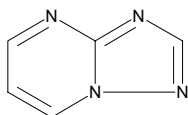




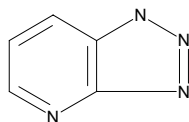
**Sigma - Aldrich  
Building Blocks**

Brand	Catalog Number	MW
Aldrich	324167	120.11



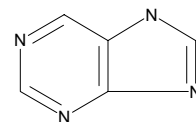
**Sigma - Aldrich  
Building Blocks**

Brand	Catalog Number	MW
Aldrich	361011	120.11



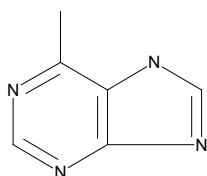
**Sigma - Aldrich  
Building Blocks**

Brand	Catalog Number	MW
Aldrich	P55805	120.11



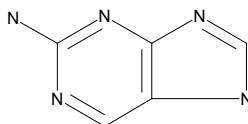
**Sigma - Aldrich  
Building Blocks**

Brand	Catalog Number	MW
Aldrich	852740	134.14



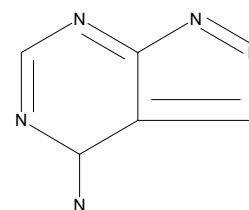
**Sigma - Aldrich  
Building Blocks**

Brand	Catalog Number	MW
Aldrich	285080	135.13



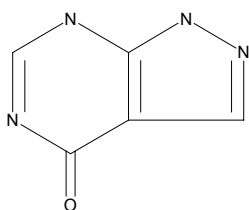
**Sigma - Aldrich  
Building Blocks**

Brand	Catalog Number	MW
Aldrich	A77806	135.13



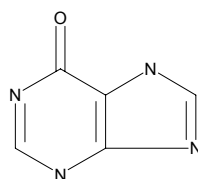
**Sigma - Aldrich  
Building Blocks**

Brand	Catalog Number	MW
Aldrich	H56606	136.11



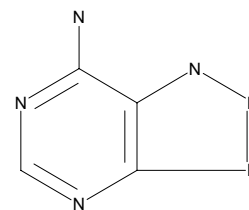
**Sigma - Aldrich  
Building Blocks**

Brand	Catalog Number	MW
Aldrich	H61200	136.11



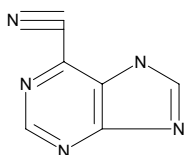
**Sigma - Aldrich  
Building Blocks**

Brand	Catalog Number	MW
Aldrich	855529	136.12



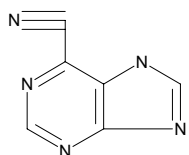
**Sigma - Aldrich  
Building Blocks**

Brand	Catalog Number	MW
Aldrich	852503	145.12



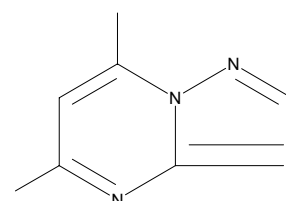
**Sigma - Aldrich  
Building Blocks**

Brand	Catalog Number	MW
Aldrich	852503	145.12



**Sigma - Aldrich  
Building Blocks**

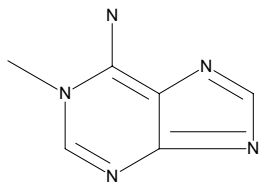
Brand	Catalog Number	MW
Aldrich	D189790	148.17





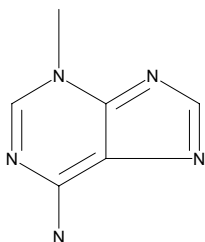
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	215325	149.16



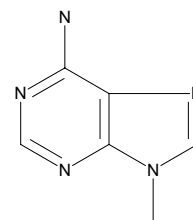
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	280879	149.16



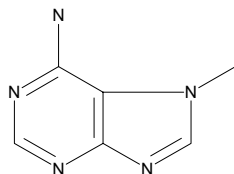
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	547328	149.16



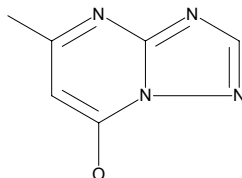
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	547336	149.16



