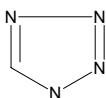




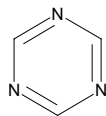
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	554049	70.05



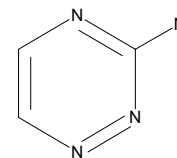
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	T46051	81.08



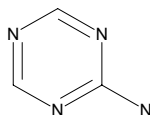
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	100625	96.09



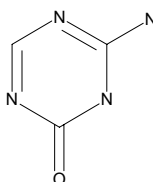
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	533823	96.09



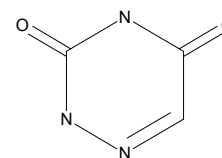
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	855049	112.09



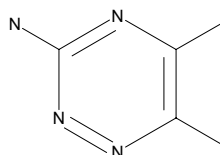
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	123293	113.08



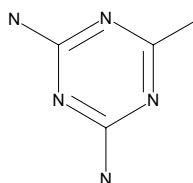
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	133469	124.15



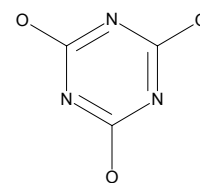
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	407801	125.13



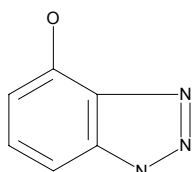
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	C95455	129.08



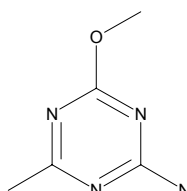
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	632791	135.13



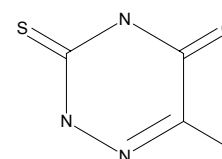
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	522996	140.15



Sigma - Aldrich Building Blocks

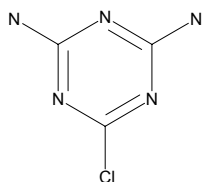
Brand	Catalog Number	MW
Aldrich	275514	143.17





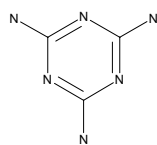
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	C33301	145.55



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	337803	146.13

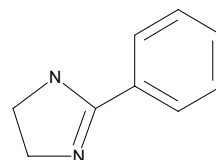


F



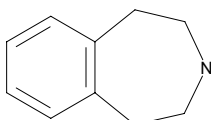
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	213039	146.19



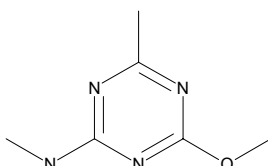
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	562483	147.22



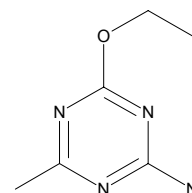
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	549843	154.17



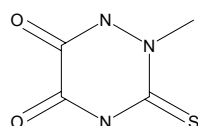
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	549851	154.17



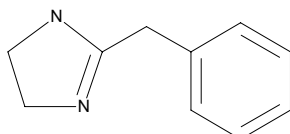
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	549746	159.17



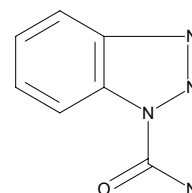
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	293490	160.22



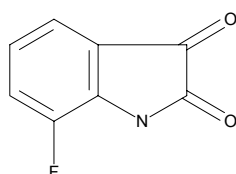
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	596485	162.15



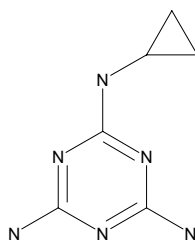
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	633194	165.12



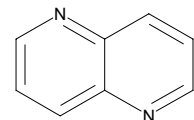
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	551295	166.19



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	594164	166.61

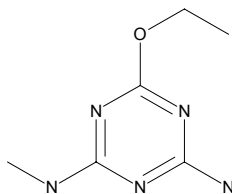


H-Cl



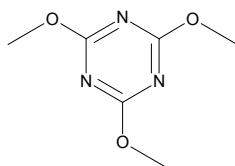
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	523003	169.19



