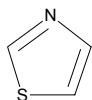




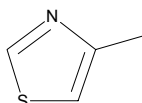
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	151645	85.13



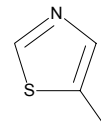
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	193925	99.16



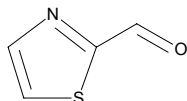
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	302740	99.16



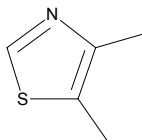
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	422460	113.14



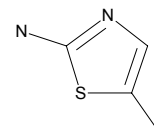
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	219177	113.18



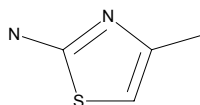
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	380563	114.17



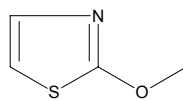
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	A66006	114.17



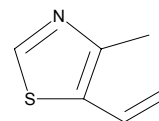
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	558532	115.15



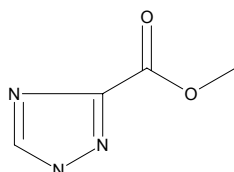
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	242004	125.19



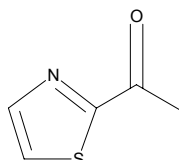
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	530352	127.10



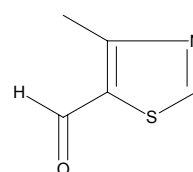
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	288411	127.17



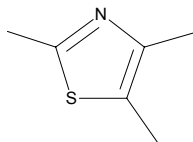
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	633208	127.17

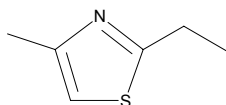


**Sigma - Aldrich Building Blocks**

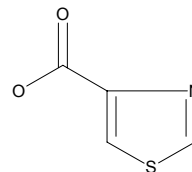
Brand	Catalog Number	MW
Aldrich	219185	127.21

**Sigma - Aldrich Building Blocks**

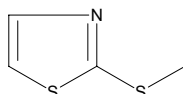
Brand	Catalog Number	MW
Aldrich	423599	127.21

**Sigma - Aldrich Building Blocks**

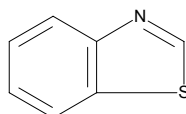
Brand	Catalog Number	MW
Aldrich	633658	129.14

**Sigma - Aldrich Building Blocks**

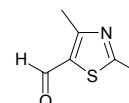
Brand	Catalog Number	MW
Aldrich	558524	131.22

**Sigma - Aldrich Building Blocks**

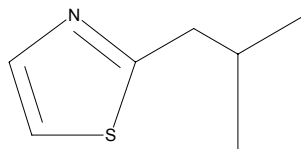
Brand	Catalog Number	MW
Aldrich	101338	135.19

**Sigma - Aldrich Building Blocks**

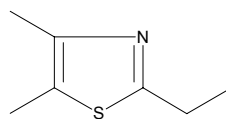
Brand	Catalog Number	MW
Aldrich	646024	141.19

**Sigma - Aldrich Building Blocks**

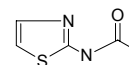
Brand	Catalog Number	MW
Aldrich	193682	141.24

**Sigma - Aldrich Building Blocks**

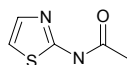
Brand	Catalog Number	MW
Aldrich	556718	141.24

**Sigma - Aldrich Building Blocks**

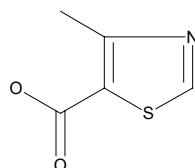
Brand	Catalog Number	MW
Aldrich	478725	142.18

**Sigma - Aldrich Building Blocks**

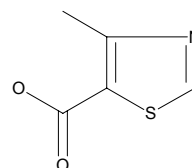
Brand	Catalog Number	MW
Aldrich	648388	142.18

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	633666	143.17

**Sigma - Aldrich Building Blocks**

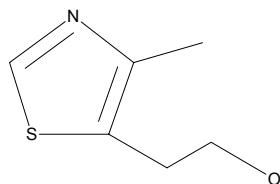
Brand	Catalog Number	MW
Aldrich	633666	143.17





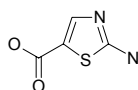
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	190675	143.21



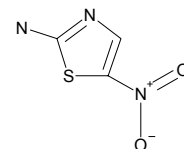
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	642479	144.15



