
Brand	Catalog Number	MW
Aldrich	221430	171.29





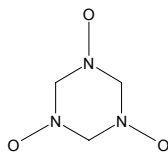
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	311294	171.29



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	238481	171.58

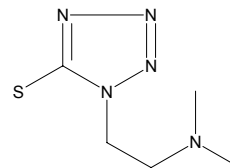


Cl



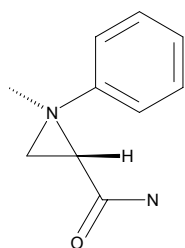
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	522333	173.24



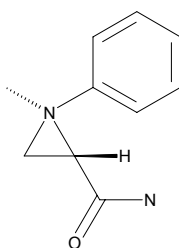
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	570524	178.24



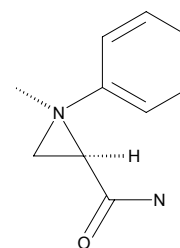
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	570524	178.24



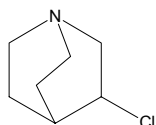
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	570559	178.24



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	125210	182.09

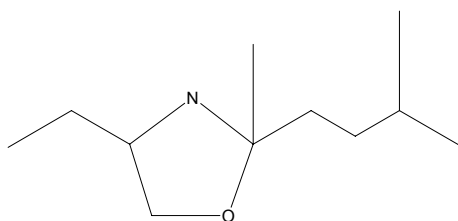


Cl



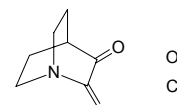
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	417815	185.31



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	M46128	191.66

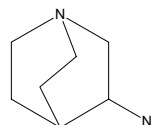


O
Cl



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	100358	199.12

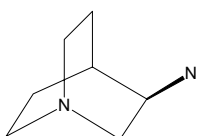


Cl
Cl



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	415715	199.12

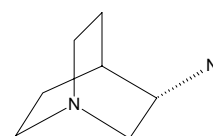


Cl
Cl



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	415723	199.12

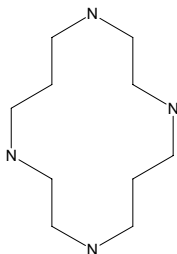


Cl
Cl



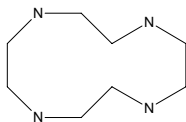
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	259160	200.33



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	433349	208.74

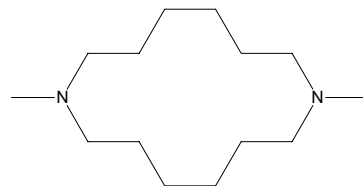


Cl



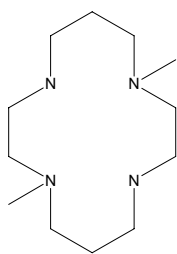
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	518859	226.41



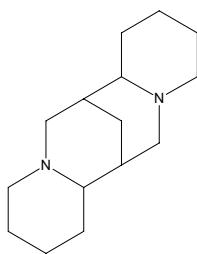
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	526169	228.38



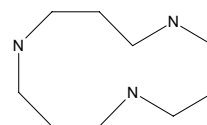
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	415316	234.39



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	323837	252.20



Br



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	473006	259.97

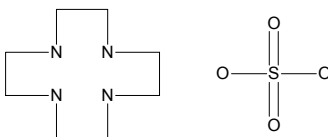


Br
Br



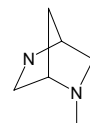
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	402621	270.35



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	476978	274.00

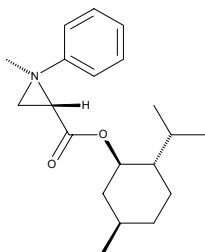


Br
Br



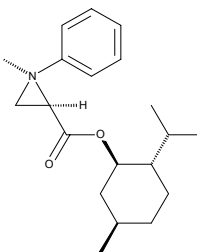
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	570516	317.48



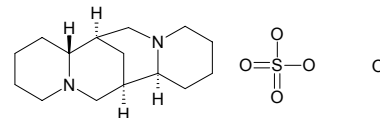
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	570540	317.48



Sigma - Aldrich Building Blocks

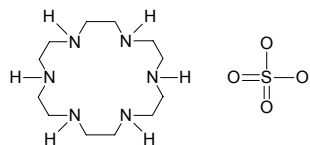
Brand	Catalog Number	MW
Aldrich	234664	350.48





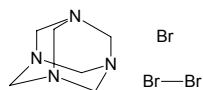
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	193933	356.49



**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	418501	380.91



**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	402508	392.96

