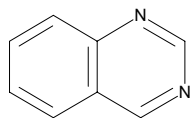




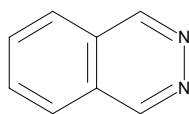
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	123323	130.15



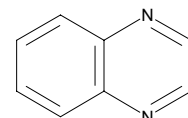
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	P38706	130.15



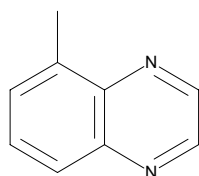
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	Q1603	130.15



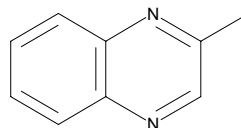
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	272310	144.18



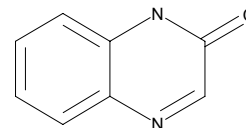
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	M80202	144.18



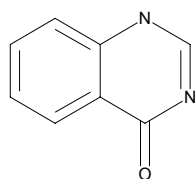
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	260517	146.15



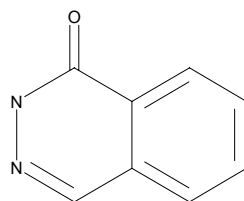
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	H57807	146.15



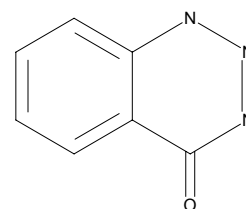
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	P39001	146.15



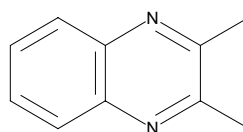
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	340545	147.14



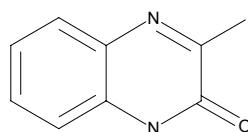
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	D184977	158.20



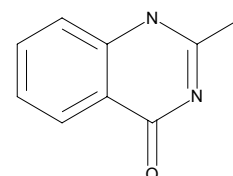
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	144460	160.18



Sigma - Aldrich Building Blocks

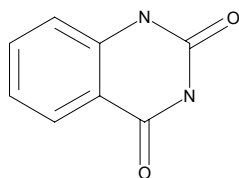
Brand	Catalog Number	MW
Aldrich	158836	160.18





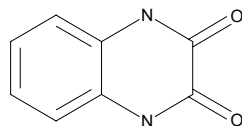
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	142026	162.15



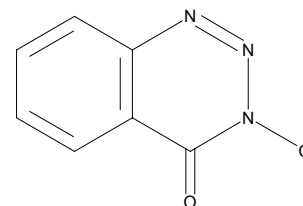
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	144789	162.15



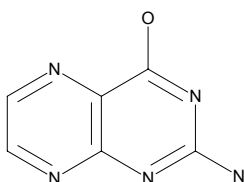
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	327964	163.14



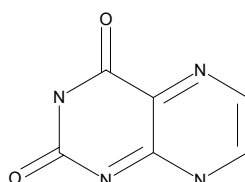
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	287083	163.14



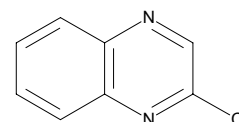
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	L3307	164.12



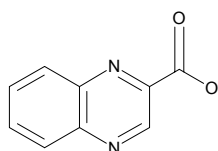
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	136301	164.60



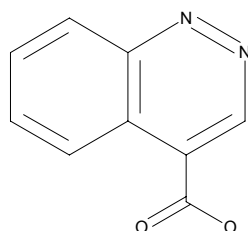
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	293407	174.16



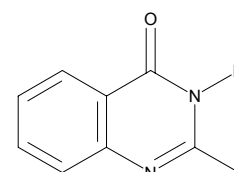
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	C82159	174.16



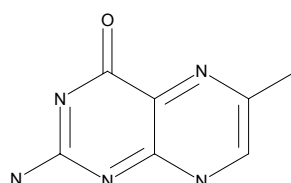
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	555797	175.19



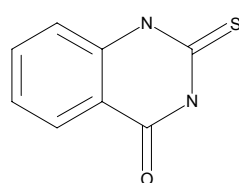
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	284793	177.17