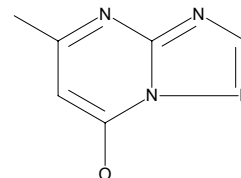
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	177679	150.14



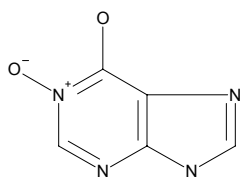
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	422509	150.14



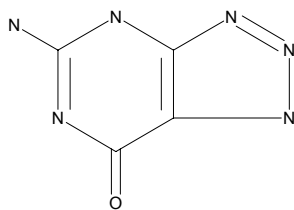
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	557889	152.11



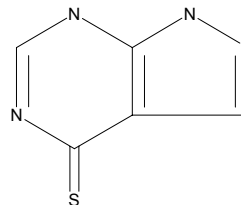
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	861448	152.12



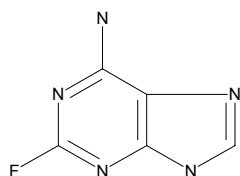
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	153060	152.18



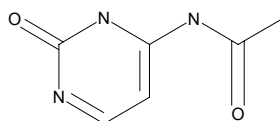
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	535087	153.12



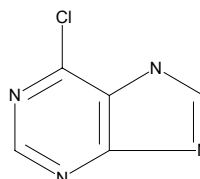
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	377910	153.14



**Sigma - Aldrich Building Blocks**

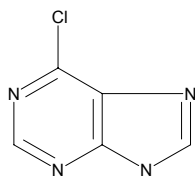
Brand	Catalog Number	MW
Aldrich	161179	154.56





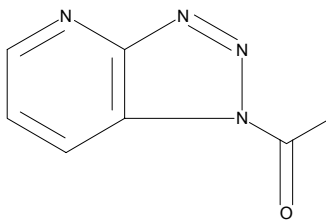
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	511617	154.56



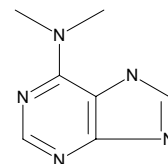
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	347272	162.15



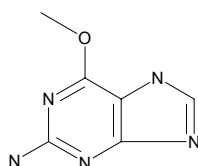
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	852554	163.18



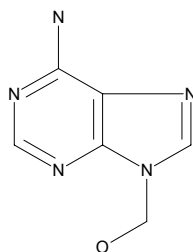
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	363057	165.16



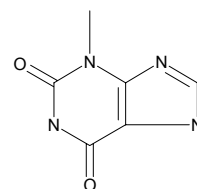
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	559385	165.16



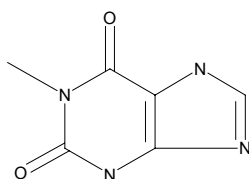
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	222526	166.14



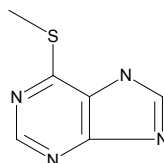
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	280984	166.14



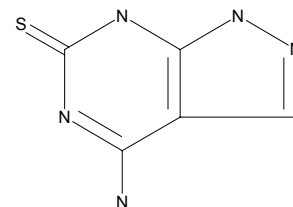
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	852732	166.21



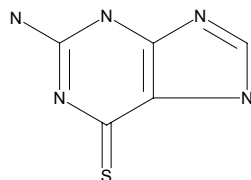
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	A60202	167.19



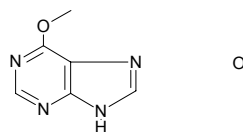
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	A76907	167.19



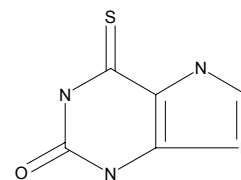
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	852708	168.16



**Sigma - Aldrich Building Blocks**

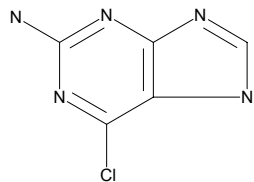
Brand	Catalog Number	MW
Aldrich	852570	168.18





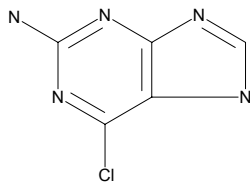
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	109789	169.57