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	407879	171.16



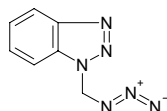
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	J1001	173.26



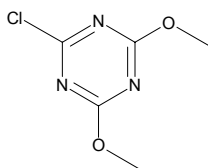
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	642576	174.17



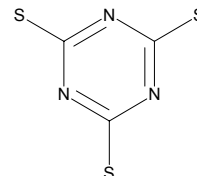
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	375217	175.58



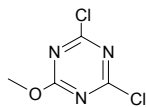
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	T88595	177.27



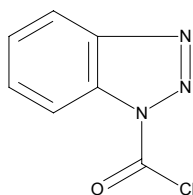
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	213535	179.99



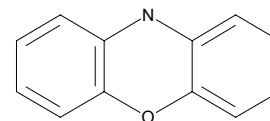
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	633917	181.58



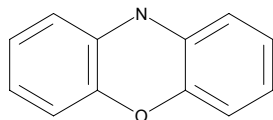
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	263893	183.21



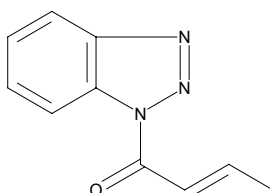
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	P14858	183.21



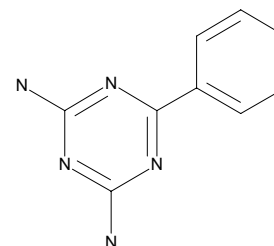
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	596361	187.20



Sigma - Aldrich Building Blocks

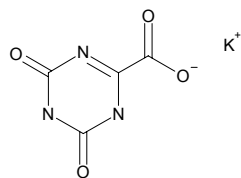
Brand	Catalog Number	MW
Aldrich	D23408	187.21





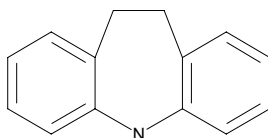
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	156124	195.18



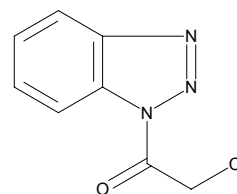
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	I1308	195.27



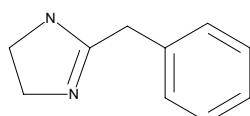
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	596922	195.61



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	T35467	196.68

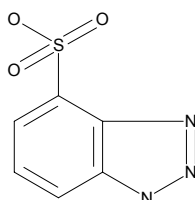


Cl



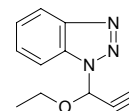
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	632805	199.19



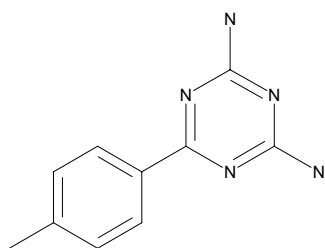
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	642584	201.23



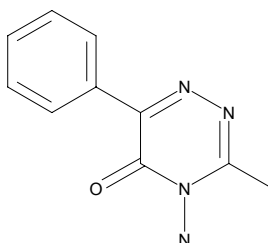
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	579521	201.23



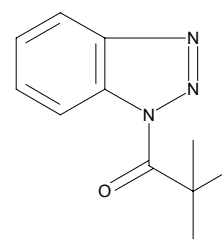
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	559806	202.22



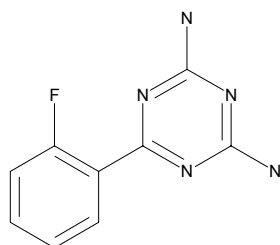
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	596264	203.25



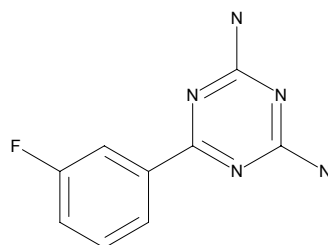
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	597341	205.20



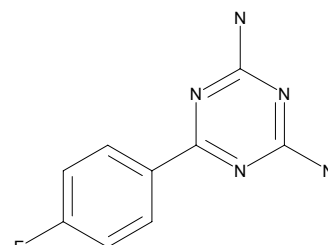
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	597457	205.20



