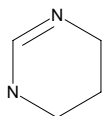




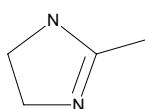
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	293334	84.12



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	375403	84.12



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	317454	98.15



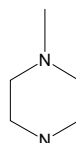
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	641065	100.12



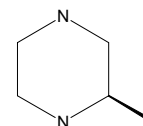
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	130001	100.16



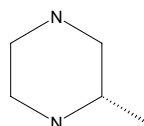
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	397164	100.16



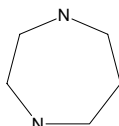
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	397172	100.16



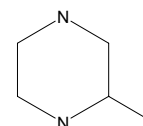
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	H16604	100.16



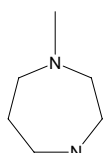
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	M72404	100.16



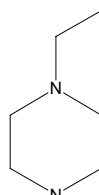
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	186090	114.19



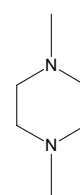
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	415308	114.19



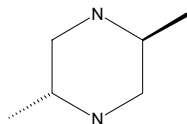
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	D179302	114.19

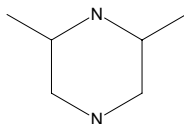


**Sigma - Aldrich Building Blocks**

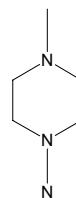
Brand	Catalog Number	MW
Aldrich	D179604	114.19

**Sigma - Aldrich Building Blocks**

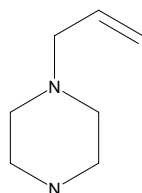
Brand	Catalog Number	MW
Aldrich	D179809	114.19

**Sigma - Aldrich Building Blocks**

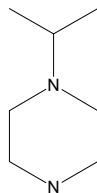
Brand	Catalog Number	MW
Aldrich	255688	115.18

**Sigma - Aldrich Building Blocks**

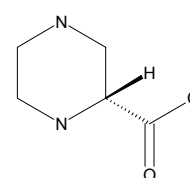
Brand	Catalog Number	MW
Aldrich	566888	126.20

**Sigma - Aldrich Building Blocks**

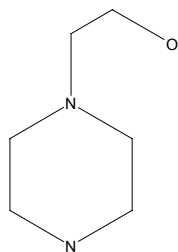
Brand	Catalog Number	MW
Aldrich	566853	128.22

**Sigma - Aldrich Building Blocks**

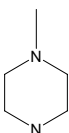
Brand	Catalog Number	MW
Aldrich	543543	130.15

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	H28807	130.19

**Sigma - Aldrich Building Blocks**

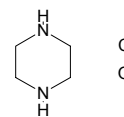
Brand	Catalog Number	MW
Aldrich	278793	136.63



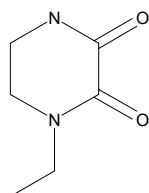
Cl

**Sigma - Aldrich Building Blocks**

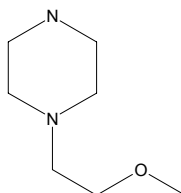
Brand	Catalog Number	MW
Aldrich	284904	140.61

**Sigma - Aldrich Building Blocks**

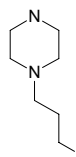
Brand	Catalog Number	MW
Aldrich	535044	142.16

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	571423	144.22

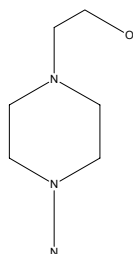
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	641642	144.22

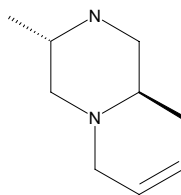


**Sigma - Aldrich Building Blocks**

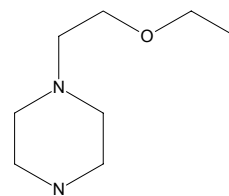
Brand	Catalog Number	MW
Aldrich	122963	145.21

**Sigma - Aldrich Building Blocks**

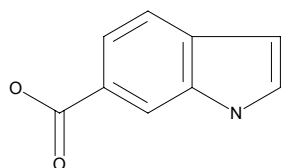
Brand	Catalog Number	MW
Aldrich	578118	154.26