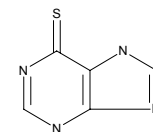
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	342300	169.57



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	852678	170.19

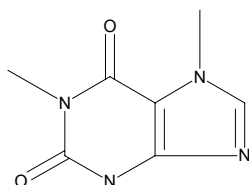


o



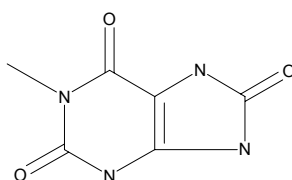
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	271519	180.17



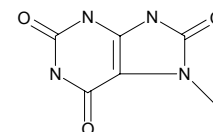
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	360236	182.14



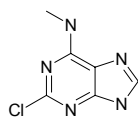
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	360244	182.14



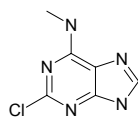
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	644595	183.60



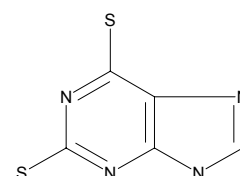
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	644595	183.60



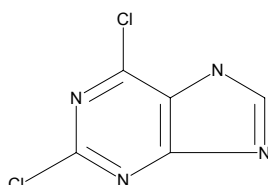
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	852627	184.24



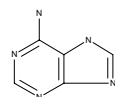
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	D73103	189.00



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	271934	189.61



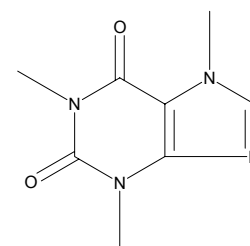
ci

o



**Sigma - Aldrich Building Blocks**

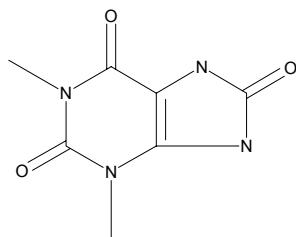
Brand	Catalog Number	MW
Aldrich	C53	194.19





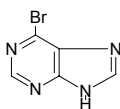
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	245097	196.17



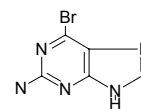
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	104981	199.01



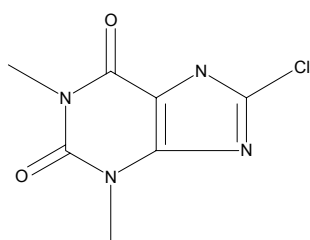
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	475254	214.03



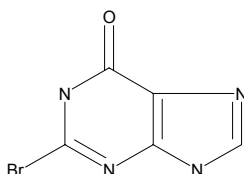
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	C71807	214.61



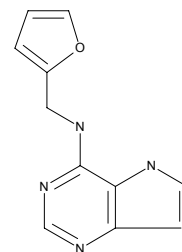
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	535125	215.01



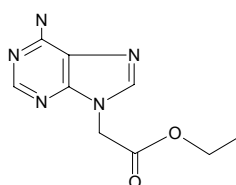
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	852643	215.22



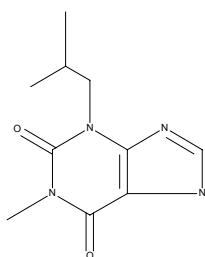
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	461768	221.22



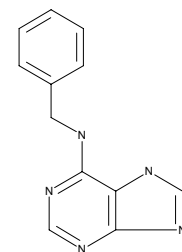
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	858455	222.25



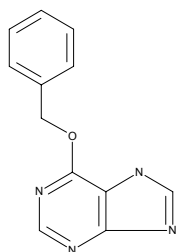
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	852430	225.25



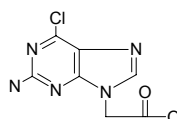
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	387606	226.24



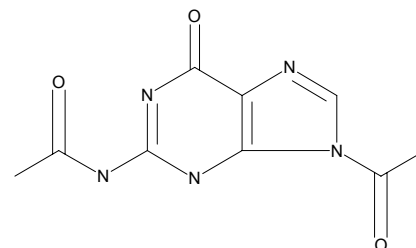
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	478563	227.61



