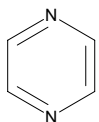




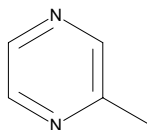
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	P56003	80.09



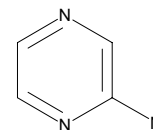
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	M75608	94.12



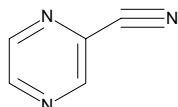
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	A76958	95.10



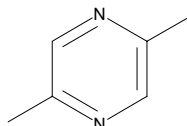
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	349496	105.10



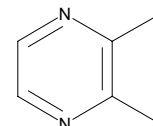
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	175420	108.14



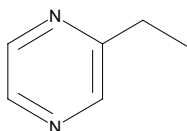
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	199400	108.14



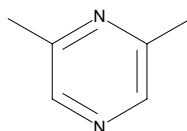
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	250384	108.14



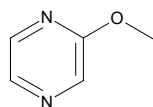
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	D181803	108.14



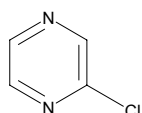
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	291420	110.12



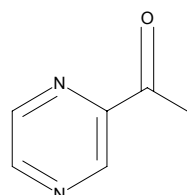
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	132489	114.53



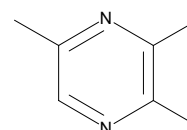
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	251801	122.13



**Sigma - Aldrich Building Blocks**

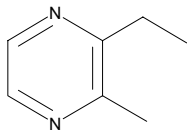
Brand	Catalog Number	MW
Aldrich	199419	122.17





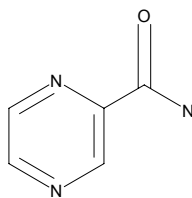
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	289582	122.17



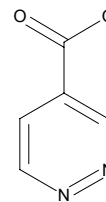
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	131571	123.12



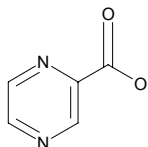
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	297763	124.10



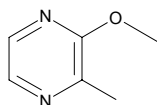
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	P56100	124.10



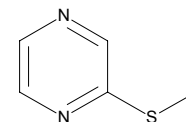
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	297941	124.14



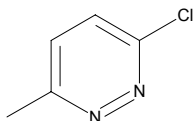
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	467677	126.18



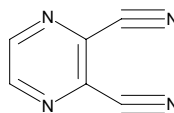
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	637637	128.56



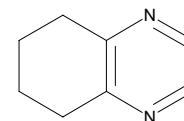
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	399620	130.11



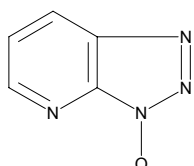
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	272701	134.18



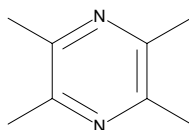
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	559156	136.11



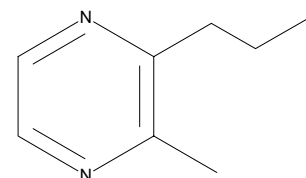
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	183938	136.20



**Sigma - Aldrich Building Blocks**

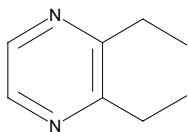
Brand	Catalog Number	MW
Aldrich	278025	136.20





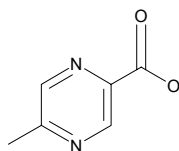
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	292982	136.20



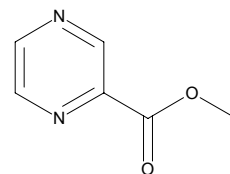
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	347647	138.13



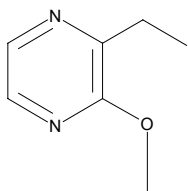
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	519677	138.13



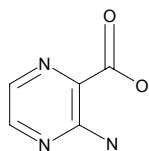
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	293806	138.17



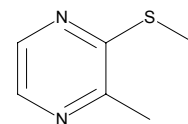
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	A76982	139.11



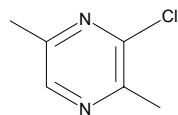
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	545791	140.21



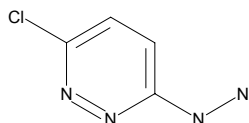
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	C38400	142.59



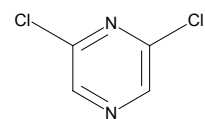
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	632619	144.56



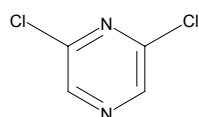
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	132497	148.98



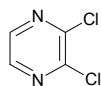
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	373575	148.98



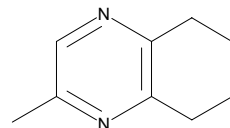
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	465208	148.98



**Sigma - Aldrich Building Blocks**

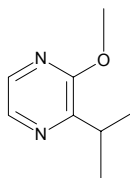
Brand	Catalog Number	MW
Aldrich	293814	150.23





