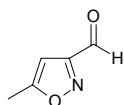




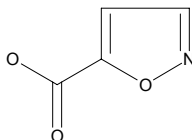
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	644684	111.10



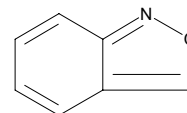
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	636258	113.07



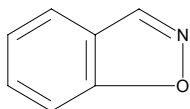
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	144517	119.12



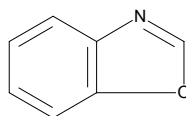
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	232300	119.12



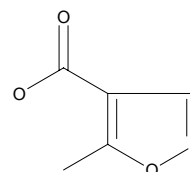
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	B11702	119.12



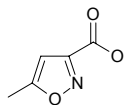
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	633771	127.10



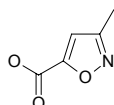
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	644676	127.10



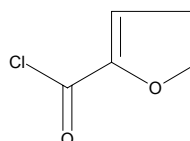
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	644900	127.10



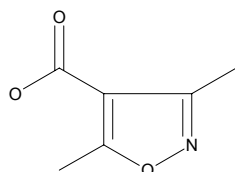
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	636231	131.52



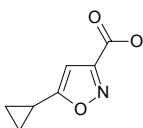
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	637653	141.13



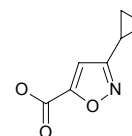
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	644803	153.14



**Sigma - Aldrich
Building Blocks**

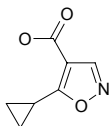
Brand	Catalog Number	MW
Aldrich	644811	153.14





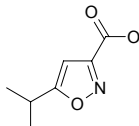
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	644838	153.14



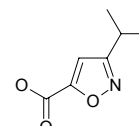
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	644730	155.15



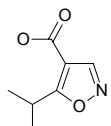
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	644854	155.15



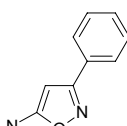
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	644862	155.15



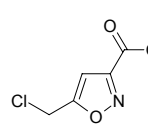
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	644978	160.18



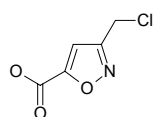
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	644749	161.55



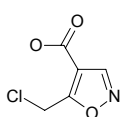
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	644757	161.55



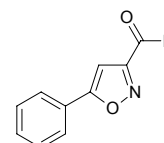
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	644765	161.55



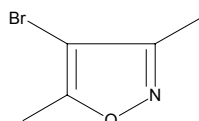
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	645060	173.17



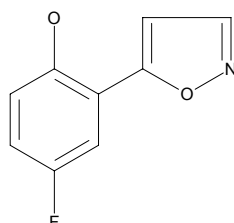
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	544299	176.01



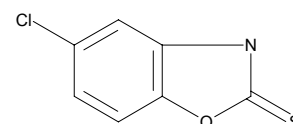
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	543683	179.15



Sigma - Aldrich Building Blocks

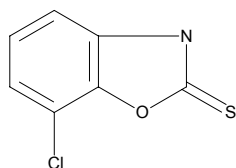
Brand	Catalog Number	MW
Aldrich	538426	185.63





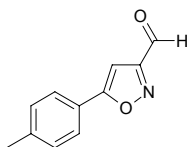
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	538442	185.63



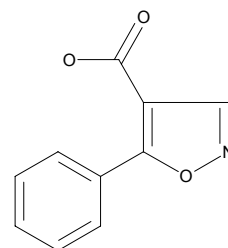
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	645044	187.20



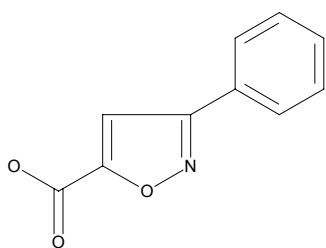
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	633682	189.17



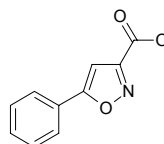
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	633690	189.17



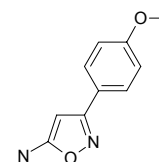
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	645087	189.17



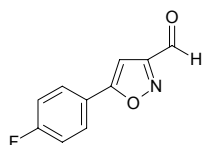
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	644951	190.20



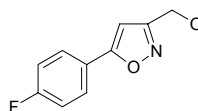
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	644668	191.16



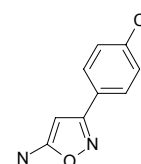
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	644641	193.18



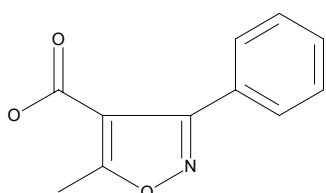
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	644943	194.62



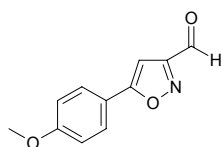
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	134198	203.20



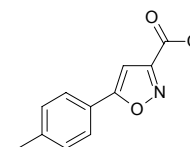
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	644617	203.20



Sigma - Aldrich Building Blocks

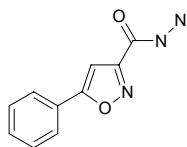
Brand	Catalog Number	MW
Aldrich	644986	203.20





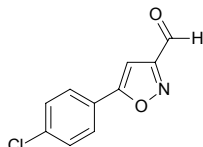
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	645079	203.20



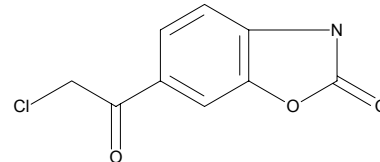
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	645036	207.62



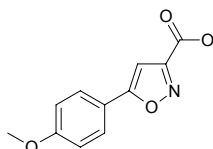
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	535400	211.61



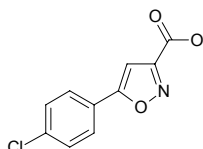
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	645028	219.20



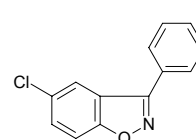
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	644994	223.62



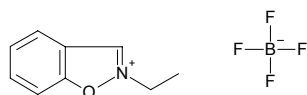
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	642657	229.67



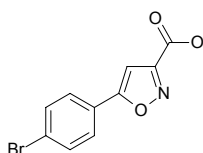
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	382647	234.99



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	645001	268.07



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	644625	282.10

