



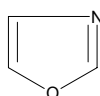
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	151637	69.06



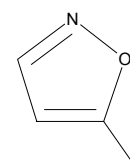
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	230138	69.06



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	153702	83.09



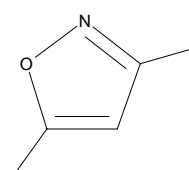
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	424218	84.08



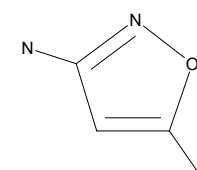
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	D167509	97.12



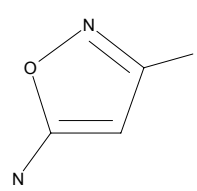
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	232270	98.11



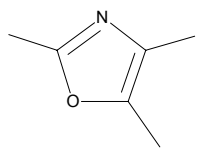
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	304271	98.11



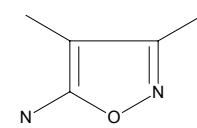
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	193704	111.14



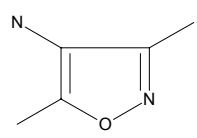
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	280240	112.13



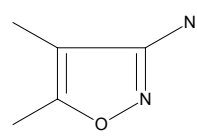
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	575069	112.13



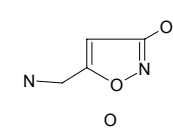
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	576050	112.13



Sigma - Aldrich Building Blocks

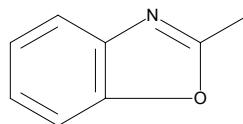
Brand	Catalog Number	MW
Aldrich	237795	132.12





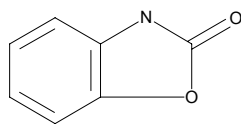
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	108936	133.15



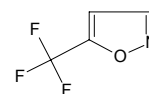
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	157058	135.12



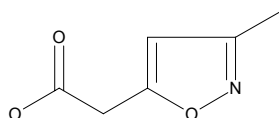
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	420638	137.06



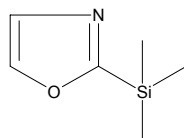
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	489689	141.13



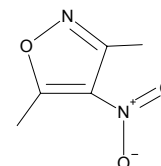
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	638609	141.25



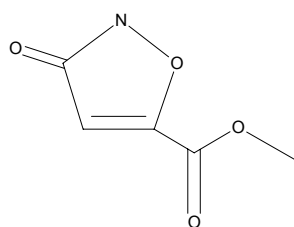
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	338249	142.12



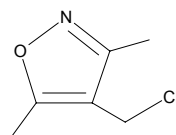
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	338605	143.10



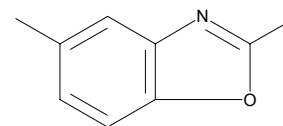
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	247316	145.59



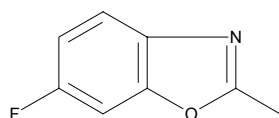
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	D150045	147.18



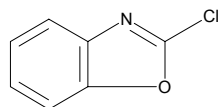
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	538434	151.14



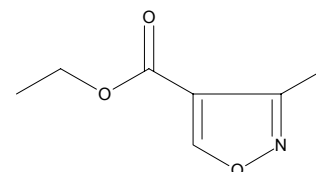
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	274089	153.57



Sigma - Aldrich Building Blocks

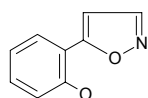
Brand	Catalog Number	MW
Aldrich	558362	155.15





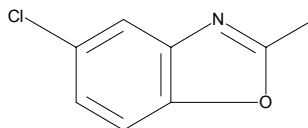
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	465429	161.16



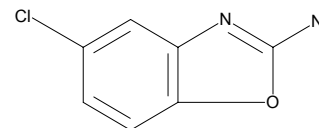
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	404500	167.60



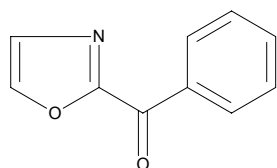
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	A45807	168.58



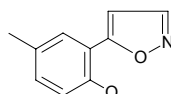
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	638595	173.17



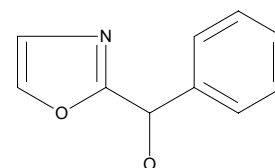
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	465453	175.19



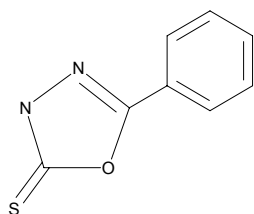
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	638544	175.19