**Sigma - Aldrich Building Blocks**

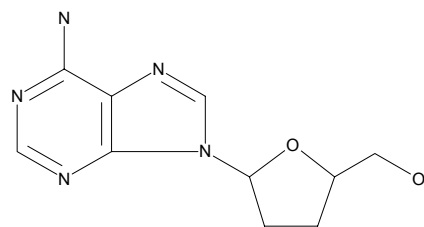
Brand	Catalog Number	MW
Aldrich	377899	235.20





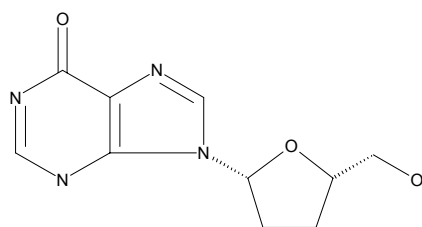
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	308331	235.25



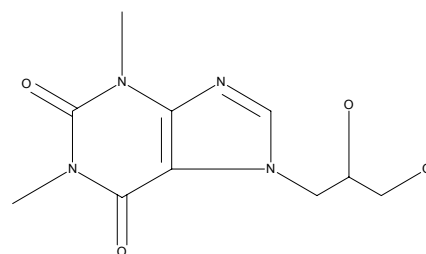
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	362646	236.23



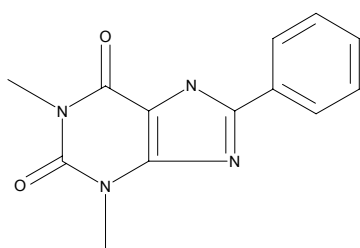
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	D119407	254.25



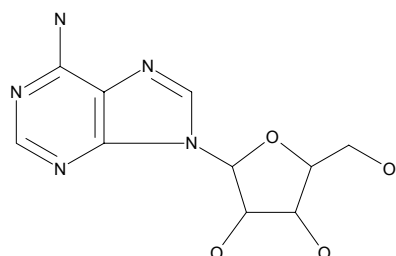
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	227161	256.27



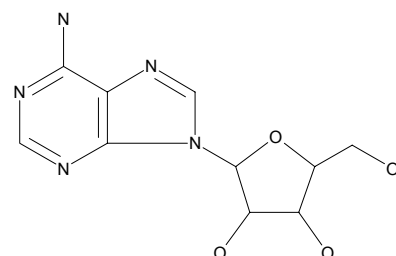
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	146595	267.25



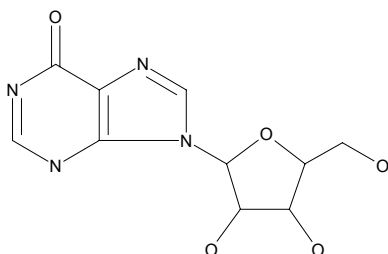
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	284971	267.25



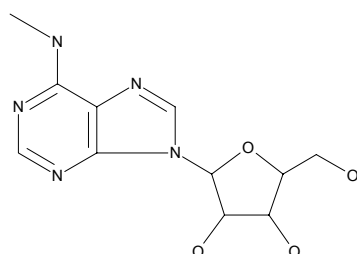
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	I6407	268.23



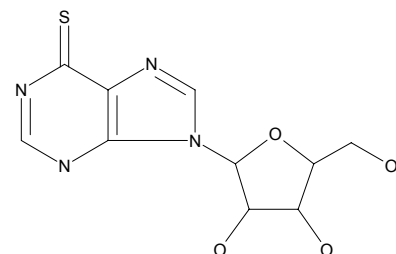
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	852724	281.27



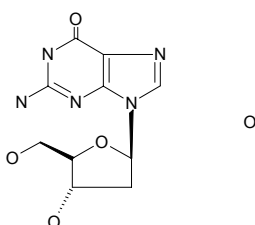
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	852686	284.30



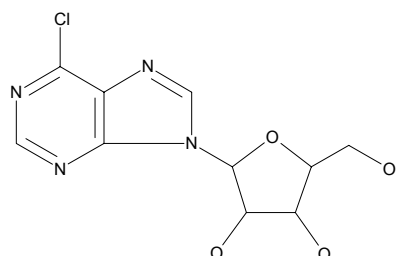
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	854999	285.26



