



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	516694	84.14



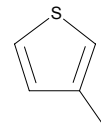
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	T31801	84.14



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	M84402	98.17



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	479209	104.17



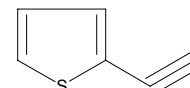
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	479217	104.17



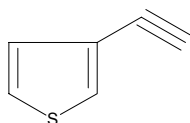
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	457469	108.16



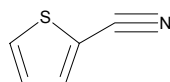
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	578797	108.16



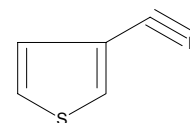
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	154385	109.15



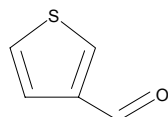
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	597449	109.15



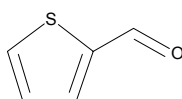
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	196282	112.15



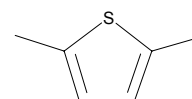
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	T32409	112.15



**Sigma - Aldrich Building Blocks**

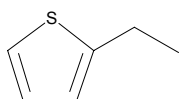
Brand	Catalog Number	MW
Aldrich	D188603	112.19





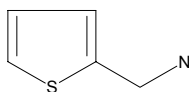
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	E49207	112.19



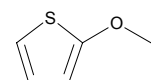
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	220884	113.18



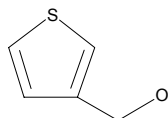
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	331597	114.17



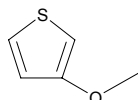
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	332399	114.17



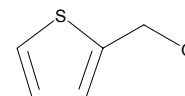
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	374024	114.17



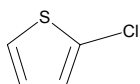
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	T32654	114.17



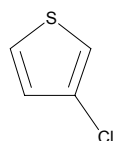
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	155667	118.59



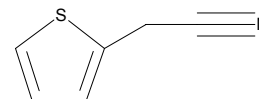
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	341061	118.59



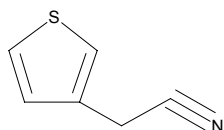
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	141682	123.18



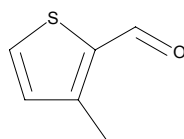
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	219193	123.18



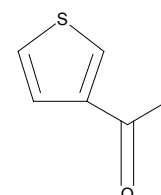
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	164135	126.18



**Sigma - Aldrich Building Blocks**

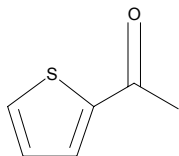
Brand	Catalog Number	MW
Aldrich	196320	126.18





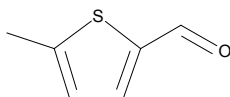
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	A22602	126.18



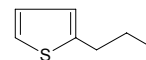
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	M84410	126.18



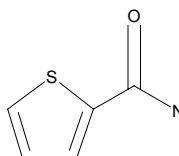
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	P54345	126.22



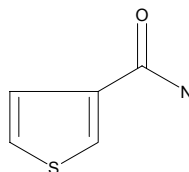
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	164453	127.17



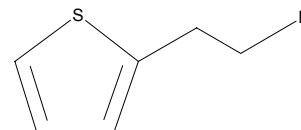
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	519502	127.17



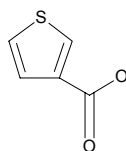
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	423270	127.21



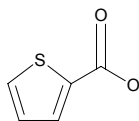
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	247766	128.15



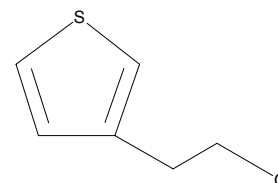
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	T32603	128.15



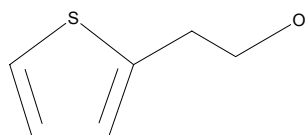
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	228796	128.19



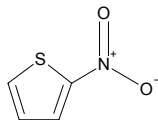
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	T27855	128.19



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	N27004	129.14



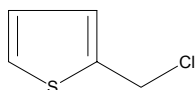
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	367141	131.04

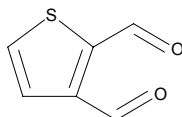


**Sigma - Aldrich Building Blocks**

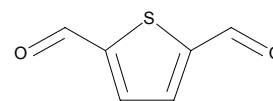
Brand	Catalog Number	MW
Aldrich	549282	132.61

