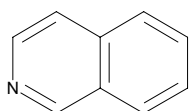
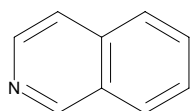


**Sigma - Aldrich Building Blocks**

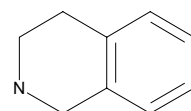
Brand	Catalog Number	MW
Aldrich	282162	129.16

**Sigma - Aldrich Building Blocks**

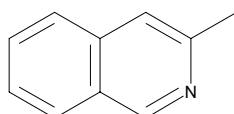
Brand	Catalog Number	MW
Aldrich	I28208	129.16

**Sigma - Aldrich Building Blocks**

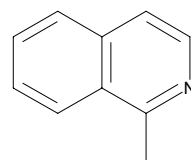
Brand	Catalog Number	MW
Aldrich	T13005	133.19

**Sigma - Aldrich Building Blocks**

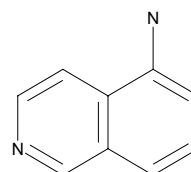
Brand	Catalog Number	MW
Aldrich	129895	143.19

**Sigma - Aldrich Building Blocks**

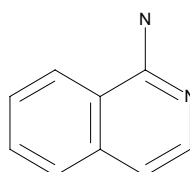
Brand	Catalog Number	MW
Aldrich	264938	143.19

**Sigma - Aldrich Building Blocks**

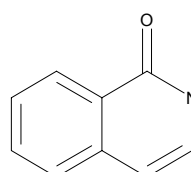
Brand	Catalog Number	MW
Aldrich	136107	144.18

**Sigma - Aldrich Building Blocks**

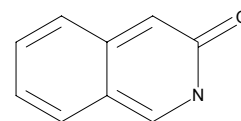
Brand	Catalog Number	MW
Aldrich	178594	144.18

**Sigma - Aldrich Building Blocks**

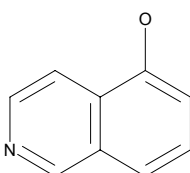
Brand	Catalog Number	MW
Aldrich	152102	145.16

**Sigma - Aldrich Building Blocks**

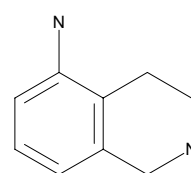
Brand	Catalog Number	MW
Aldrich	368954	145.16

**Sigma - Aldrich Building Blocks**

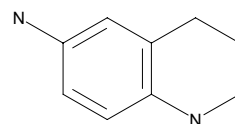
Brand	Catalog Number	MW
Aldrich	H33207	145.16

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	596825	148.21

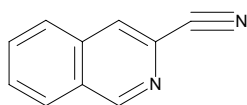
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	630276	148.21

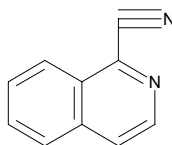


**Sigma - Aldrich Building Blocks**

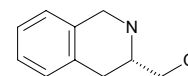
Brand	Catalog Number	MW
Aldrich	338532	154.17

**Sigma - Aldrich Building Blocks**

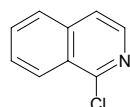
Brand	Catalog Number	MW
Aldrich	357952	154.17

**Sigma - Aldrich Building Blocks**

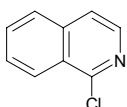
Brand	Catalog Number	MW
Aldrich	494186	163.22

**Sigma - Aldrich Building Blocks**

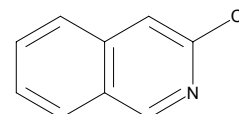
Brand	Catalog Number	MW
Aldrich	156744	163.61

**Sigma - Aldrich Building Blocks**

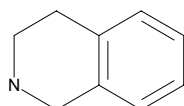
Brand	Catalog Number	MW
Aldrich	156744	163.61

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	595233	163.61

**Sigma - Aldrich Building Blocks**

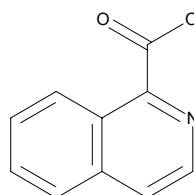
Brand	Catalog Number	MW
Aldrich	307548	169.66



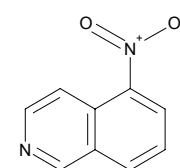
Cl

**Sigma - Aldrich Building Blocks**

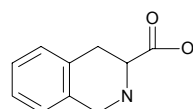
Brand	Catalog Number	MW
Aldrich	150134	173.17

**Sigma - Aldrich Building Blocks**

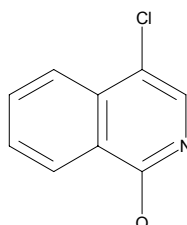
Brand	Catalog Number	MW
Aldrich	130222	174.16

**Sigma - Aldrich Building Blocks**

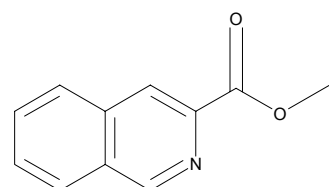
Brand	Catalog Number	MW
Aldrich	485039	177.20

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	156736	179.61

**Sigma - Aldrich Building Blocks**

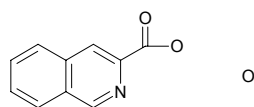
Brand	Catalog Number	MW
Aldrich	338559	187.20





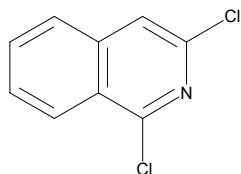
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	338540	191.19



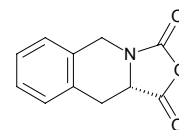
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	368830	198.05