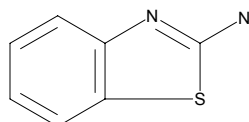
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	133507	145.14



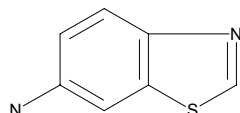
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	108812	150.20



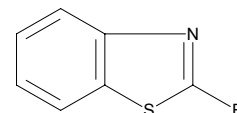
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	562440	150.20



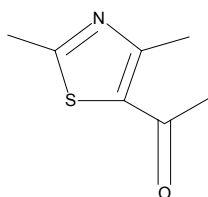
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	553085	153.18



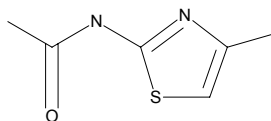
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	298085	155.22



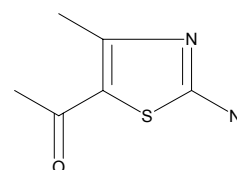
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	301922	156.21



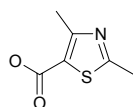
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	586145	156.21



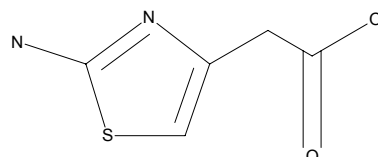
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	642509	157.19



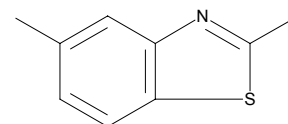
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	249696	158.18



**Sigma - Aldrich Building Blocks**

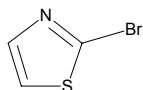
Brand	Catalog Number	MW
Aldrich	108839	163.24





**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	160474	164.02



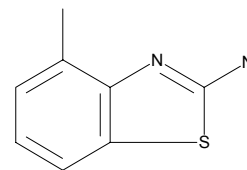
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	642517	164.02



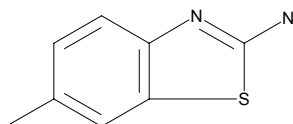
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	193224	164.23



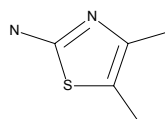
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	288373	164.23



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	174408	164.66

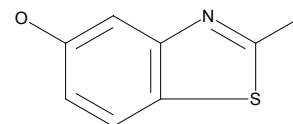


Cl



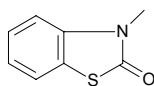
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	413259	165.22



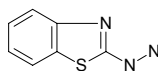
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	460281	165.22



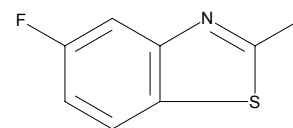
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	642436	165.22



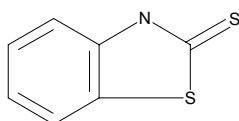
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	304476	167.21



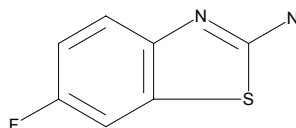
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	M3302	167.25



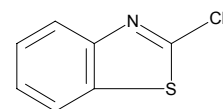
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	324213	168.19



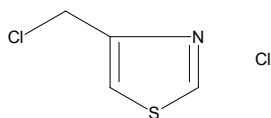
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	167576	169.63

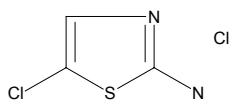


**Sigma - Aldrich Building Blocks**

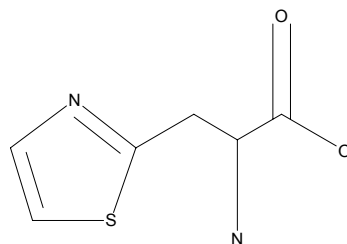
Brand	Catalog Number	MW
Aldrich	534277	170.06

**Sigma - Aldrich Building Blocks**

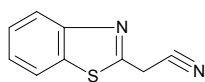
Brand	Catalog Number	MW
Aldrich	143456	171.05

**Sigma - Aldrich Building Blocks**

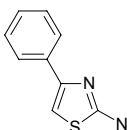
Brand	Catalog Number	MW
Aldrich	862193	172.21