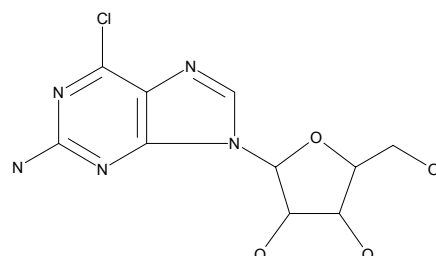
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	852481	286.68



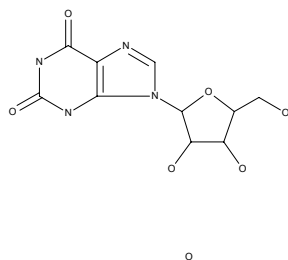
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	852392	301.69

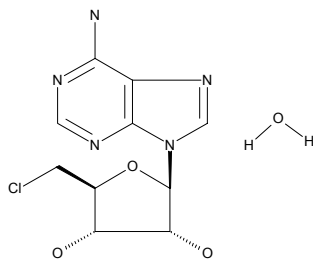


**Sigma - Aldrich Building Blocks**

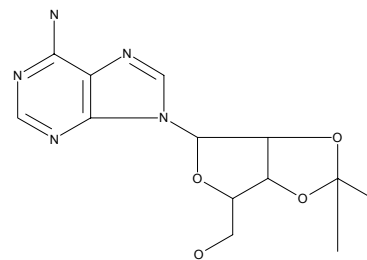
Brand	Catalog Number	MW
Aldrich	223344	302.25

**Sigma - Aldrich Building Blocks**

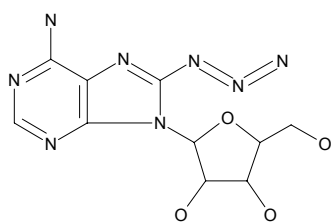
Brand	Catalog Number	MW
Aldrich	548488	303.71

**Sigma - Aldrich Building Blocks**

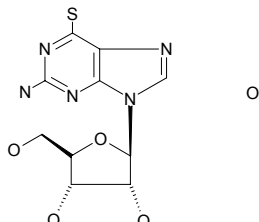
Brand	Catalog Number	MW
Aldrich	122404	307.31

**Sigma - Aldrich Building Blocks**

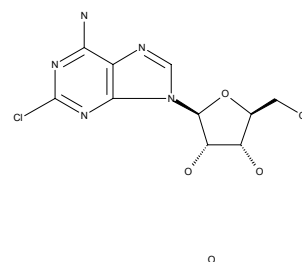
Brand	Catalog Number	MW
Aldrich	263788	310.27

**Sigma - Aldrich Building Blocks**

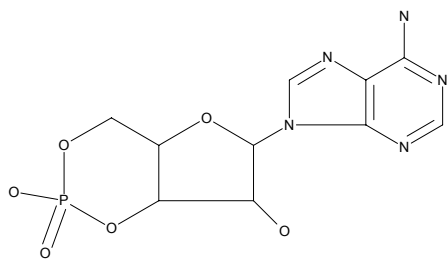
Brand	Catalog Number	MW
Aldrich	858412	317.33

**Sigma - Aldrich Building Blocks**

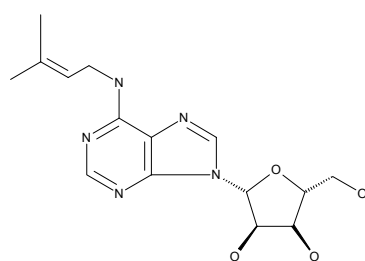
Brand	Catalog Number	MW
Aldrich	861863	319.71

**Sigma - Aldrich Building Blocks**

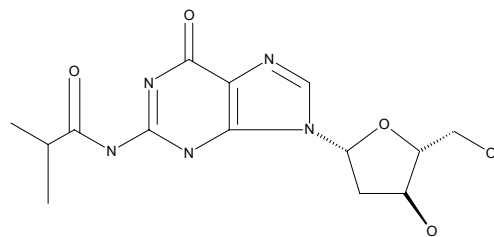
Brand	Catalog Number	MW
Aldrich	851205	329.21