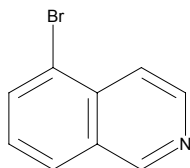
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	493759	203.20



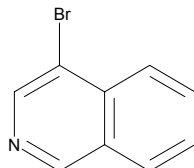
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	594261	208.06



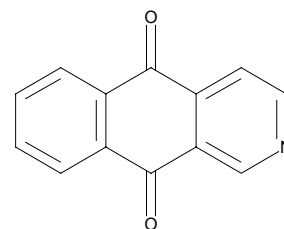
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	B70202	208.06



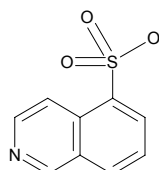
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	282219	209.21



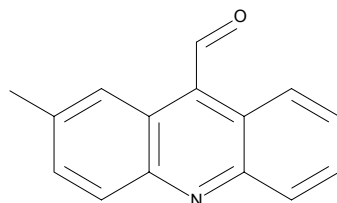
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	246123	209.23



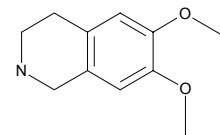
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	220604	221.26



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	291919	229.71

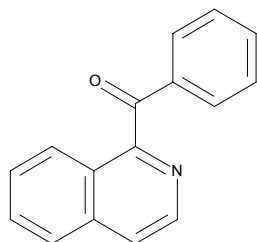


Cl



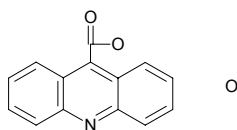
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	357960	233.27



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	246344	241.25

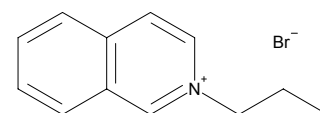


O



**Sigma - Aldrich Building Blocks**

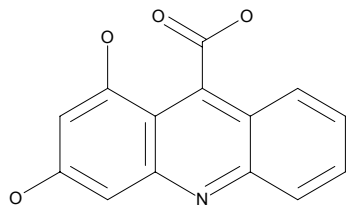
Brand	Catalog Number	MW
Aldrich	412872	252.16





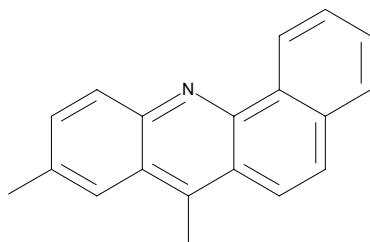
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	344990	255.23



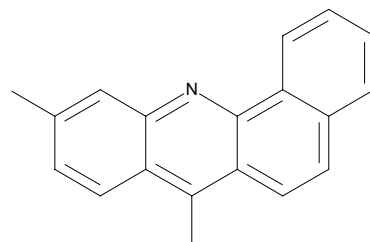
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	386332	257.34



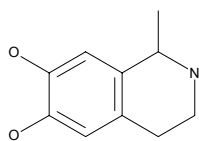
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	386677	257.34



### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	244201	260.13

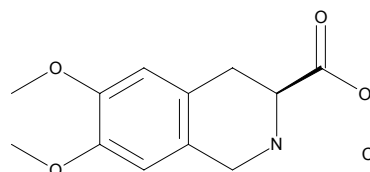


Br



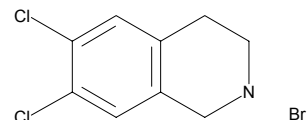
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	543527	273.72



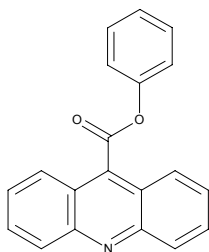
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	543616	283.00



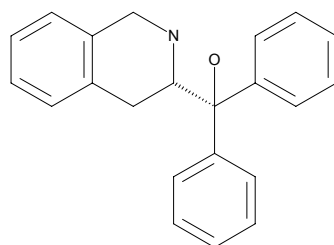
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	344168	299.33



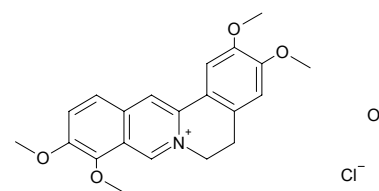
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	496057	315.42



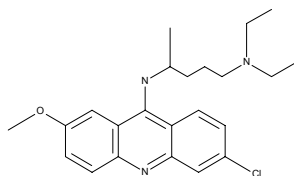
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	361615	405.88



### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	222992	436.43

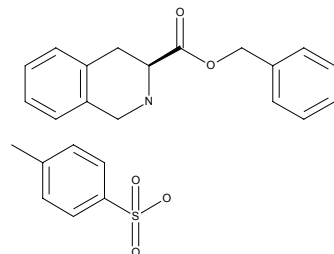


Cl



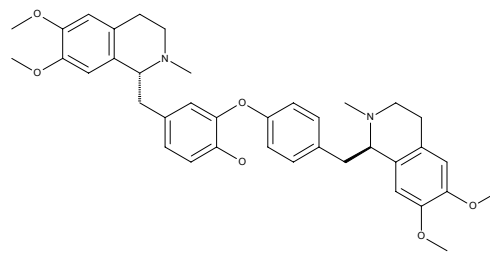
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	465062	439.53



### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	547255	624.78





**Sigma - Aldrich  
Building Blocks**

Brand	Catalog Number	MW
Aldrich	547190	681.66