**Sigma - Aldrich Building Blocks**

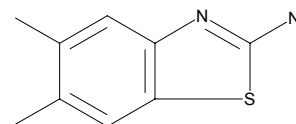
Brand	Catalog Number	MW
Aldrich	443042	174.23

**Sigma - Aldrich Building Blocks**

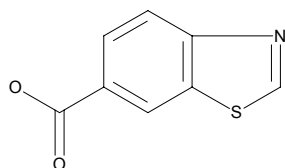
Brand	Catalog Number	MW
Aldrich	642754	176.24

**Sigma - Aldrich Building Blocks**

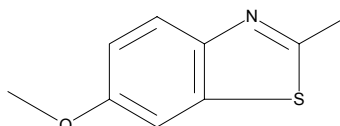
Brand	Catalog Number	MW
Aldrich	A51408	178.26

**Sigma - Aldrich Building Blocks**

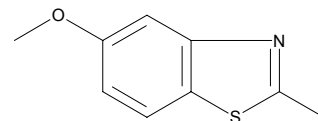
Brand	Catalog Number	MW
Aldrich	530336	179.20

**Sigma - Aldrich Building Blocks**

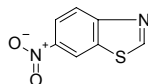
Brand	Catalog Number	MW
Aldrich	129682	179.24

**Sigma - Aldrich Building Blocks**

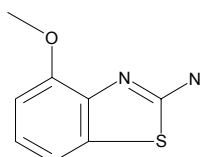
Brand	Catalog Number	MW
Aldrich	545767	179.24

**Sigma - Aldrich Building Blocks**

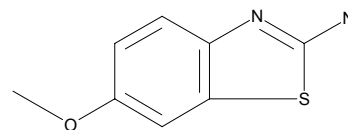
Brand	Catalog Number	MW
Aldrich	469114	180.19

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	138215	180.23

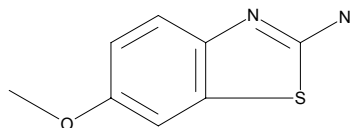
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	162590	180.23

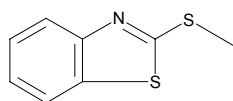


**Sigma - Aldrich Building Blocks**

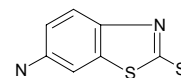
Brand	Catalog Number	MW
Aldrich	A60806	180.23

**Sigma - Aldrich Building Blocks**

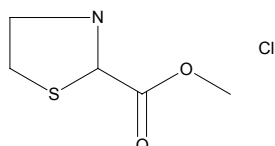
Brand	Catalog Number	MW
Aldrich	168653	181.28

**Sigma - Aldrich Building Blocks**

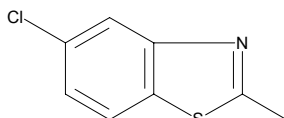
Brand	Catalog Number	MW
Aldrich	454370	182.27

**Sigma - Aldrich Building Blocks**

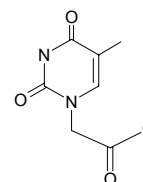
Brand	Catalog Number	MW
Aldrich	548758	183.66

**Sigma - Aldrich Building Blocks**

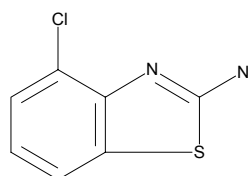
Brand	Catalog Number	MW
Aldrich	545775	183.66

**Sigma - Aldrich Building Blocks**

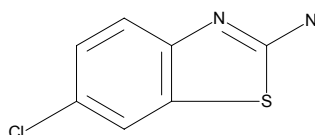
Brand	Catalog Number	MW
Aldrich	448958	184.15

**Sigma - Aldrich Building Blocks**

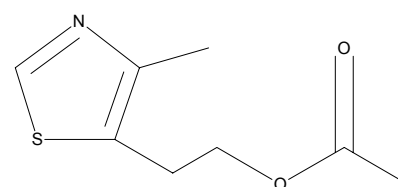
Brand	Catalog Number	MW
Aldrich	133434	184.65

**Sigma - Aldrich Building Blocks**

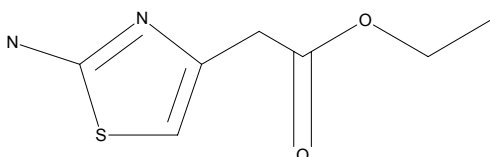
Brand	Catalog Number	MW
Aldrich	136085	184.65