**Sigma - Aldrich Building Blocks**

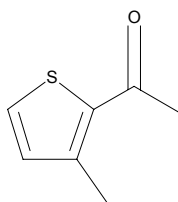
Brand	Catalog Number	MW
Aldrich	429872	140.16

**Sigma - Aldrich Building Blocks**

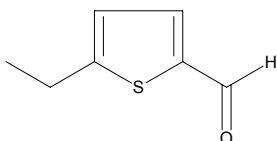
Brand	Catalog Number	MW
Aldrich	429880	140.16

**Sigma - Aldrich Building Blocks**

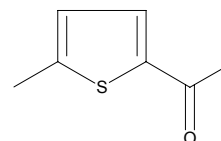
Brand	Catalog Number	MW
Aldrich	249645	140.21

**Sigma - Aldrich Building Blocks**

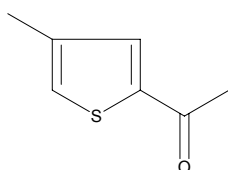
Brand	Catalog Number	MW
Aldrich	492809	140.21

**Sigma - Aldrich Building Blocks**

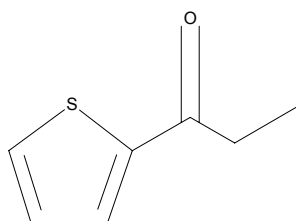
Brand	Catalog Number	MW
Aldrich	533610	140.21

**Sigma - Aldrich Building Blocks**

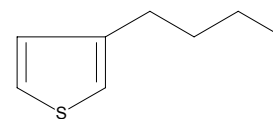
Brand	Catalog Number	MW
Aldrich	633046	140.21

**Sigma - Aldrich Building Blocks**

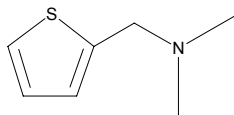
Brand	Catalog Number	MW
Aldrich	T27952	140.21

**Sigma - Aldrich Building Blocks**

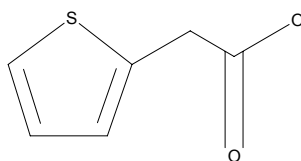
Brand	Catalog Number	MW
Aldrich	510424	140.25

**Sigma - Aldrich Building Blocks**

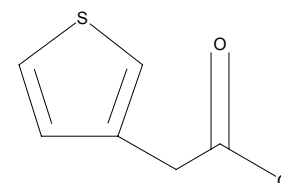
Brand	Catalog Number	MW
Aldrich	283460	141.24

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	195944	142.18

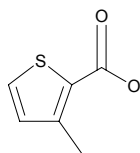
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	220639	142.18

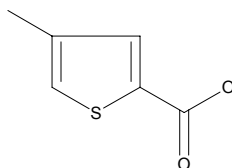


**Sigma - Aldrich Building Blocks**

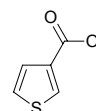
Brand	Catalog Number	MW
Aldrich	248207	142.18

**Sigma - Aldrich Building Blocks**

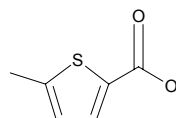
Brand	Catalog Number	MW
Aldrich	633550	142.18

**Sigma - Aldrich Building Blocks**

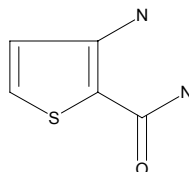
Brand	Catalog Number	MW
Aldrich	645559	142.18

**Sigma - Aldrich Building Blocks**

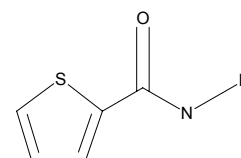
Brand	Catalog Number	MW
Aldrich	M84429	142.18

**Sigma - Aldrich Building Blocks**

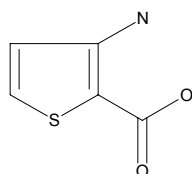
Brand	Catalog Number	MW
Aldrich	632813	142.18

**Sigma - Aldrich Building Blocks**

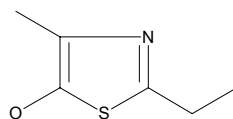
Brand	Catalog Number	MW
Aldrich	T32611	142.18

**Sigma - Aldrich Building Blocks**

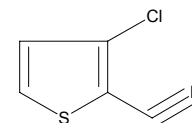
Brand	Catalog Number	MW
Aldrich	499099	143.17

