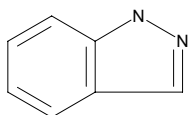




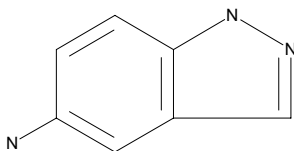
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	I2401	118.14



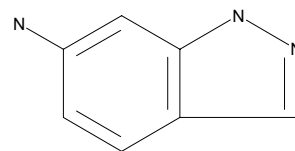
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	A59557	133.15



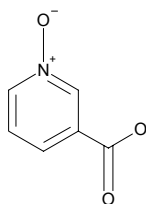
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	A59565	133.15



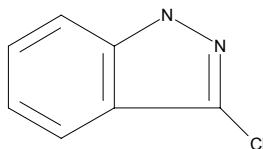
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	106194	139.11



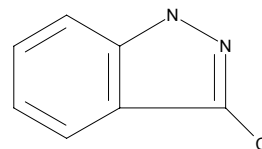
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	335207	152.58



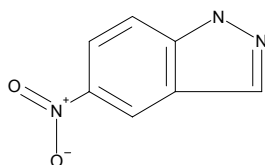
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	335207	152.58



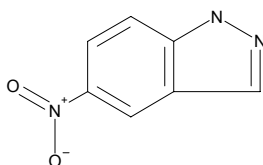
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	216755	163.14



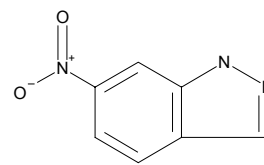
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	N17505	163.14



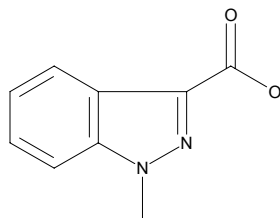
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	N17513	163.14



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	637645	176.18



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	561312	226.64