**Sigma - Aldrich Building Blocks**

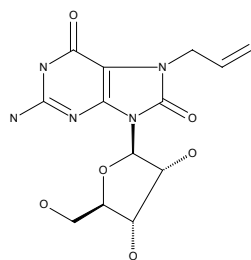
Brand	Catalog Number	MW
Aldrich	852619	335.37

**Sigma - Aldrich Building Blocks**

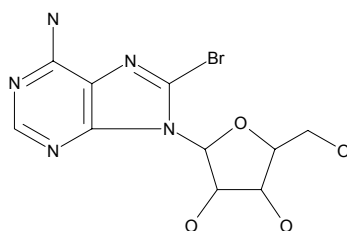
Brand	Catalog Number	MW
Aldrich	258482	337.34

**Sigma - Aldrich Building Blocks**

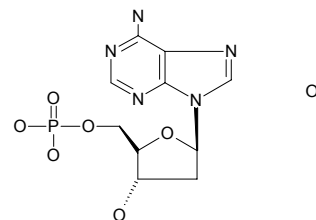
Brand	Catalog Number	MW
Aldrich	496812	339.31

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	127507	346.14

**Sigma - Aldrich Building Blocks**

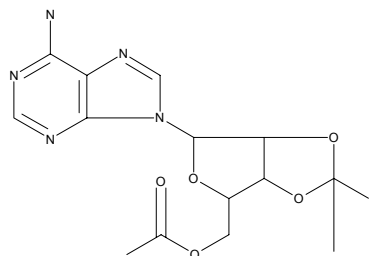
Brand	Catalog Number	MW
Aldrich	852198	349.24





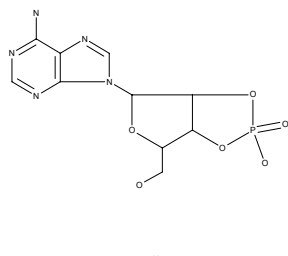
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	858013	349.35



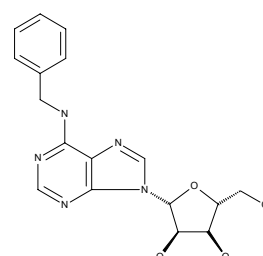
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	857920	353.21



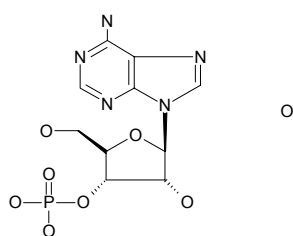
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	852449	357.37



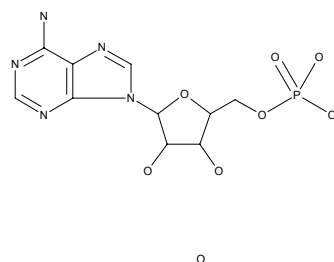
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	851949	365.24



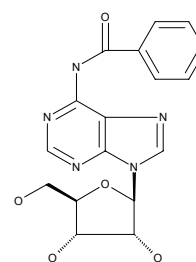
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	A25008	365.24



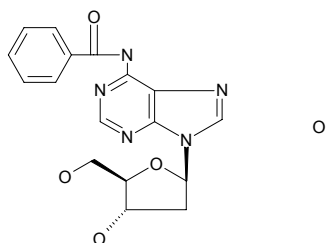
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	586846	371.36



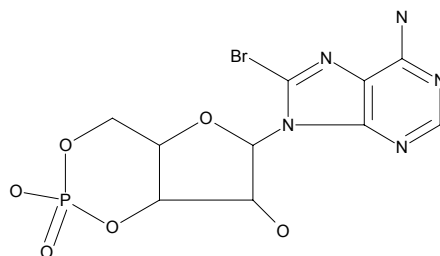
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	258466	373.37



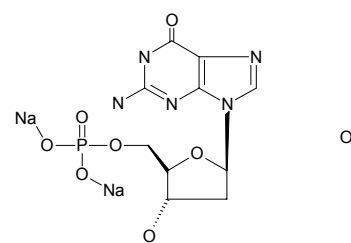
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	372447	408.11