**Sigma - Aldrich Building Blocks**

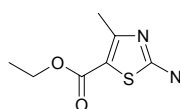
Brand	Catalog Number	MW
Aldrich	311316	185.25

**Sigma - Aldrich Building Blocks**

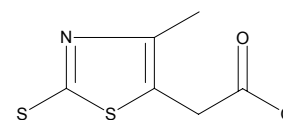
Brand	Catalog Number	MW
Aldrich	220558	186.23

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	646016	186.23

**Sigma - Aldrich Building Blocks**

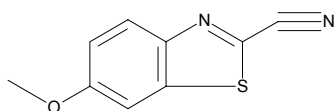
Brand	Catalog Number	MW
Aldrich	522317	189.26





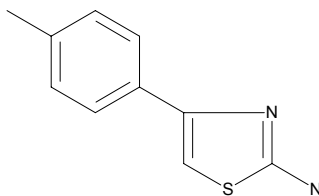
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	261858	190.22



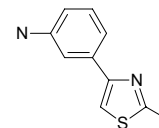
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	534161	190.27



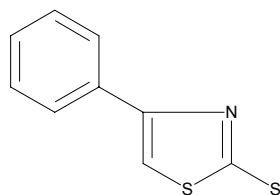
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	642630	190.27



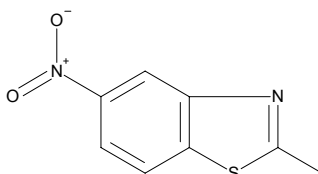
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	632570	193.29



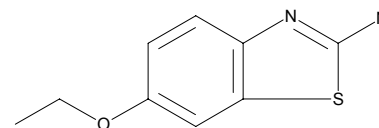
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	545821	194.21



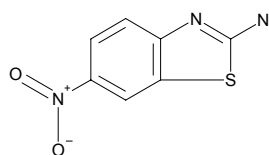
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	A54008	194.26



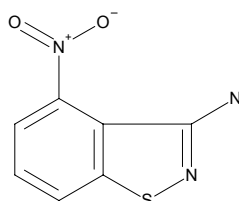
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	190241	195.20



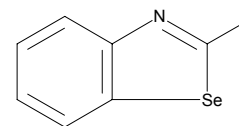
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	550884	195.20



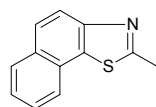
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	M30604	196.11



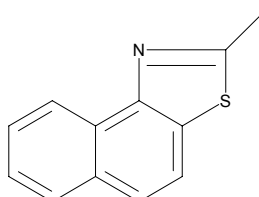
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	641847	199.28



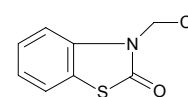
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	M57804	199.28



**Sigma - Aldrich Building Blocks**

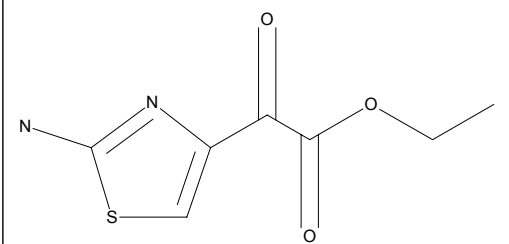
Brand	Catalog Number	MW
Aldrich	460303	199.66





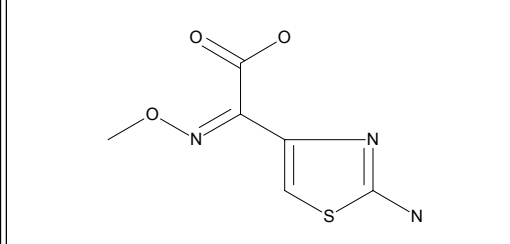
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	280062	200.22



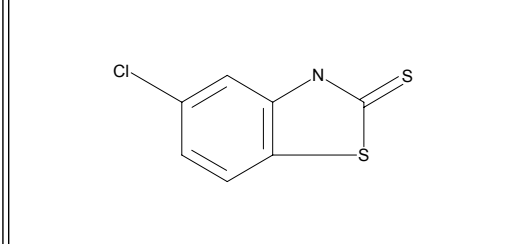
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	280143	201.20