**Sigma - Aldrich Building Blocks**

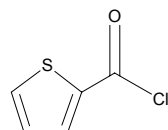
Brand	Catalog Number	MW
Aldrich	556726	143.21

**Sigma - Aldrich Building Blocks**

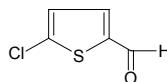
Brand	Catalog Number	MW
Aldrich	630047	143.60

**Sigma - Aldrich Building Blocks**

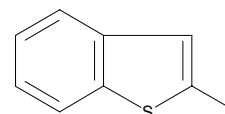
Brand	Catalog Number	MW
Aldrich	288985	146.60

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	443239	146.60

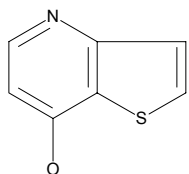
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	467456	148.23

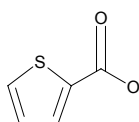


**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	498246	151.19

**Sigma - Aldrich Building Blocks**

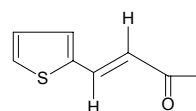
Brand	Catalog Number	MW
Aldrich	288977	152.15



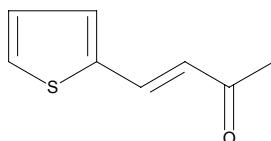
Na

**Sigma - Aldrich Building Blocks**

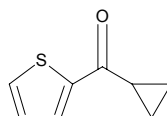
Brand	Catalog Number	MW
Aldrich	456225	152.22

**Sigma - Aldrich Building Blocks**

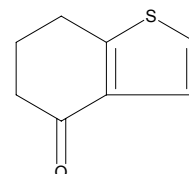
Brand	Catalog Number	MW
Aldrich	574457	152.22

**Sigma - Aldrich Building Blocks**

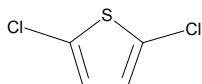
Brand	Catalog Number	MW
Aldrich	C121207	152.22

**Sigma - Aldrich Building Blocks**

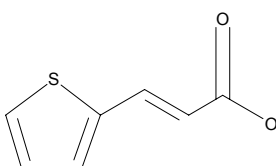
Brand	Catalog Number	MW
Aldrich	K3603	152.22

**Sigma - Aldrich Building Blocks**

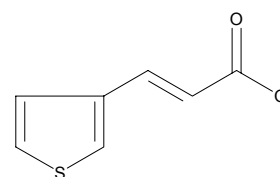
Brand	Catalog Number	MW
Aldrich	D75807	153.03

**Sigma - Aldrich Building Blocks**

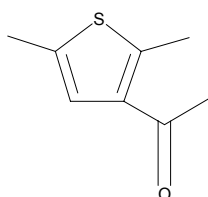
Brand	Catalog Number	MW
Aldrich	130583	154.19

**Sigma - Aldrich Building Blocks**

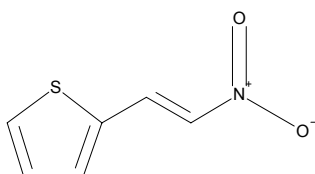
Brand	Catalog Number	MW
Aldrich	467987	154.19

**Sigma - Aldrich Building Blocks**

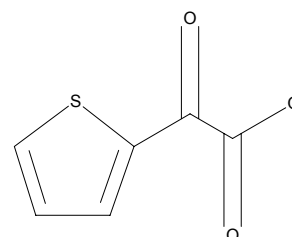
Brand	Catalog Number	MW
Aldrich	225797	154.23

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	387584	155.18

**Sigma - Aldrich Building Blocks**

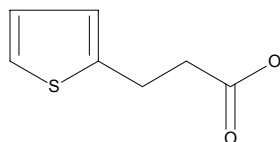
Brand	Catalog Number	MW
Aldrich	222275	156.16





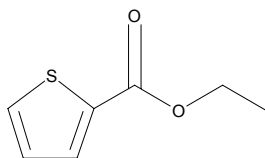
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	595535	156.20



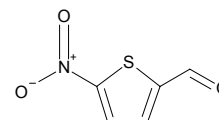
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	E49258	156.20



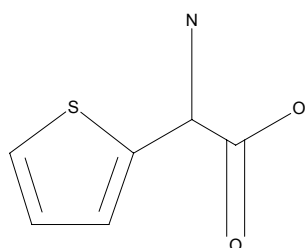
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	302295	157.15



