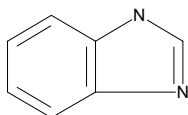




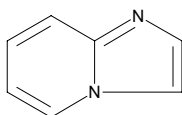
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	194123	118.14



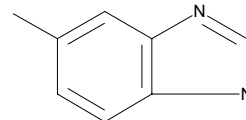
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	275778	118.14



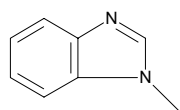
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	305235	132.17



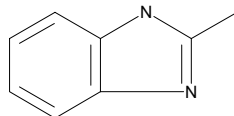
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	399353	132.17



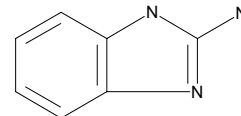
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	M29800	132.17



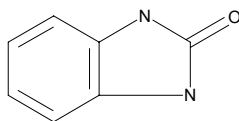
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	171778	133.15



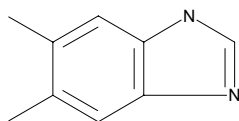
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	H19859	134.14



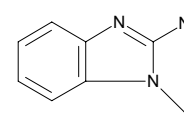
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	D147206	146.19



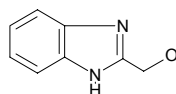
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	412546	147.18



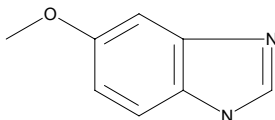
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	158356	148.17



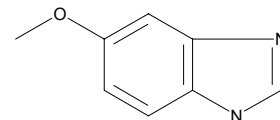
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	515108	148.17



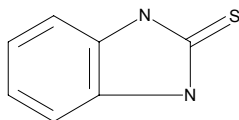
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	515108	148.17

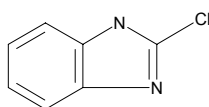


**Sigma - Aldrich Building Blocks**

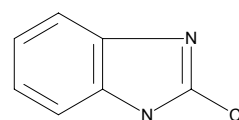
Brand	Catalog Number	MW
Aldrich	M3205	150.20

**Sigma - Aldrich Building Blocks**

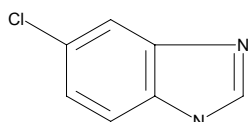
Brand	Catalog Number	MW
Aldrich	412538	152.58

**Sigma - Aldrich Building Blocks**

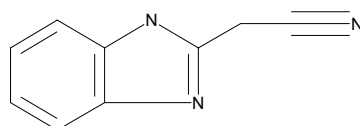
Brand	Catalog Number	MW
Aldrich	592277	152.58

**Sigma - Aldrich Building Blocks**

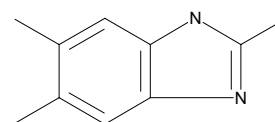
Brand	Catalog Number	MW
Aldrich	631248	152.58

**Sigma - Aldrich Building Blocks**

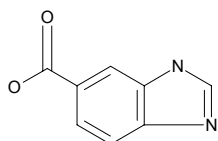
Brand	Catalog Number	MW
Aldrich	B5291	157.18

**Sigma - Aldrich Building Blocks**

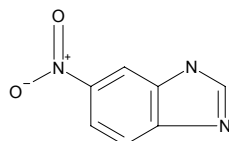
Brand	Catalog Number	MW
Aldrich	T74004	160.22

**Sigma - Aldrich Building Blocks**

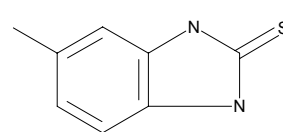
Brand	Catalog Number	MW
Aldrich	296783	162.15

**Sigma - Aldrich Building Blocks**

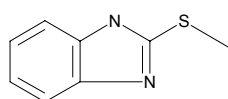
Brand	Catalog Number	MW
Aldrich	N11604	163.14

**Sigma - Aldrich Building Blocks**

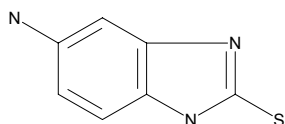
Brand	Catalog Number	MW
Aldrich	305243	164.23

**Sigma - Aldrich Building Blocks**

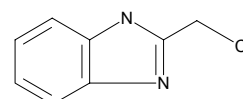
Brand	Catalog Number	MW
Aldrich	M54600	164.23

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	536067	165.22

**Sigma - Aldrich Building Blocks**

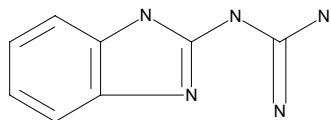
Brand	Catalog Number	MW
Aldrich	C52055	166.61





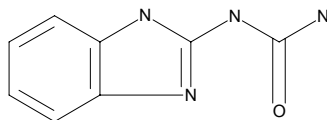
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	G11802	175.19



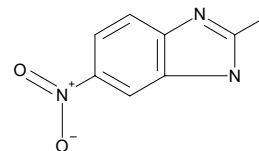
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	B5453	176.18



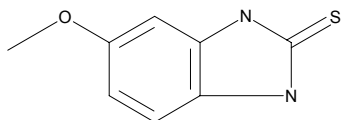
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	178225	177.16



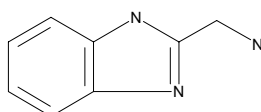
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	382485	180.23



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	165638	183.64

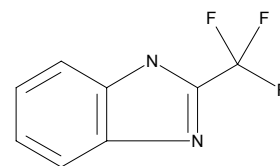


Cl



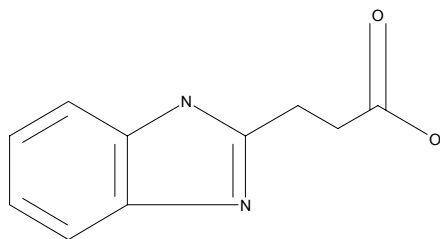
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	105694	186.14