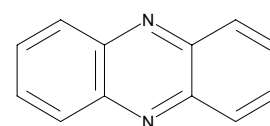
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	158828	178.21



Sigma - Aldrich Building Blocks

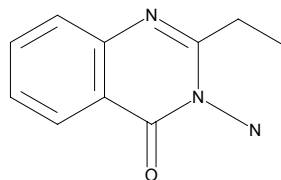
Brand	Catalog Number	MW
Aldrich	P13207	180.21





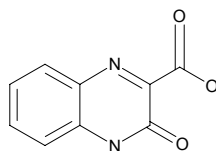
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	386324	189.22



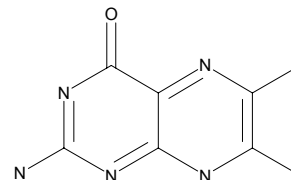
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	108340	190.16



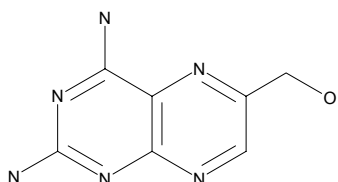
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	A51602	191.19



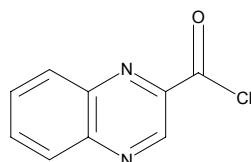
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	861634	192.18



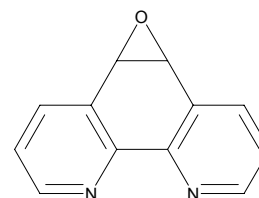
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	145998	192.61



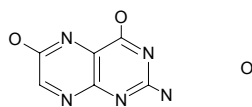
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	578789	196.21



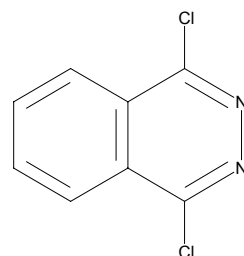
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	X708	197.15



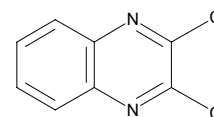
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	126020	199.04



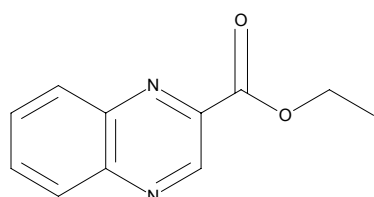
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	144525	199.04



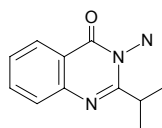
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	289051	202.21



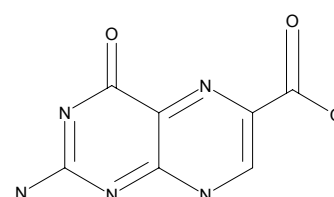
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	450642	203.25



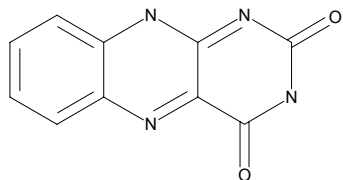
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	100080	207.15

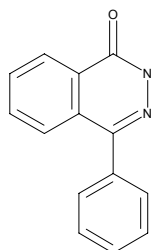


**Sigma - Aldrich Building Blocks**

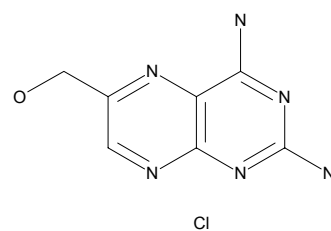
Brand	Catalog Number	MW
Aldrich	A28651	214.18

**Sigma - Aldrich Building Blocks**

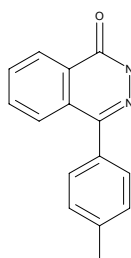
Brand	Catalog Number	MW
Aldrich	525553	222.25

**Sigma - Aldrich Building Blocks**

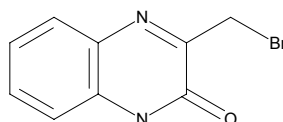
Brand	Catalog Number	MW
Aldrich	541575	228.64