Sigma - Aldrich Building Blocks

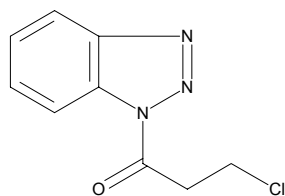
Brand	Catalog Number	MW
Aldrich	597570	205.20





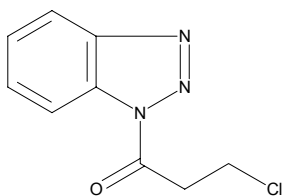
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	596817	209.64



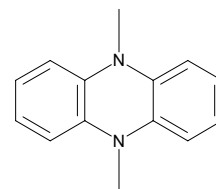
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	596817	209.64



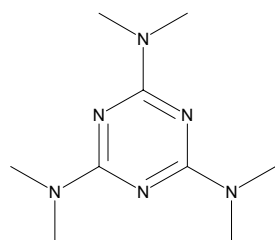
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	395013	210.28



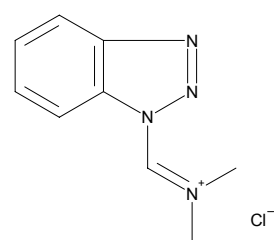
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	549835	210.28



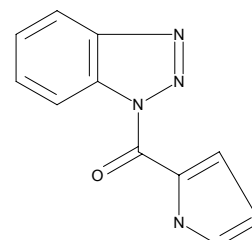
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	633763	210.67



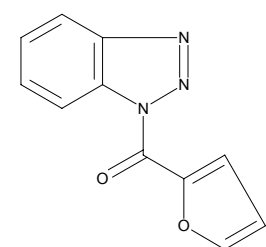
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	631892	212.21



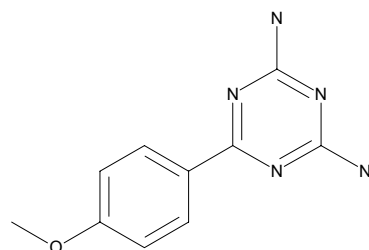
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	630799	213.20



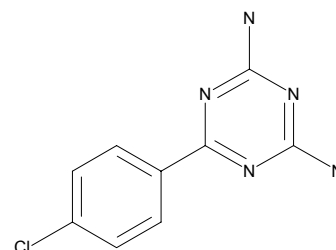
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	579939	217.23



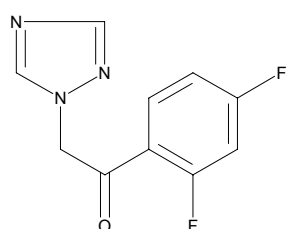
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	592188	221.65



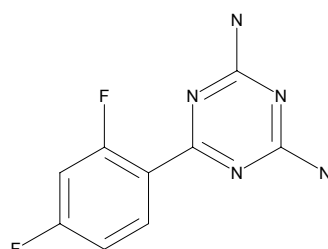
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	636886	223.18



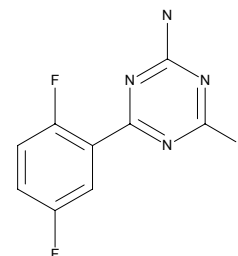
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	597686	223.19



Sigma - Aldrich Building Blocks

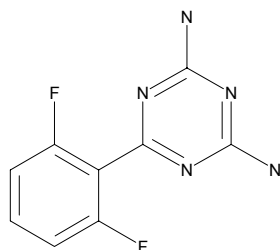
Brand	Catalog Number	MW
Aldrich	597783	223.19





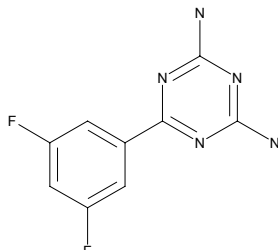
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	597899	223.19