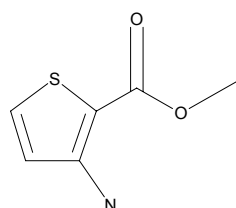
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	188808	157.19



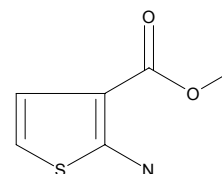
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	232904	157.19



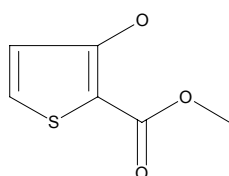
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	563080	157.19



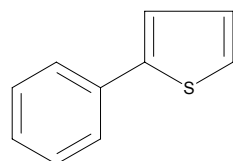
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	566977	158.18



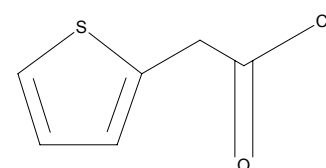
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	520578	160.24



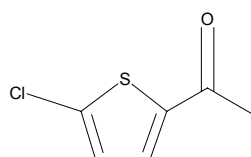
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	195995	160.62



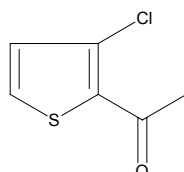
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	247073	160.62



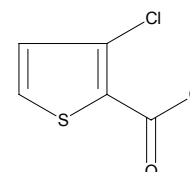
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	633038	160.62



**Sigma - Aldrich Building Blocks**

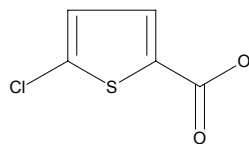
Brand	Catalog Number	MW
Aldrich	586994	162.60





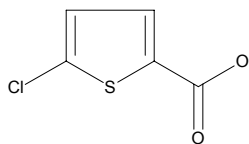
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	633003	162.60



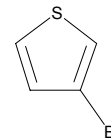
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	633003	162.60



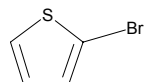
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	106224	163.04



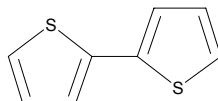
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	124168	163.04



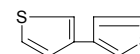
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	241636	166.27



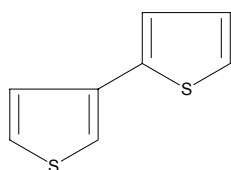
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	472913	166.27



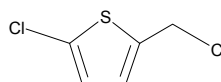
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	577812	166.27



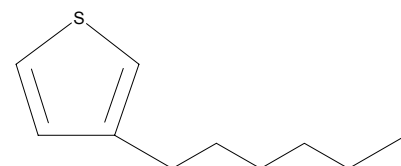
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	241822	167.06



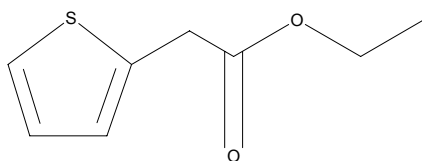
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	399051	168.30



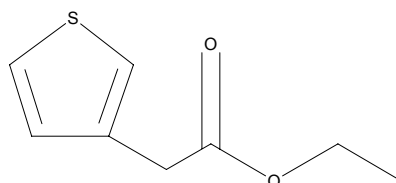
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	257648	170.23



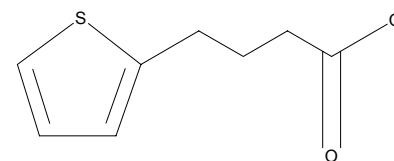
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	257656	170.23



**Sigma - Aldrich Building Blocks**

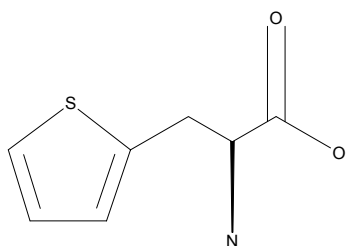
Brand	Catalog Number	MW
Aldrich	T27804	170.23





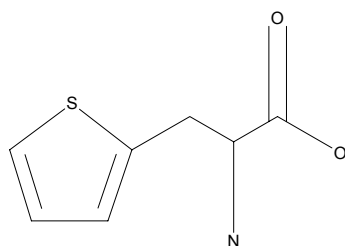
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	282154	171.22