**Sigma - Aldrich Building Blocks**

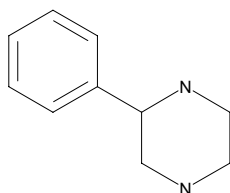
Brand	Catalog Number	MW
Aldrich	566861	158.25

**Sigma - Aldrich Building Blocks**

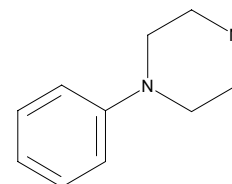
Brand	Catalog Number	MW
Aldrich	576662	161.16

**Sigma - Aldrich Building Blocks**

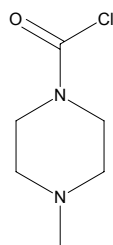
Brand	Catalog Number	MW
Aldrich	638641	162.24

**Sigma - Aldrich Building Blocks**

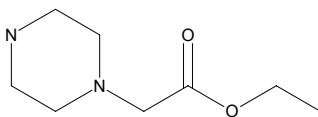
Brand	Catalog Number	MW
Aldrich	P30004	162.24

**Sigma - Aldrich Building Blocks**

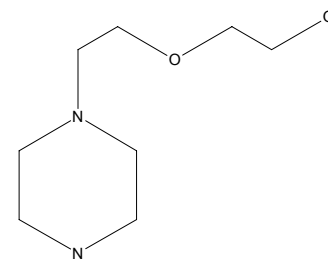
Brand	Catalog Number	MW
Aldrich	563250	162.62

**Sigma - Aldrich Building Blocks**

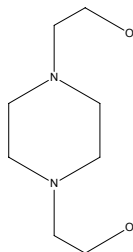
Brand	Catalog Number	MW
Aldrich	568732	172.23

**Sigma - Aldrich Building Blocks**

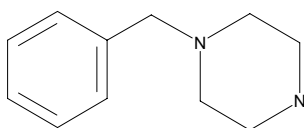
Brand	Catalog Number	MW
Aldrich	331260	174.24

**Sigma - Aldrich Building Blocks**

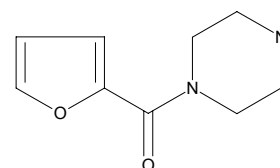
Brand	Catalog Number	MW
Aldrich	B45402	174.24

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	136832	176.26

**Sigma - Aldrich Building Blocks**

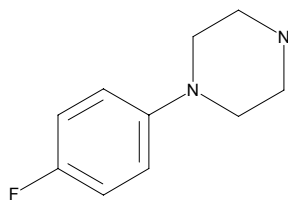
Brand	Catalog Number	MW
Aldrich	558966	180.21





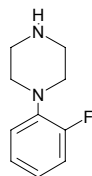
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	191337	180.23



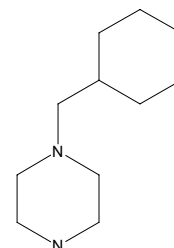
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	444804	180.23



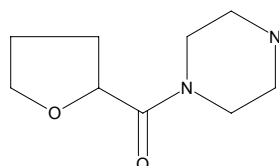
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	566535	182.31



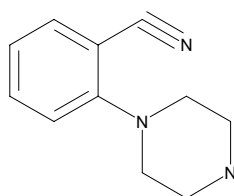
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	561304	184.24



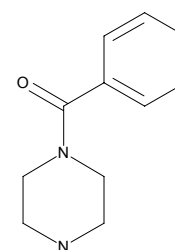
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	567183	187.25



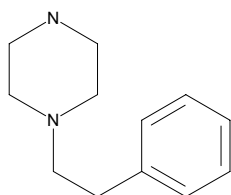
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	571083	190.25



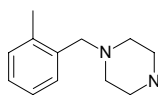
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	571458	190.29



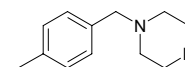
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	644226	190.29



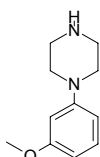
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	646172	190.29