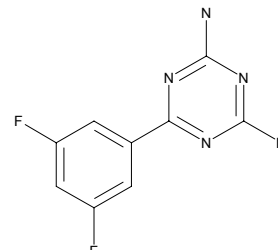
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	598003	223.19



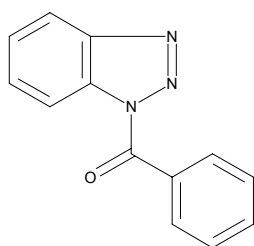
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	598003	223.19



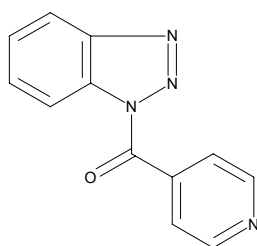
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	596590	223.24



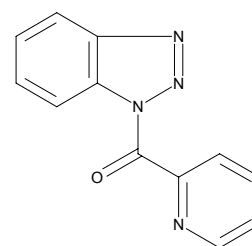
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	631906	224.22



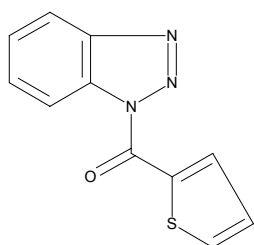
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	631914	224.22



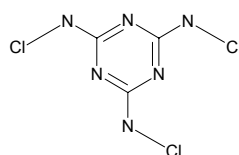
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	630772	229.26



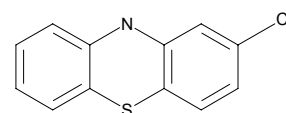
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	106240	229.46



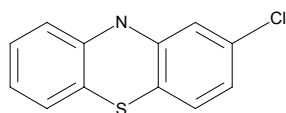
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	373559	233.72



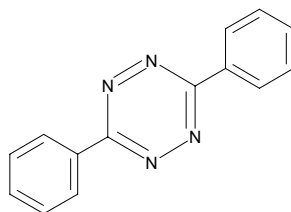
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	C63006	233.72



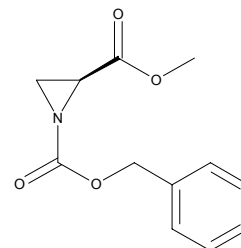
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	403555	234.26



Sigma - Aldrich Building Blocks

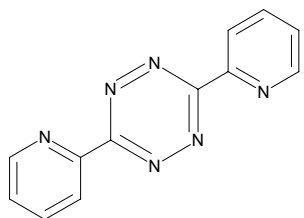
Brand	Catalog Number	MW
Aldrich	516031	235.24





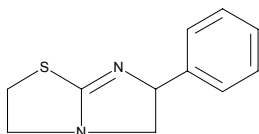
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	403547	236.24



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	196134	240.76

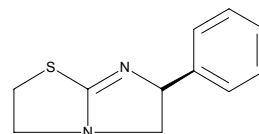


Cl



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	196142	240.76

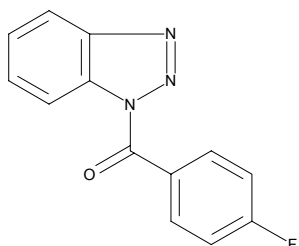


Cl



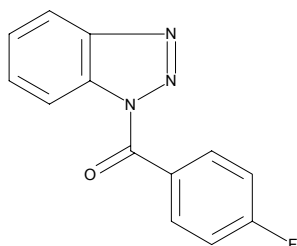
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	630780	241.23



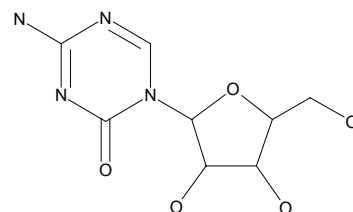
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	630780	241.23



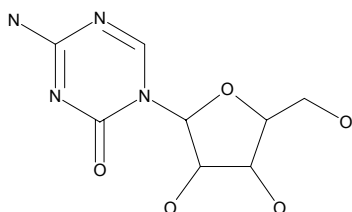
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	852880	244.21