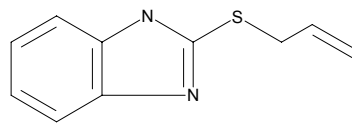
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	361933	190.20



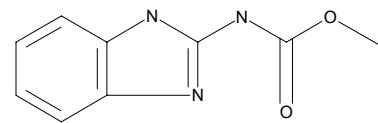
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	392065	190.27



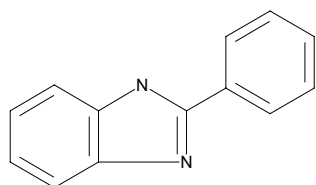
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	378674	191.19



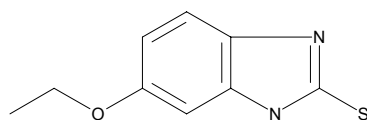
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	P19809	194.24



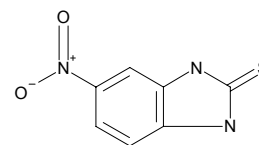
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	536075	194.26



Sigma - Aldrich Building Blocks

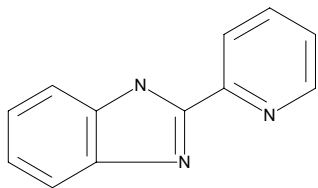
Brand	Catalog Number	MW
Aldrich	263575	195.20





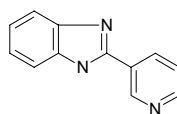
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	155101	195.23



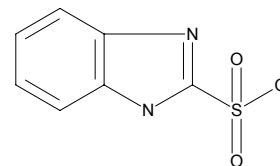
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	642711	195.23



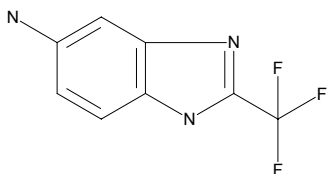
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	530646	198.20



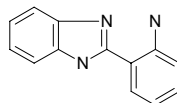
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	579483	201.15



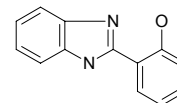
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	642681	209.25



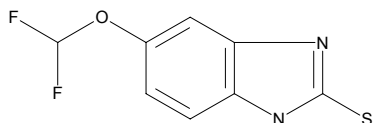
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	642738	210.24



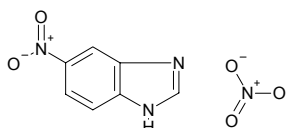
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	542210	216.21



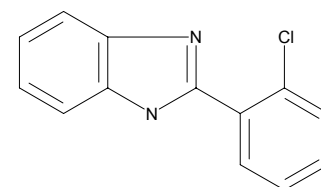
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	266507	226.15



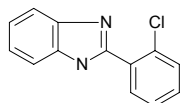
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	632600	228.68



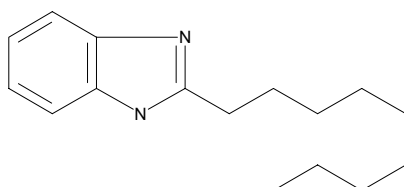
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	642703	228.68



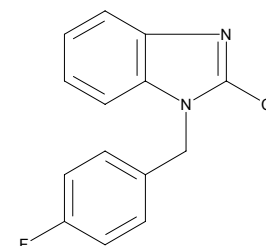
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	538353	244.38



Sigma - Aldrich Building Blocks

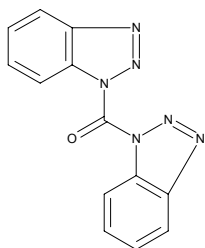
Brand	Catalog Number	MW
Aldrich	568694	260.70





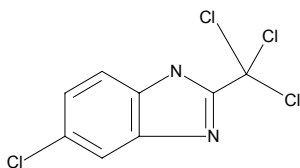
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	512974	264.25



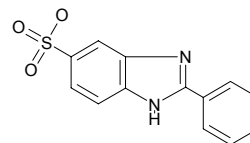
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	192538	269.95



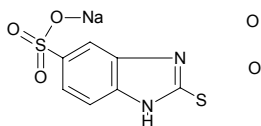
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	437166	274.30



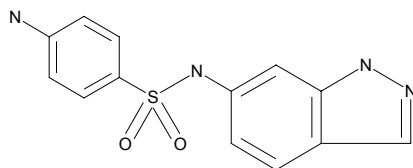
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	385700	288.28



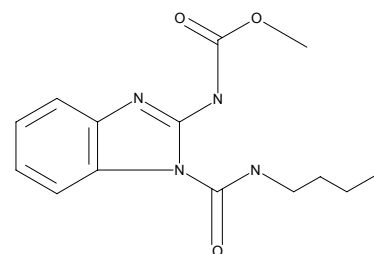
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	155306	288.33



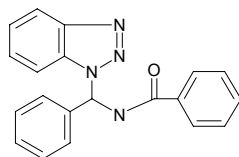
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	381586	290.32



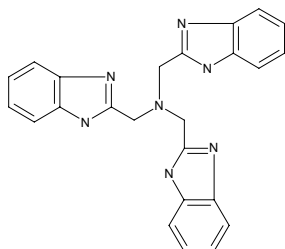
Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	452726	328.38



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	368520	407.48



Sigma - Aldrich Building Blocks

Brand	Catalog Number	MW
Aldrich	226084	442.29