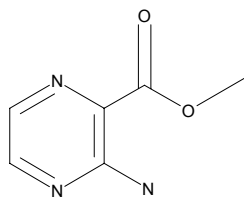
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	243132	152.20



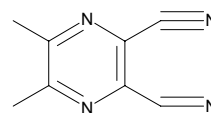
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	276154	153.14



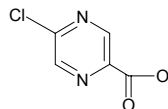
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	399639	158.16



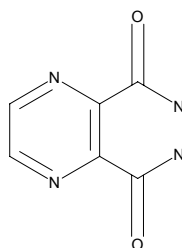
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	439576	158.54



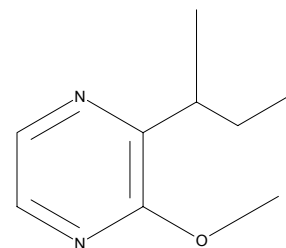
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	P56151	166.14



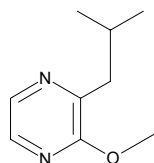
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	243116	166.22



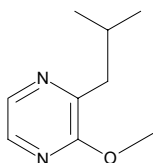
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	243124	166.22



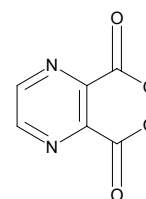
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	297666	166.22



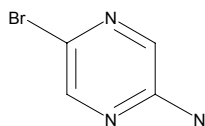
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	P56208	168.11



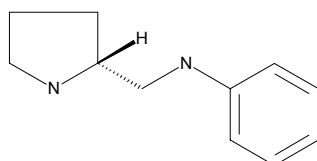
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	636320	174.00



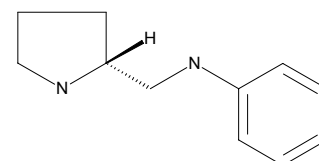
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	510505	176.26



**Sigma - Aldrich Building Blocks**

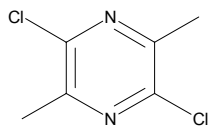
Brand	Catalog Number	MW
Aldrich	510505	176.26





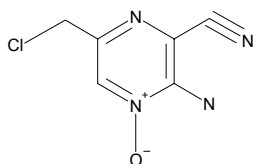
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	630055	177.03



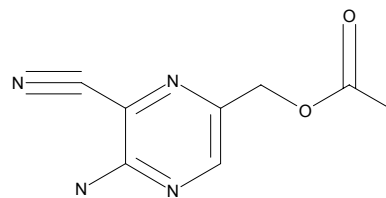
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	221139	184.59



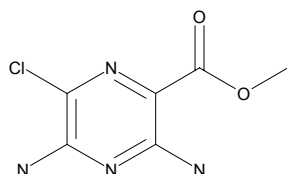
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	861626	192.18



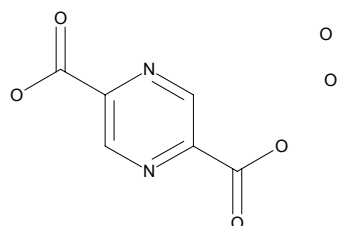
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	596833	202.60



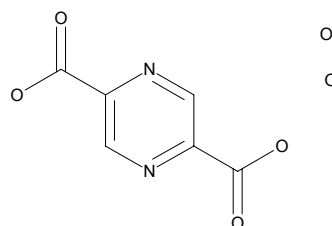
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	540498	204.14



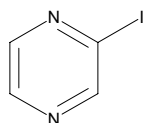
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	540498	204.14



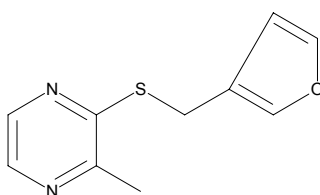
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	513164	205.99



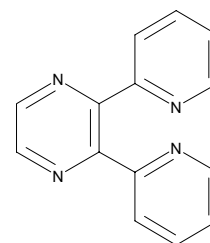
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	545783	206.27



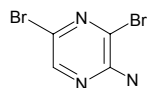
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	281646	234.26



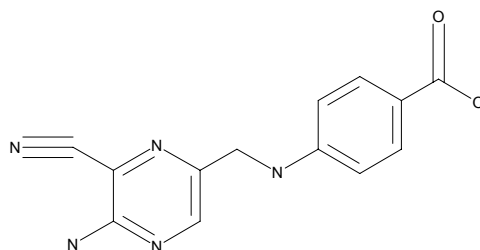
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	494364	252.90



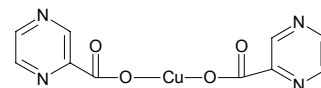
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	861596	269.26



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	300993	309.73





**Sigma - Aldrich  
Building Blocks**

Brand	Catalog Number	MW
Aldrich	495557	388.44