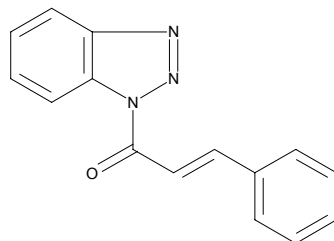
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	852880	244.21



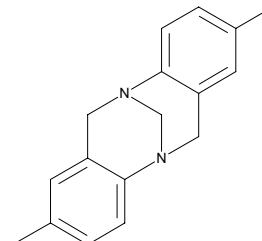
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	630616	249.27



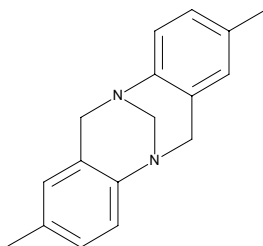
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	132225	250.35



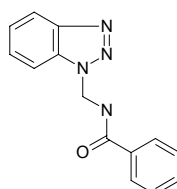
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	374644	250.35



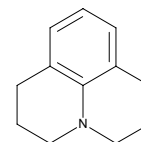
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	642851	252.28



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	407518	254.17

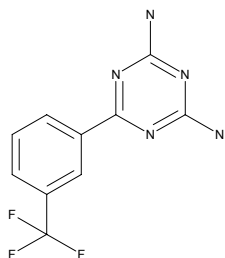


Br



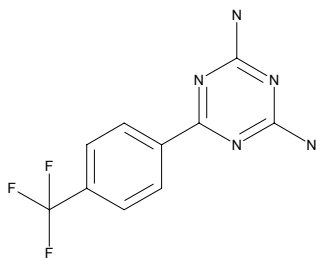
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	579580	255.20



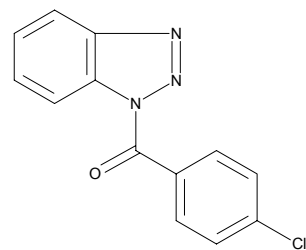
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	579807	255.20



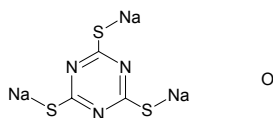
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	596701	257.68



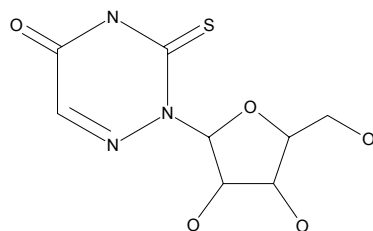
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	382922	261.23



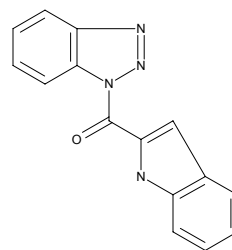
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	209406	261.26



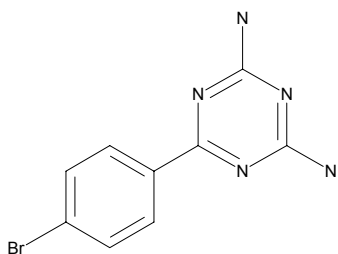
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	630608	262.27



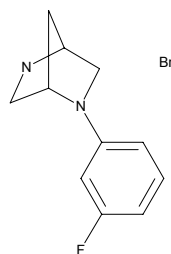
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	597244	266.10



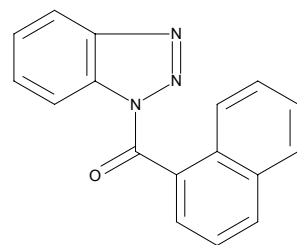
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	473405	273.15



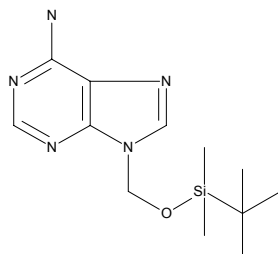
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	632163	273.30



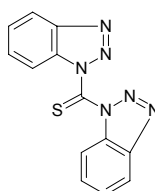
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	637734	279.42