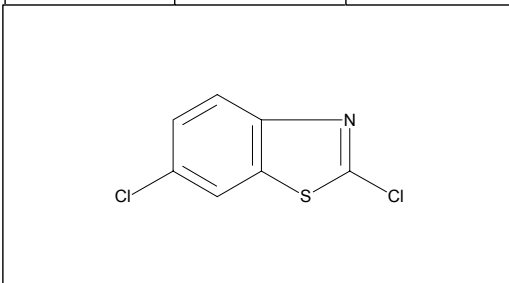
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	125571	201.70



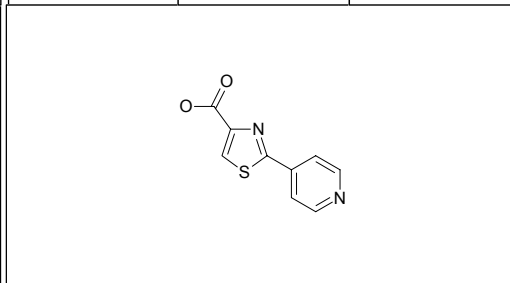
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	553670	204.08



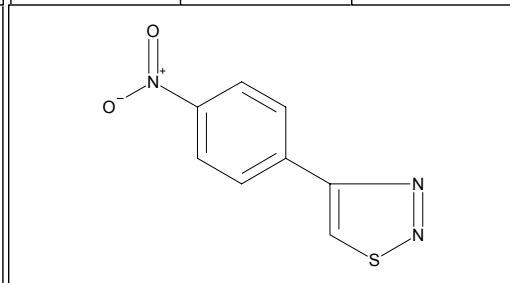
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	646032	206.22



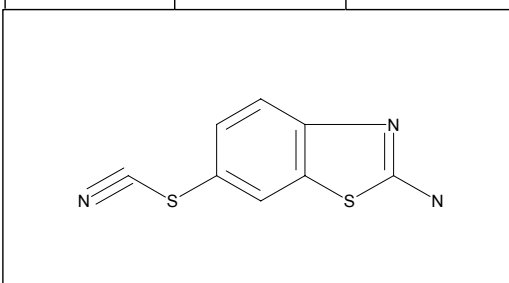
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	558494	207.21



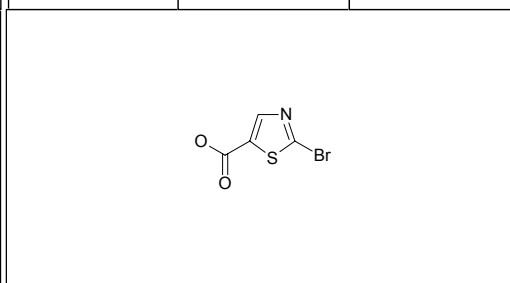
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	546887	207.28



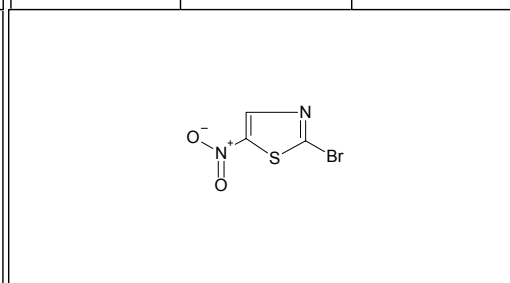
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	642487	208.03



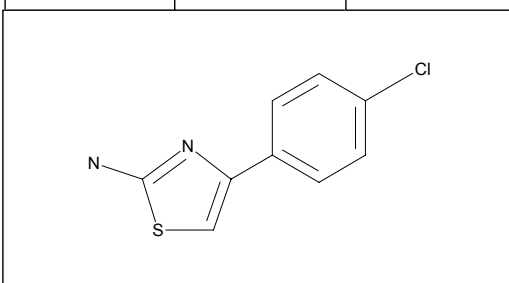
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	B74372	209.02