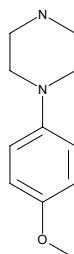
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	471682	192.26



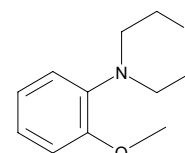
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	571415	192.26



Sigma - Aldrich Building Blocks

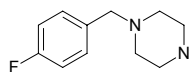
Brand	Catalog Number	MW
Aldrich	M22601	192.26





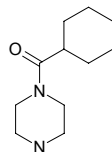
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	646164	194.25



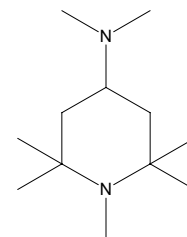
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	647128	196.29



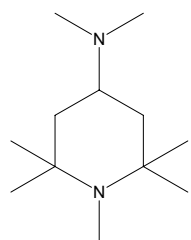
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	302201	198.35



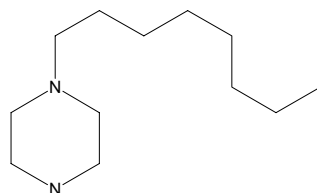
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	302201	198.35



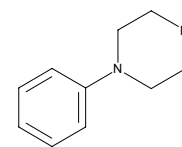
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	566896	198.35



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	281891	198.70

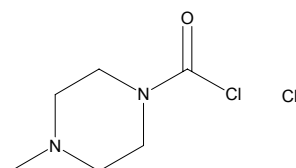


Cl



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	515027	199.08

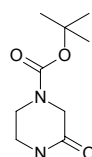


Cl



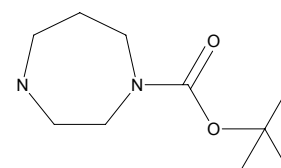
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	641057	200.24



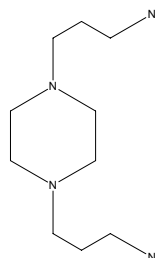
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	511382	200.28



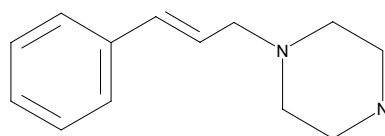
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	239488	200.33



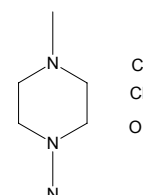
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	404470	202.30



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	A65131	206.12

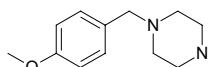


Cl
Cl
O



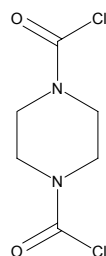
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	646156	206.29



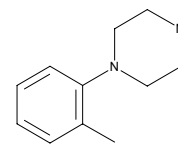
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	518530	211.05



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	T42005	212.72

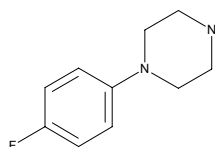


Cl



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	301280	216.69

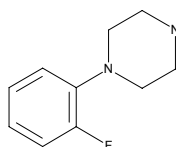


Cl



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	375306	216.69

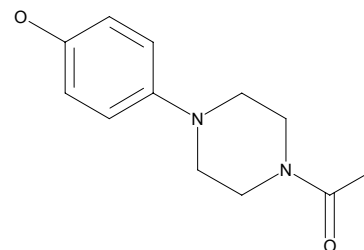


Cl



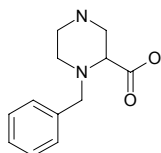
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	511943	220.27



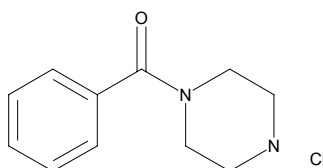
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	641987	220.27