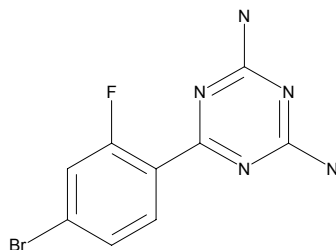
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	643610	280.31



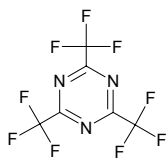
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	598119	284.09

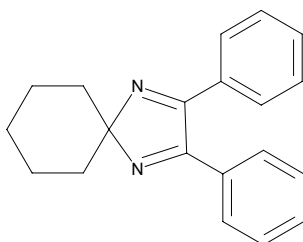


**Sigma - Aldrich Building Blocks**

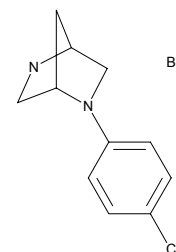
Brand	Catalog Number	MW
Aldrich	443980	285.07

**Sigma - Aldrich Building Blocks**

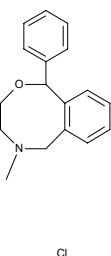
Brand	Catalog Number	MW
Aldrich	371467	288.40

**Sigma - Aldrich Building Blocks**

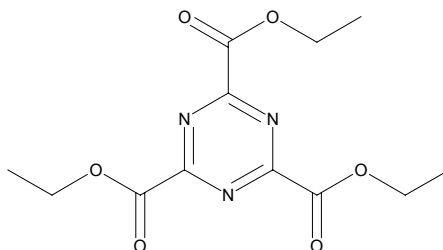
Brand	Catalog Number	MW
Aldrich	473391	289.60

**Sigma - Aldrich Building Blocks**

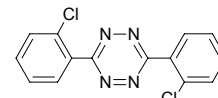
Brand	Catalog Number	MW
Aldrich	257745	289.81

**Sigma - Aldrich Building Blocks**

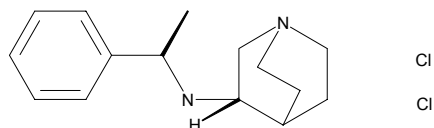
Brand	Catalog Number	MW
Aldrich	576697	297.27

**Sigma - Aldrich Building Blocks**

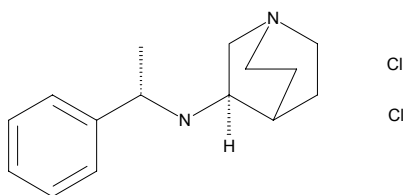
Brand	Catalog Number	MW
Aldrich	476838	303.15

**Sigma - Aldrich Building Blocks**

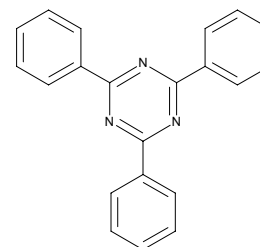
Brand	Catalog Number	MW
Aldrich	487244	303.28

**Sigma - Aldrich Building Blocks**

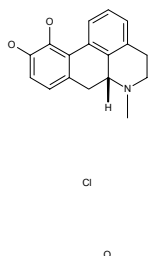
Brand	Catalog Number	MW
Aldrich	510343	303.28

**Sigma - Aldrich Building Blocks**

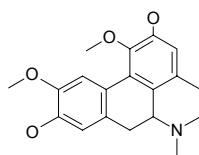
Brand	Catalog Number	MW
Aldrich	259810	309.37

**Sigma - Aldrich Building Blocks**

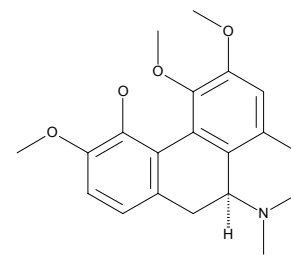
Brand	Catalog Number	MW
Aldrich	363944	321.81