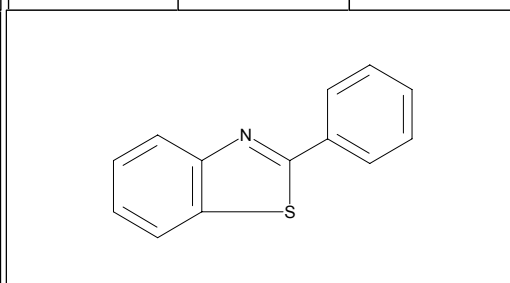
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	339369	210.69



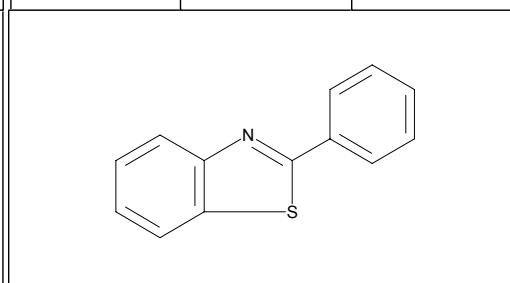
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	225444	211.29



**Sigma - Aldrich Building Blocks**

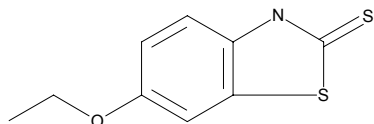
Brand	Catalog Number	MW
Aldrich	268437	211.29





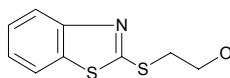
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	142174	211.31



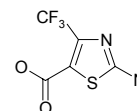
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	454451	211.31



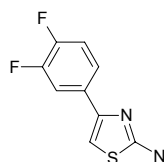
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	646059	212.15



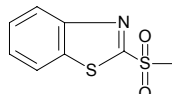
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	642444	212.22



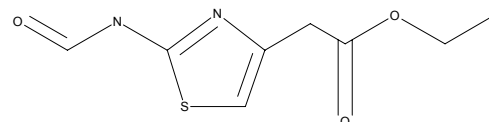
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	454354	213.28



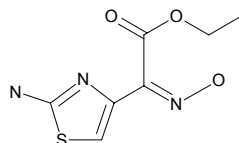
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	279757	214.24



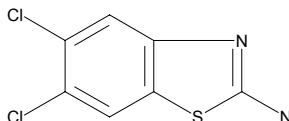
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	280178	215.23



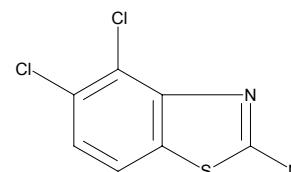
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	546852	219.09



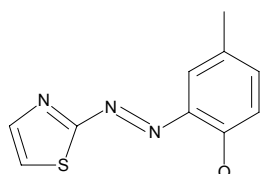
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	550876	219.09



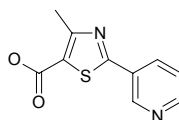
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	294543	219.27



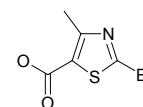
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	646040	220.25



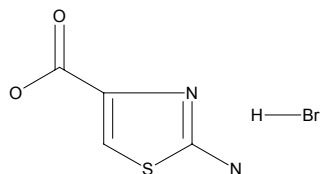
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	646008	222.06

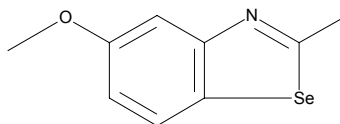


**Sigma - Aldrich Building Blocks**

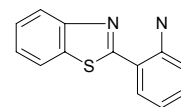
Brand	Catalog Number	MW
Aldrich	636282	225.06

**Sigma - Aldrich Building Blocks**

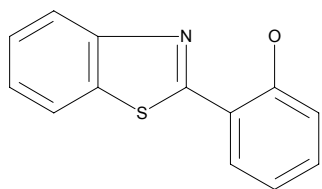
Brand	Catalog Number	MW
Aldrich	M15206	226.14

**Sigma - Aldrich Building Blocks**

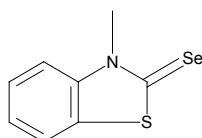
Brand	Catalog Number	MW
Aldrich	642428	226.30

**Sigma - Aldrich Building Blocks**

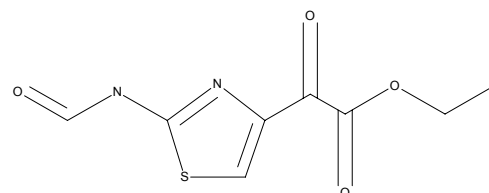
Brand	Catalog Number	MW
Aldrich	632589	227.29