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	592714	226.71

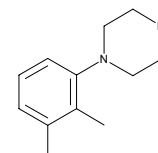


Cl



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	275182	226.75

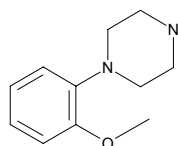


Cl



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	270067	228.72

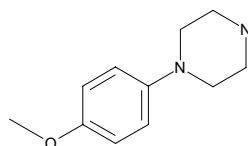


Cl



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	M23004	228.72

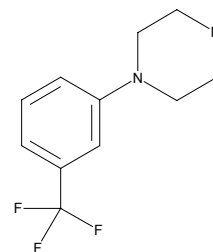


Cl



Sigma - Aldrich Building Blocks

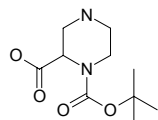
Brand	Catalog Number	MW
Aldrich	149853	230.23





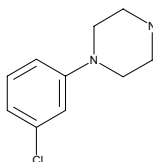
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	640131	230.27



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	125180	233.14

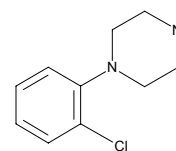


Cl



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	C67605	233.14

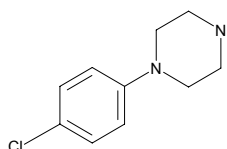


Cl



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	C68008	233.14

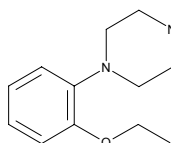


Cl



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	275174	242.75

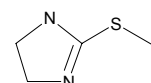


Cl



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	158844	244.10

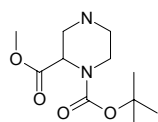


I



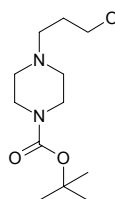
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	640158	244.29



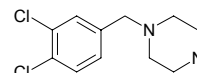
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	643467	244.34



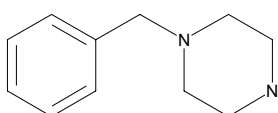
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	644269	245.15



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	539945	249.19

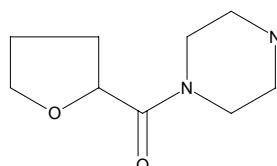


Cl
Cl



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	570990	265.15

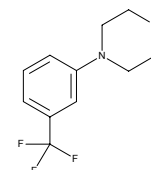


Br



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	270075	266.70

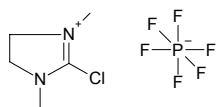


Cl



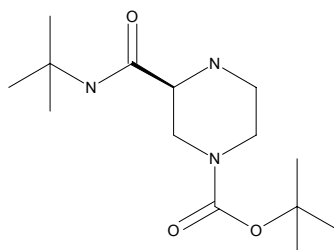
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	420336	278.57



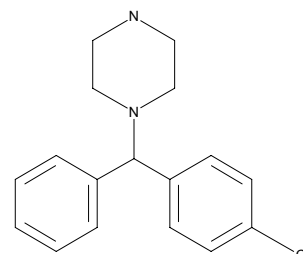
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	534382	285.39



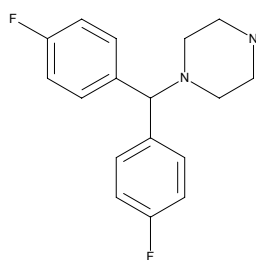
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	C24507	286.81



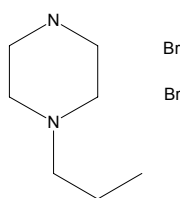
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	552402	288.34



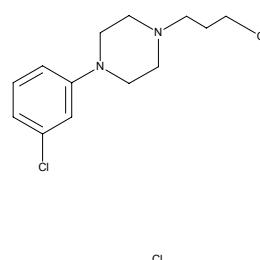
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	141674	290.04



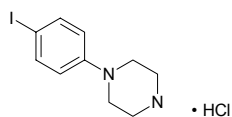
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	363464	309.67



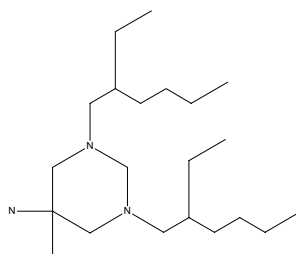
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	647268	324.59



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	259187	339.61



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	473588	452.51

