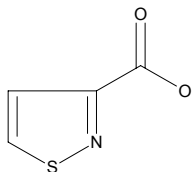




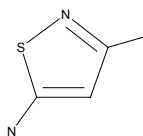
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	578258	129.14



**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	155640	150.63

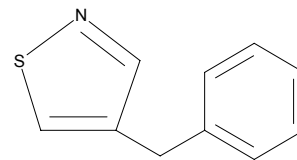


Cl



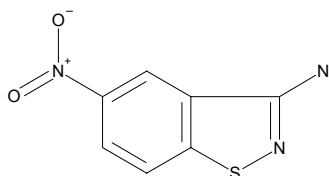
**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	305448	175.25



**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	522457	195.20



**Sigma - Aldrich
Building Blocks**

Brand	Catalog Number	MW
Aldrich	586366	528.43