**Sigma - Aldrich Building Blocks**

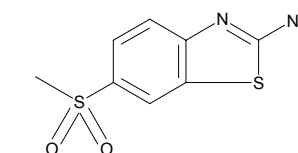
Brand	Catalog Number	MW
Aldrich	308978	228.18

**Sigma - Aldrich Building Blocks**

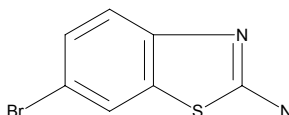
Brand	Catalog Number	MW
Aldrich	280054	228.23

**Sigma - Aldrich Building Blocks**

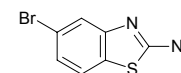
Brand	Catalog Number	MW
Aldrich	419478	228.29

**Sigma - Aldrich Building Blocks**

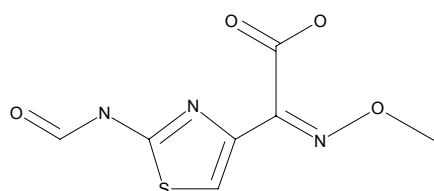
Brand	Catalog Number	MW
Aldrich	538507	229.10

**Sigma - Aldrich Building Blocks**

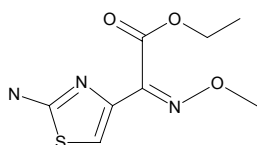
Brand	Catalog Number	MW
Aldrich	647683	229.10

**Sigma - Aldrich Building Blocks**

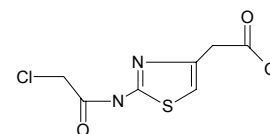
Brand	Catalog Number	MW
Aldrich	280194	229.22

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	280151	229.26

**Sigma - Aldrich Building Blocks**

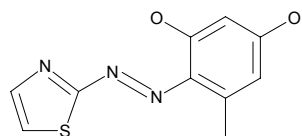
Brand	Catalog Number	MW
Aldrich	448117	234.66





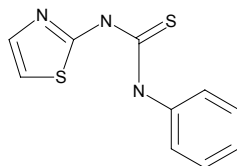
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	323179	235.27



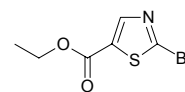
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	157961	235.33



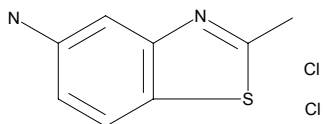
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	642495	236.09



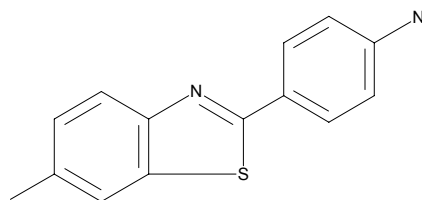
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	A63309	237.15



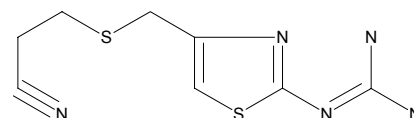
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	A72995	240.33



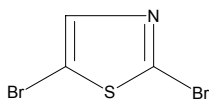
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	546879	241.34



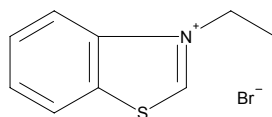
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	524182	242.92



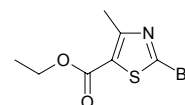
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	558443	244.15



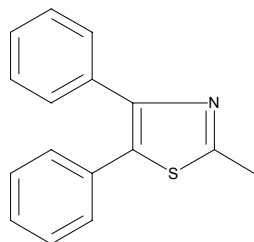
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	645990	250.12



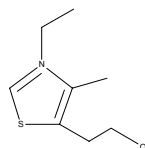
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	536059	251.35



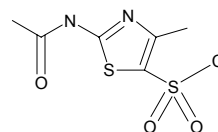
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	331244	254.19



**Sigma - Aldrich Building Blocks**

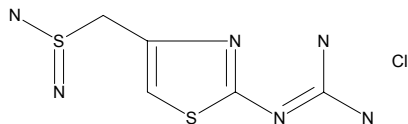
Brand	Catalog Number	MW
Aldrich	107859	254.72





