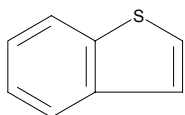




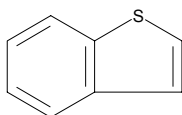
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	357790	134.20



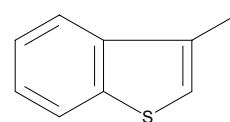
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	T27405	134.20



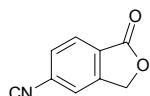
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	638587	148.23



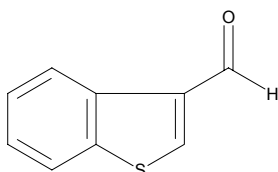
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	647195	159.15



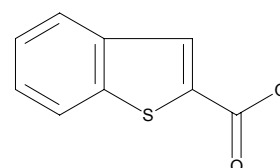
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	494968	162.21



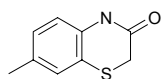
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	467464	178.21



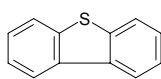
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	647519	179.24



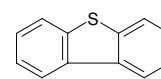
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	347833	184.26



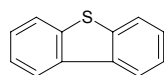
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	433306	184.26



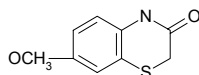
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	D32202	184.26



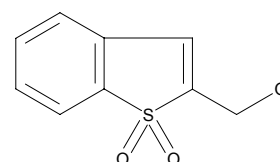
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	647500	195.24



**Sigma - Aldrich Building Blocks**

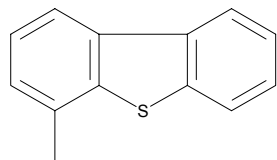
Brand	Catalog Number	MW
Aldrich	559164	196.23





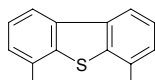
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	496901	198.29



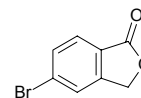
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	479411	212.32



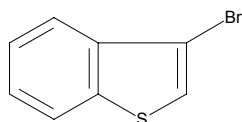
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	647187	213.03



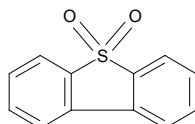
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	494976	213.10



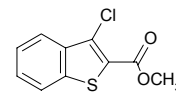
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	D32407	216.26



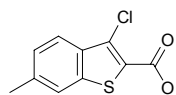
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	646741	226.68



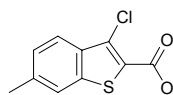
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	646814	226.68



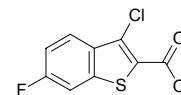
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	646814	226.68



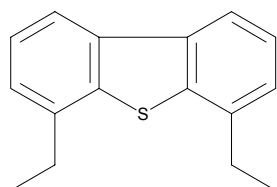
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	646806	230.65



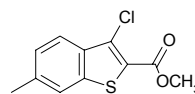
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	597996	240.37



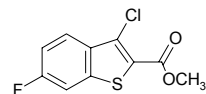
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	646784	240.71



**Sigma - Aldrich Building Blocks**

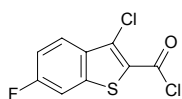
Brand	Catalog Number	MW
Aldrich	646776	244.67





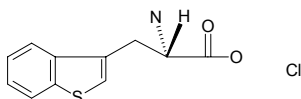
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	646822	249.09



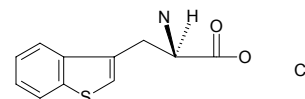
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	468827	257.74



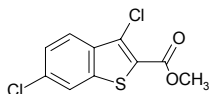
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	468835	257.74



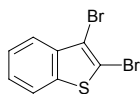
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	646792	261.13



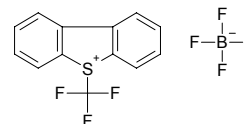
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	647373	291.99



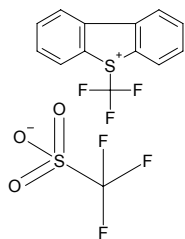
### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	483869	340.07



### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	483877	402.34



### Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	218057	445.41