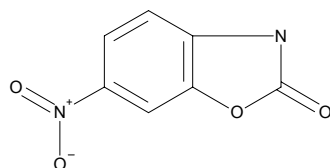
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	391808	178.21



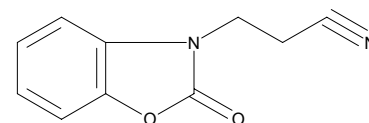
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	535605	180.12



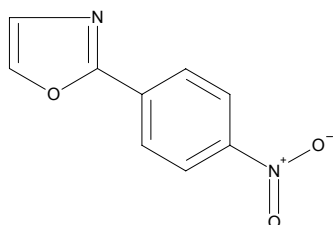
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	538531	188.19



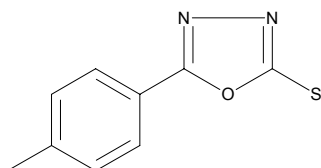
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	638633	190.16



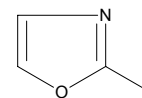
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	548952	192.24



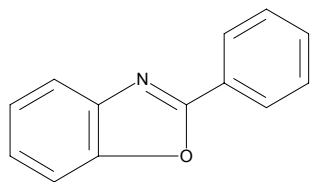
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	638552	194.96

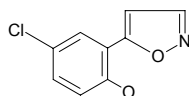


**Sigma - Aldrich Building Blocks**

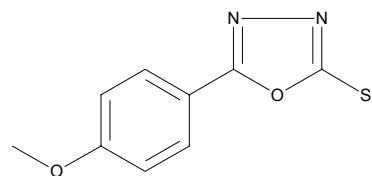
Brand	Catalog Number	MW
Aldrich	310565	195.22

**Sigma - Aldrich Building Blocks**

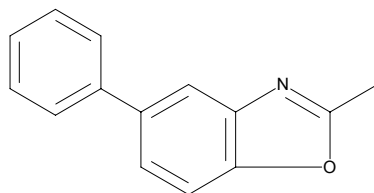
Brand	Catalog Number	MW
Aldrich	465437	195.61

**Sigma - Aldrich Building Blocks**

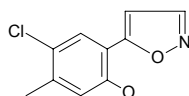
Brand	Catalog Number	MW
Aldrich	548960	208.24

**Sigma - Aldrich Building Blocks**

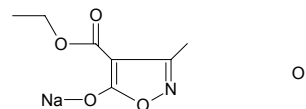
Brand	Catalog Number	MW
Aldrich	222259	209.25

**Sigma - Aldrich Building Blocks**

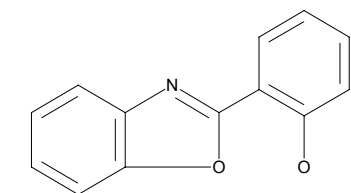
Brand	Catalog Number	MW
Aldrich	465488	209.63

**Sigma - Aldrich Building Blocks**

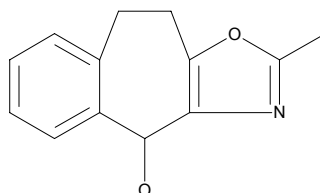
Brand	Catalog Number	MW
Aldrich	383643	211.15

**Sigma - Aldrich Building Blocks**

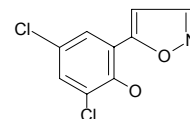
Brand	Catalog Number	MW
Aldrich	H50802	211.22

**Sigma - Aldrich Building Blocks**

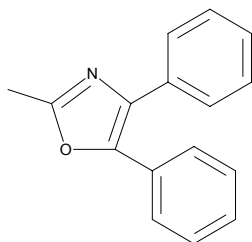
Brand	Catalog Number	MW
Aldrich	214329	215.25

**Sigma - Aldrich Building Blocks**

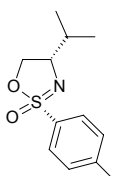
Brand	Catalog Number	MW
Aldrich	465461	230.05

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	225592	235.29

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	648345	239.34

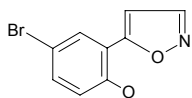
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	B96953	239.66

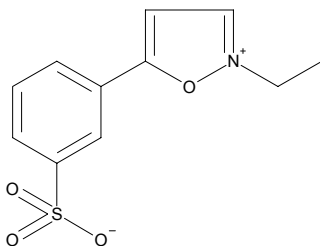


**Sigma - Aldrich Building Blocks**

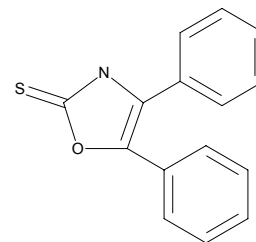
Brand	Catalog Number	MW
Aldrich	465496	240.06