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	469580	327.38

**Sigma - Aldrich Building Blocks**

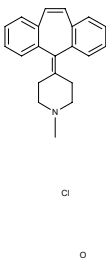
Brand	Catalog Number	MW
Aldrich	545368	341.41





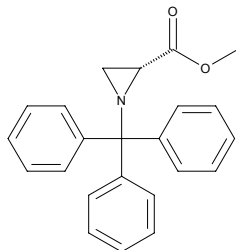
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	279072	341.88



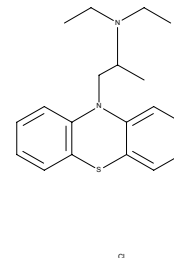
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	516007	343.43



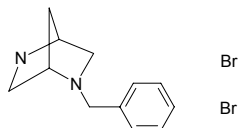
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	285838	348.94



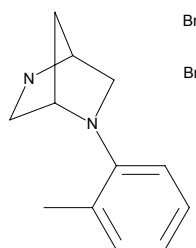
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	473030	350.10



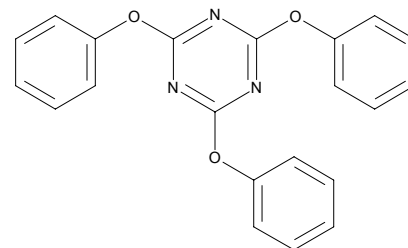
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	473413	350.10



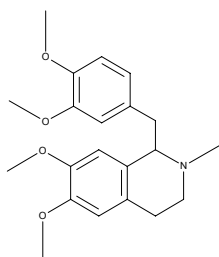
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	153664	357.37



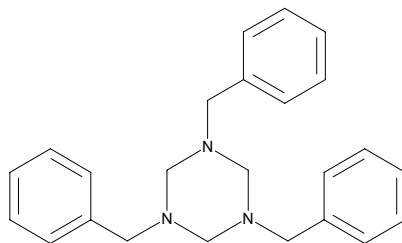
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	104671	357.45



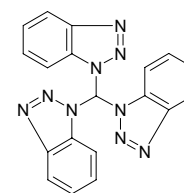
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	337722	357.50



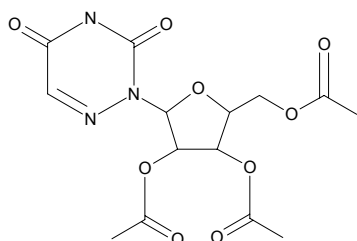
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	642568	367.38



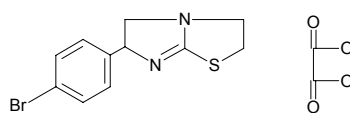
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	115185	371.31



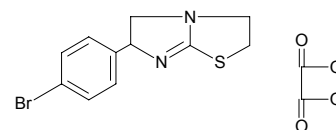
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	190047	373.23



Sigma - Aldrich Building Blocks

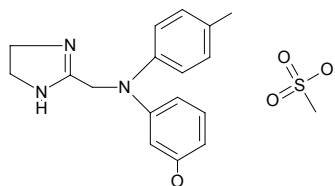
Brand	Catalog Number	MW
Aldrich	190055	373.23





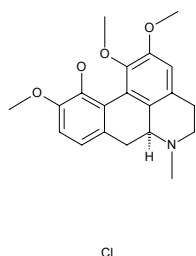
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	361658	377.47



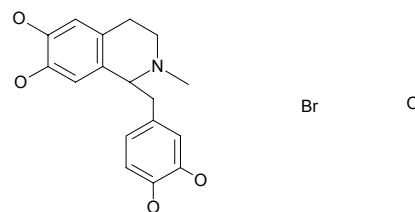
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	361607	377.87



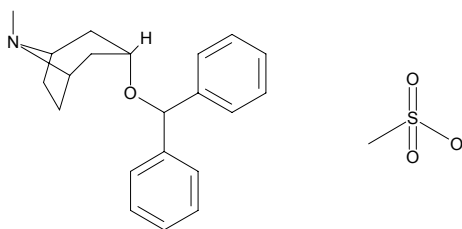
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	115770	400.27



