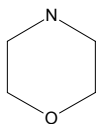




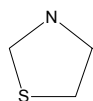
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	394467	87.12



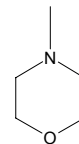
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	149691	89.16



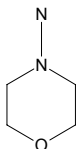
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	407704	101.15



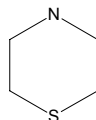
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	A66308	102.14



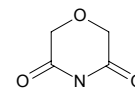
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	196274	103.19



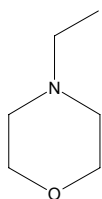
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	493724	115.09



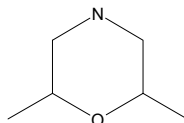
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	109932	115.18



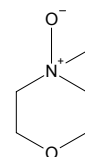
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	126527	115.18



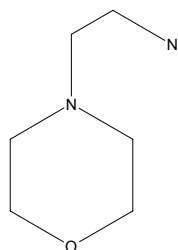
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	224286	117.15



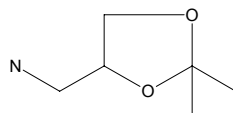
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	A55004	130.19



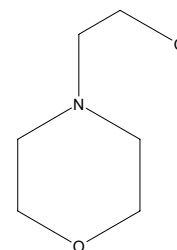
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	483117	131.18



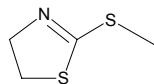
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	H28203	131.18

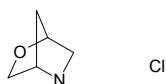


**Sigma - Aldrich Building Blocks**

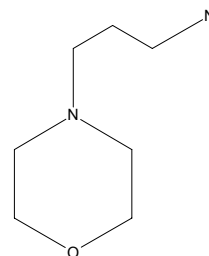
Brand	Catalog Number	MW
Aldrich	250406	133.24

**Sigma - Aldrich Building Blocks**

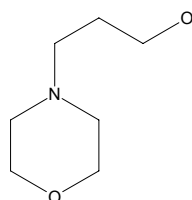
Brand	Catalog Number	MW
Aldrich	473022	135.59

**Sigma - Aldrich Building Blocks**

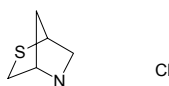
Brand	Catalog Number	MW
Aldrich	123099	144.22

**Sigma - Aldrich Building Blocks**

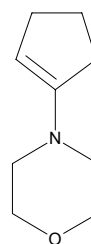
Brand	Catalog Number	MW
Aldrich	567280	145.20

**Sigma - Aldrich Building Blocks**

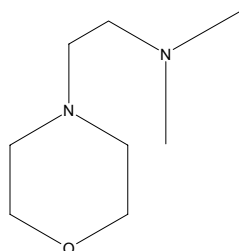
Brand	Catalog Number	MW
Aldrich	473014	151.66

**Sigma - Aldrich Building Blocks**

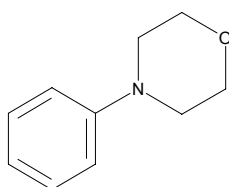
Brand	Catalog Number	MW
Aldrich	C114901	153.23

**Sigma - Aldrich Building Blocks**

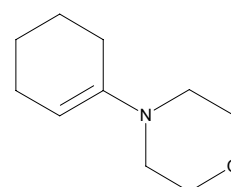
Brand	Catalog Number	MW
Aldrich	278432	158.25

**Sigma - Aldrich Building Blocks**

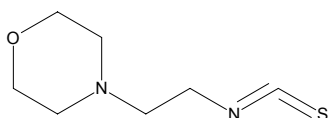
Brand	Catalog Number	MW
Aldrich	211338	163.22

**Sigma - Aldrich Building Blocks**

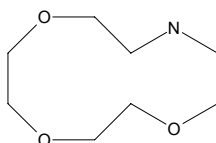
Brand	Catalog Number	MW
Aldrich	M87800	167.25

**Sigma - Aldrich Building Blocks**

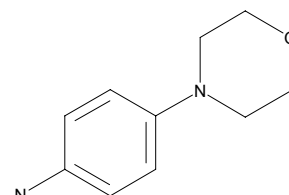
Brand	Catalog Number	MW
Aldrich	560472	172.25

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	364088	175.23

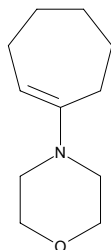
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	197157	178.24

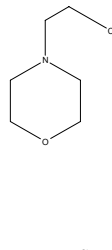


**Sigma - Aldrich Building Blocks**

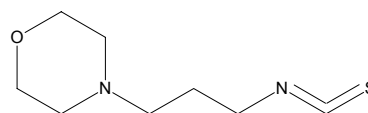
Brand	Catalog Number	MW
Aldrich	385573	181.28

**Sigma - Aldrich Building Blocks**

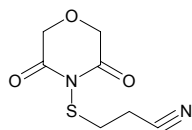
Brand	Catalog Number	MW
Aldrich	C42203	186.08

**Sigma - Aldrich Building Blocks**

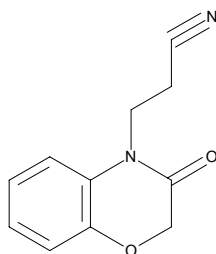
Brand	Catalog Number	MW
Aldrich	560464	186.28

**Sigma - Aldrich Building Blocks**

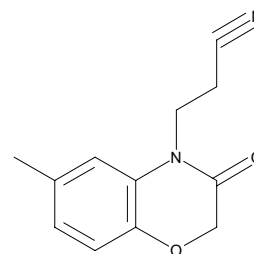
Brand	Catalog Number	MW
Aldrich	493635	200.22