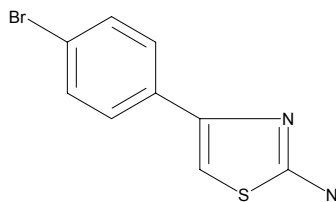
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	540269	254.76



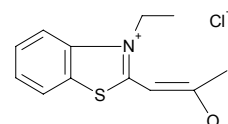
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	534129	255.14



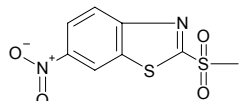
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	293652	255.77



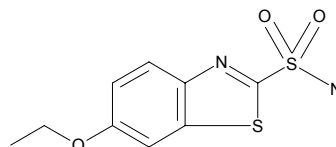
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	454362	258.28



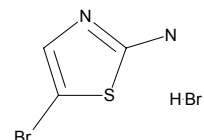
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	333328	258.32



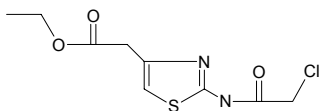
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	128023	259.95



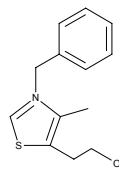
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	445436	262.72



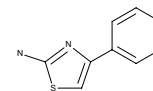
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	256234	271.81



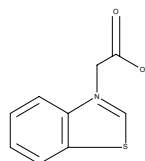
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	A75005	275.17



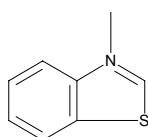
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	371637	276.15



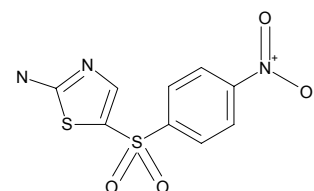
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	385972	279.14



**Sigma - Aldrich Building Blocks**

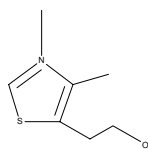
Brand	Catalog Number	MW
Aldrich	A70755	285.30





**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	257826	287.16

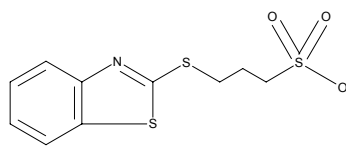


i



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	251690	313.39

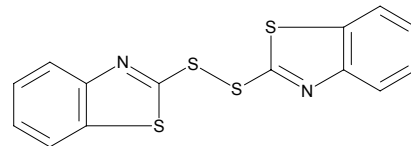


Na



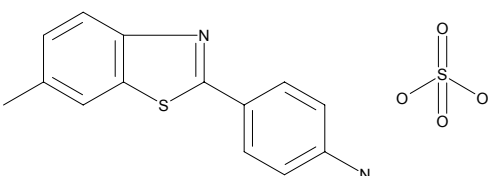
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	D218154	332.49



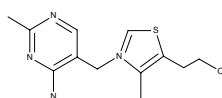
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	555789	338.41



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	109177	339.29



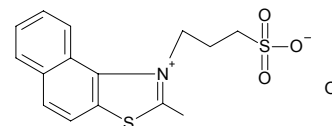
Cl

Cl



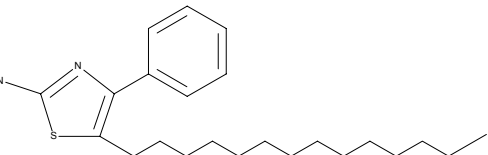
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	381233	339.44



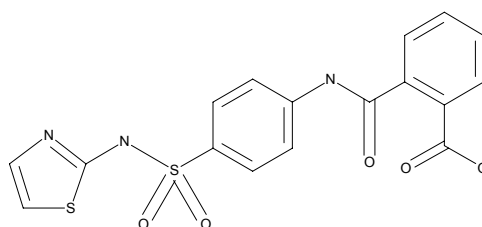
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	141054	372.62



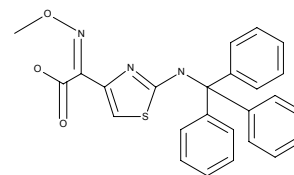
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	211834	403.44



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	280186	479.99



Cl