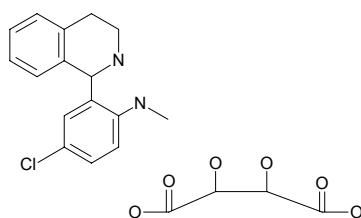
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	484628	403.54



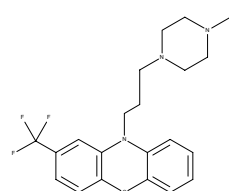
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	214345	422.87



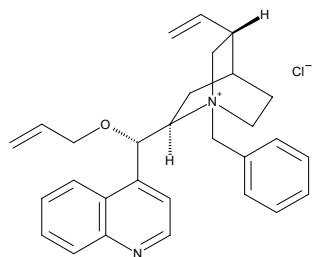
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	283886	443.97



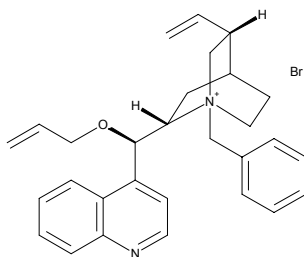
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	514268	461.05



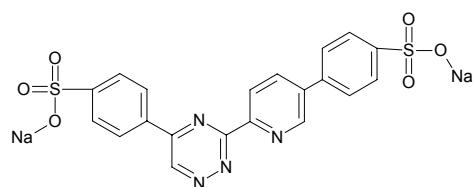
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	514276	505.50



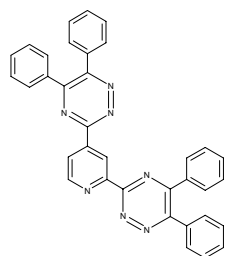
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	378143	514.45



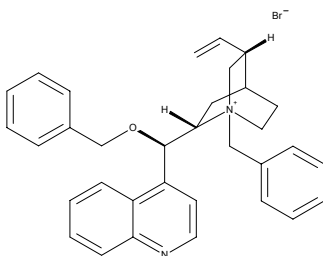
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	301752	541.62



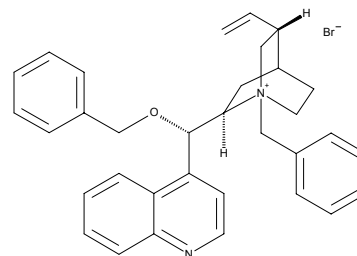
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	511552	555.56



Sigma - Aldrich Building Blocks

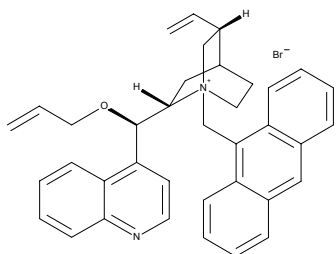
Brand	Catalog Number	MW
Aldrich	511811	555.56





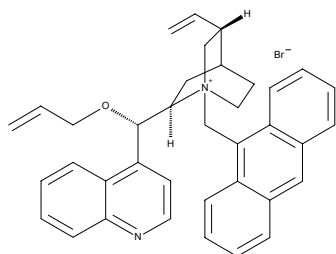
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	499617	605.62



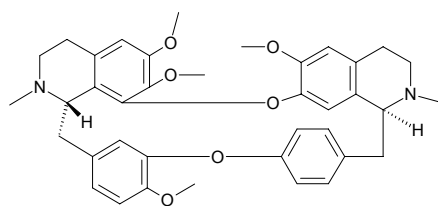
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	511838	605.62



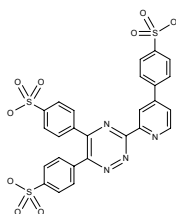
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	365629	622.77



**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	301876	650.64

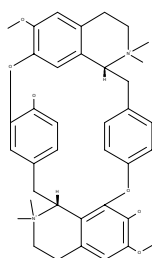


Na



**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	317330	713.71



**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	358797	1185.22