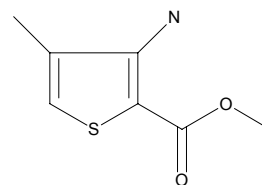
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	287288	171.22



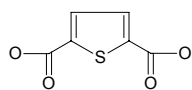
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	546658	171.22



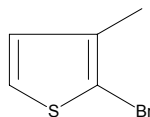
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	405191	172.16



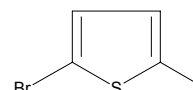
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	337021	177.06



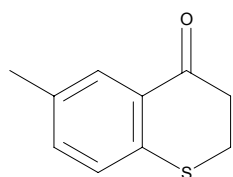
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	515485	177.06



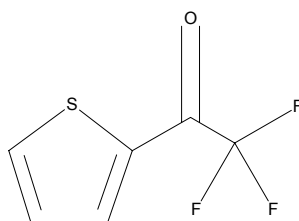
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	630683	178.25



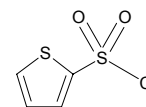
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	422215	180.15



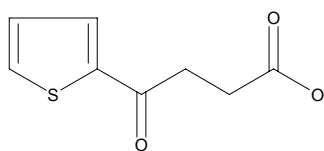
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	255335	182.65



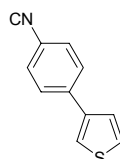
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	574406	184.22



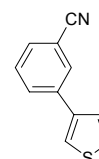
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	646520	185.25



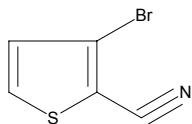
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	646539	185.25

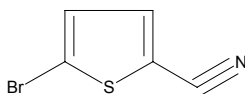


**Sigma - Aldrich Building Blocks**

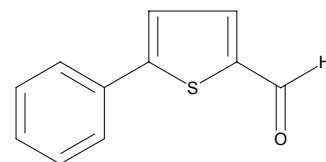
Brand	Catalog Number	MW
Aldrich	574392	188.05

**Sigma - Aldrich Building Blocks**

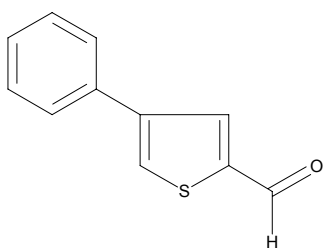
Brand	Catalog Number	MW
Aldrich	574430	188.05

**Sigma - Aldrich Building Blocks**

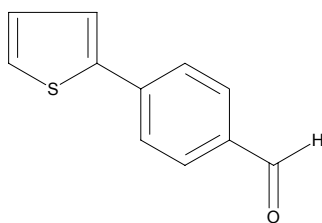
Brand	Catalog Number	MW
Aldrich	565938	188.25

**Sigma - Aldrich Building Blocks**

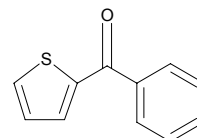
Brand	Catalog Number	MW
Aldrich	569526	188.25

**Sigma - Aldrich Building Blocks**

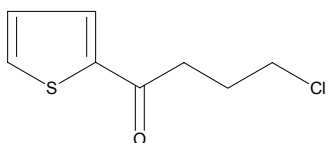
Brand	Catalog Number	MW
Aldrich	597325	188.25

**Sigma - Aldrich Building Blocks**

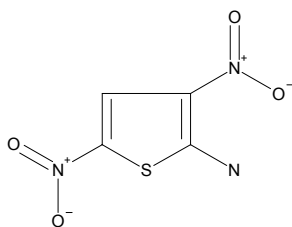
Brand	Catalog Number	MW
Aldrich	B14809	188.25

**Sigma - Aldrich Building Blocks**

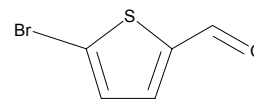
Brand	Catalog Number	MW
Aldrich	125016	188.68

**Sigma - Aldrich Building Blocks**

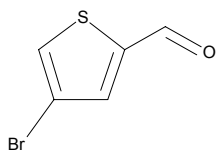
Brand	Catalog Number	MW
Aldrich	521620	189.15

**Sigma - Aldrich Building Blocks**

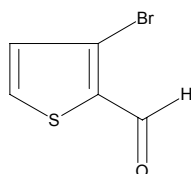
Brand	Catalog Number	MW
Aldrich	152625	191.05