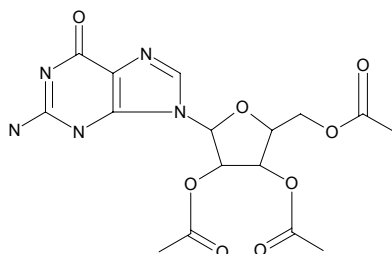
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	852228	409.21



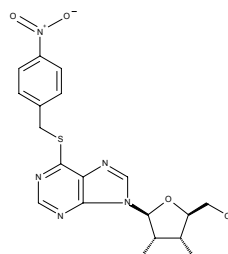
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	850926	409.36



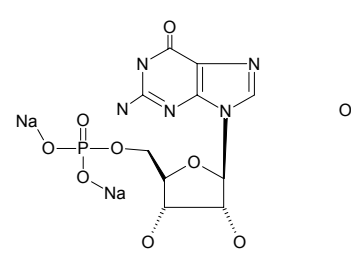
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	861499	419.42



### Sigma - Aldrich Building Blocks

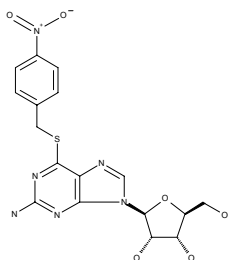
Brand	Catalog Number	MW
Aldrich	852856	425.20





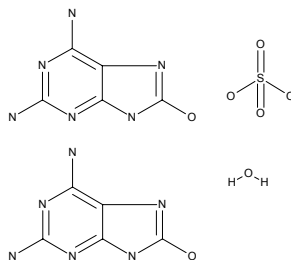
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	861669	434.43



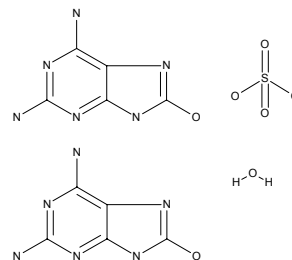
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	530662	448.38



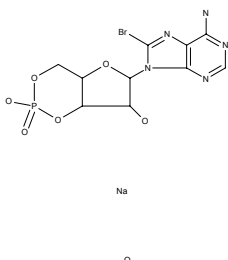
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	530662	448.38



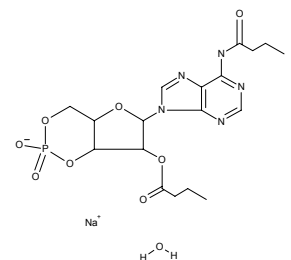
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	858463	450.12



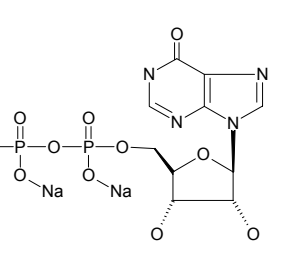
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	851965	509.39



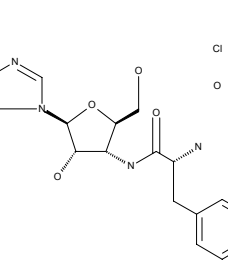
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	852074	512.15



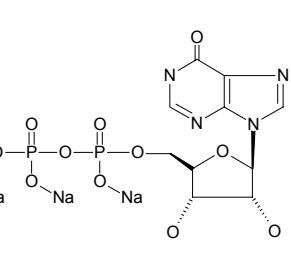
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	856339	526.00



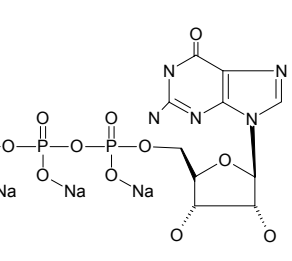
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	367184	592.13



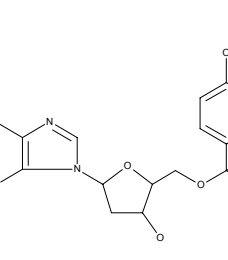
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	852058	607.15



### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	258512	639.71



### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	258490	657.73