**Sigma - Aldrich Building Blocks**

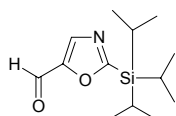
Brand	Catalog Number	MW
Aldrich	E45260	253.28

**Sigma - Aldrich Building Blocks**

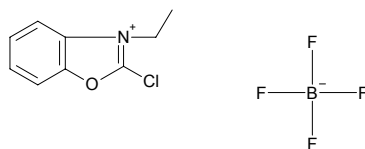
Brand	Catalog Number	MW
Aldrich	385719	253.33

**Sigma - Aldrich Building Blocks**

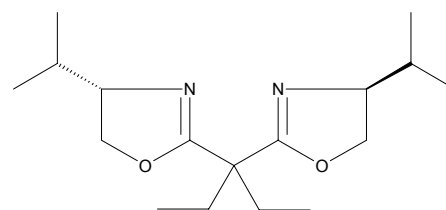
Brand	Catalog Number	MW
Aldrich	641103	253.42

**Sigma - Aldrich Building Blocks**

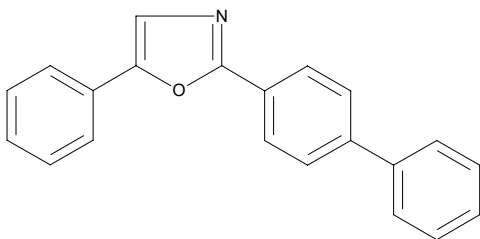
Brand	Catalog Number	MW
Aldrich	232556	269.44

**Sigma - Aldrich Building Blocks**

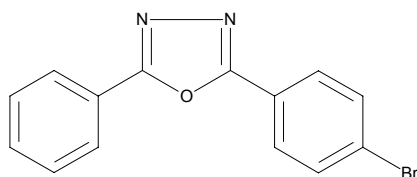
Brand	Catalog Number	MW
Aldrich	554138	294.44

**Sigma - Aldrich Building Blocks**

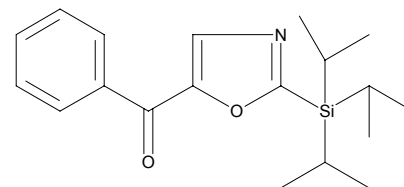
Brand	Catalog Number	MW
Aldrich	216984	297.36

**Sigma - Aldrich Building Blocks**

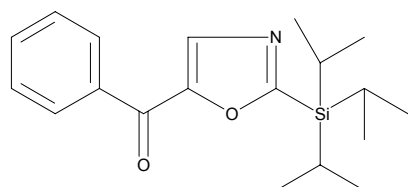
Brand	Catalog Number	MW
Aldrich	553646	301.14

**Sigma - Aldrich Building Blocks**

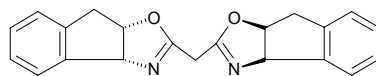
Brand	Catalog Number	MW
Aldrich	638625	329.52

**Sigma - Aldrich Building Blocks**

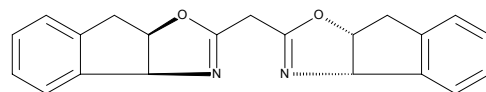
Brand	Catalog Number	MW
Aldrich	638625	329.52

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	464155	330.39

**Sigma - Aldrich Building Blocks**

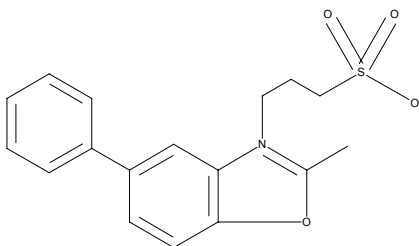
Brand	Catalog Number	MW
Aldrich	467073	330.39





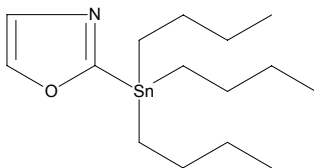
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	386081	333.41



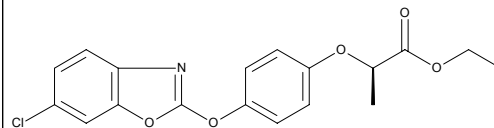
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	638617	358.09



**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	542776	361.78



**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	368598	414.47