**Sigma - Aldrich Building Blocks**

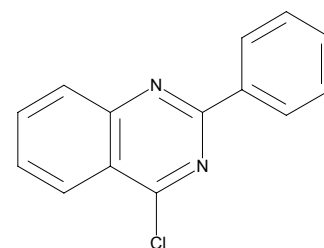
Brand	Catalog Number	MW
Aldrich	528080	236.28

**Sigma - Aldrich Building Blocks**

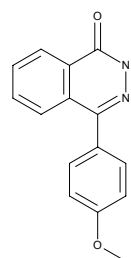
Brand	Catalog Number	MW
Aldrich	534218	239.07

**Sigma - Aldrich Building Blocks**

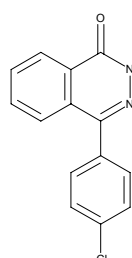
Brand	Catalog Number	MW
Aldrich	162434	240.69

**Sigma - Aldrich Building Blocks**

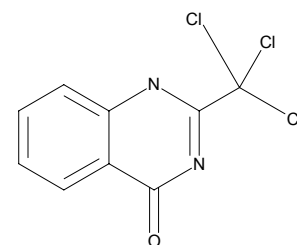
Brand	Catalog Number	MW
Aldrich	548839	252.28

**Sigma - Aldrich Building Blocks**

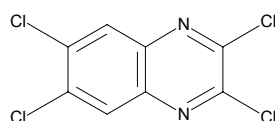
Brand	Catalog Number	MW
Aldrich	548820	256.69

**Sigma - Aldrich Building Blocks**

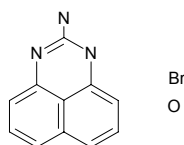
Brand	Catalog Number	MW
Aldrich	144592	263.51

**Sigma - Aldrich Building Blocks**

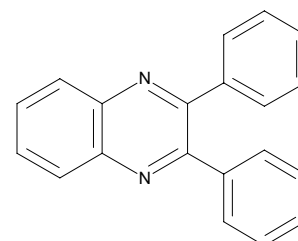
Brand	Catalog Number	MW
Aldrich	307149	267.93

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	232122	282.14

**Sigma - Aldrich Building Blocks**

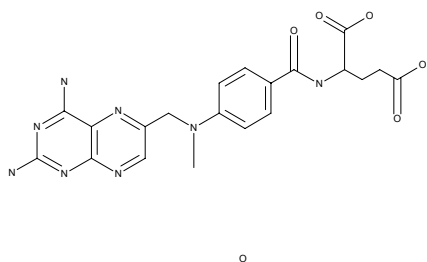
Brand	Catalog Number	MW
Aldrich	145483	282.35





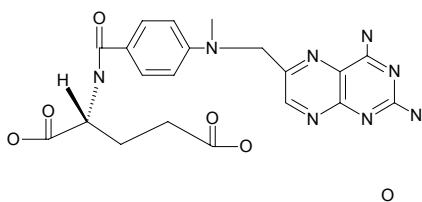
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	223948	472.46



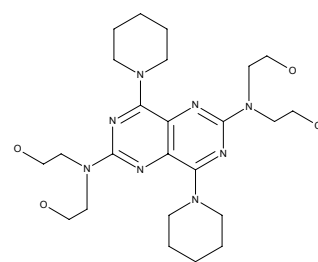
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	223956	472.46



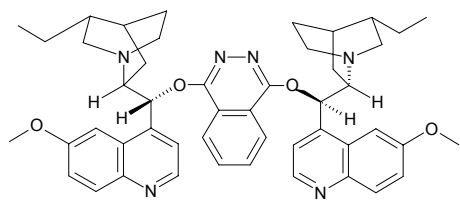
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	285676	504.64



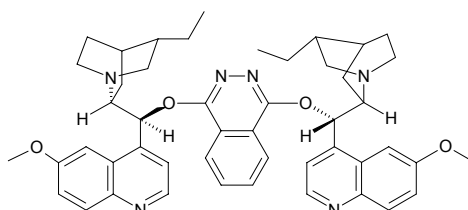
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	392723	779.00



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	392731	779.00



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	856258	1255.45