**Sigma - Aldrich Building Blocks**

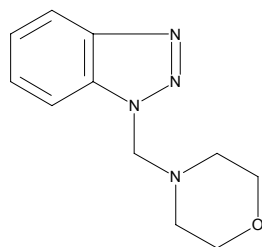
Brand	Catalog Number	MW
Aldrich	535621	202.21

**Sigma - Aldrich Building Blocks**

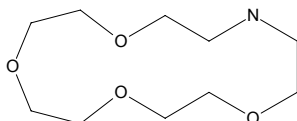
Brand	Catalog Number	MW
Aldrich	548901	216.24

**Sigma - Aldrich Building Blocks**

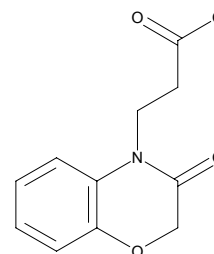
Brand	Catalog Number	MW
Aldrich	467502	218.26

**Sigma - Aldrich Building Blocks**

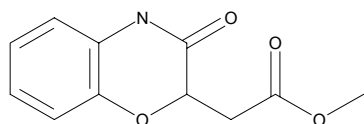
Brand	Catalog Number	MW
Aldrich	364096	219.28

**Sigma - Aldrich Building Blocks**

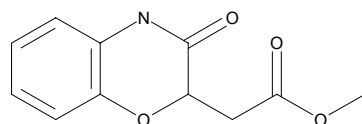
Brand	Catalog Number	MW
Aldrich	548898	221.21

**Sigma - Aldrich Building Blocks**

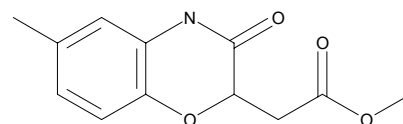
Brand	Catalog Number	MW
Aldrich	550094	221.21

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	550094	221.21

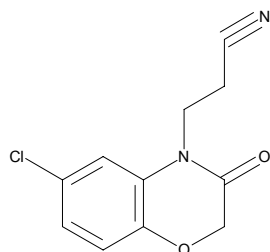
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	550108	235.24

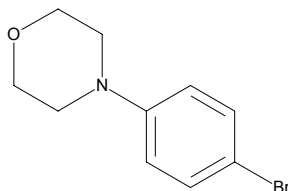


**Sigma - Aldrich Building Blocks**

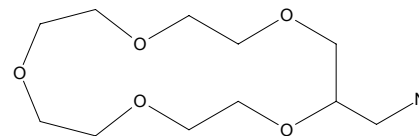
Brand	Catalog Number	MW
Aldrich	548928	236.66

**Sigma - Aldrich Building Blocks**

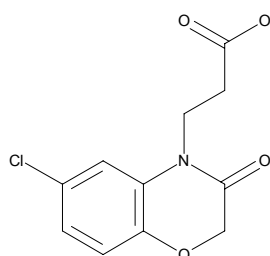
Brand	Catalog Number	MW
Aldrich	591777	242.12

**Sigma - Aldrich Building Blocks**

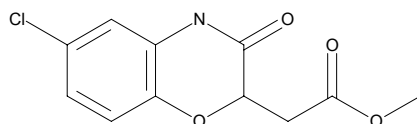
Brand	Catalog Number	MW
Aldrich	388416	249.31

**Sigma - Aldrich Building Blocks**

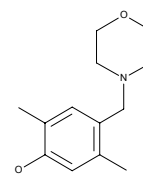
Brand	Catalog Number	MW
Aldrich	535613	255.66

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	550116	255.66

**Sigma - Aldrich Building Blocks**

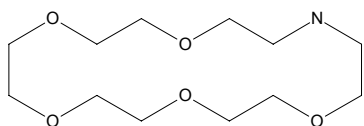
Brand	Catalog Number	MW
Aldrich	186716	257.76



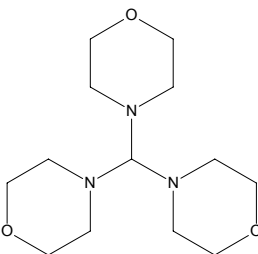
Cl

**Sigma - Aldrich Building Blocks**

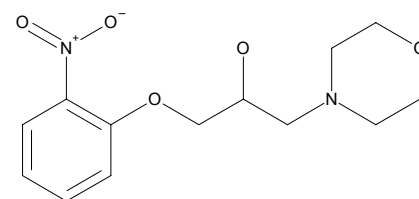
Brand	Catalog Number	MW
Aldrich	364118	263.34

**Sigma - Aldrich Building Blocks**

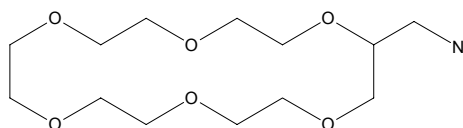
Brand	Catalog Number	MW
Aldrich	513288	271.36

**Sigma - Aldrich Building Blocks**

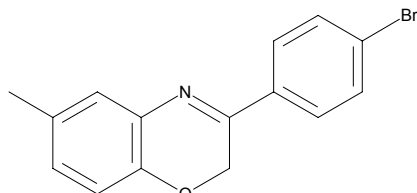
Brand	Catalog Number	MW
Aldrich	630500	282.30

**Sigma - Aldrich Building Blocks**

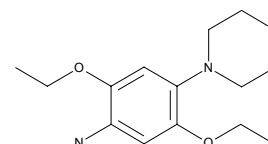
Brand	Catalog Number	MW
Aldrich	388432	293.36

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	551023	302.17

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	439363	302.80



Cl



Sigma - Aldrich
Building Blocks

Brand	Catalog Number	MW
Aldrich	551031	322.59