**Sigma - Aldrich Building Blocks**

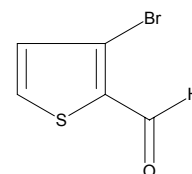
Brand	Catalog Number	MW
Aldrich	283452	191.05

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	529834	191.05

**Sigma - Aldrich Building Blocks**

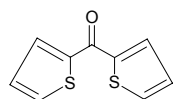
Brand	Catalog Number	MW
Aldrich	573484	191.05





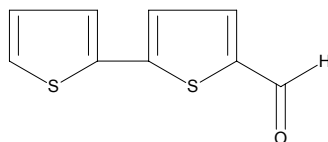
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	449377	194.28



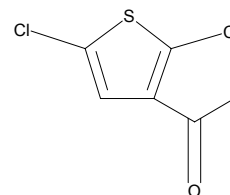
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	576700	194.28



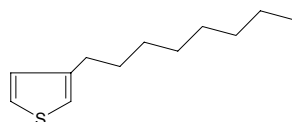
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	439509	195.07



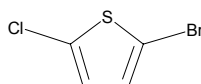
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	424285	196.36



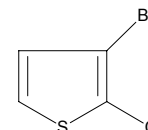
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	329762	197.48



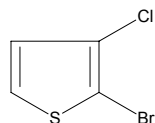
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	512389	197.48



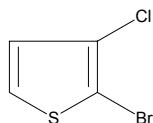
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	586978	197.48



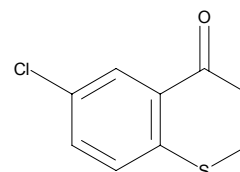
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	638188	197.48



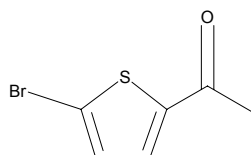
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	630691	198.67



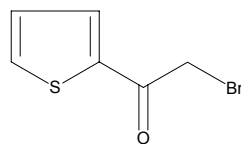
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	329738	205.07



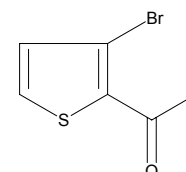
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	595594	205.07



**Sigma - Aldrich Building Blocks**

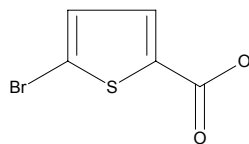
Brand	Catalog Number	MW
Aldrich	633054	205.07





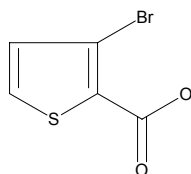
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	467944	207.05



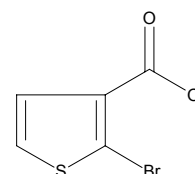
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	633011	207.05



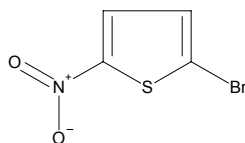
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	638129	207.05



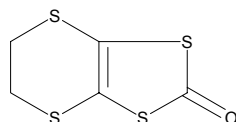
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	544531	208.03



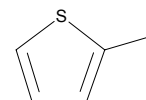
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	533467	208.34



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	196150	210.04



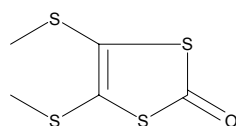
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	215023	210.04



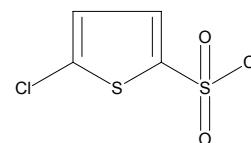
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	533440	210.36



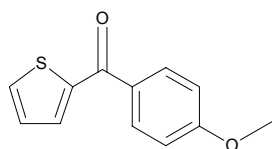
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	544272	217.09



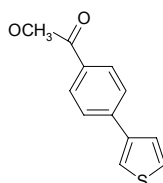
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	188336	218.28



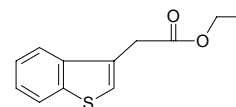
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	646512	218.28



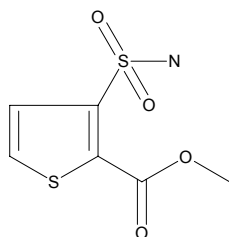
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	480843	220.29

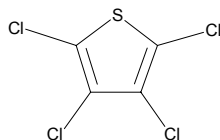


**Sigma - Aldrich Building Blocks**

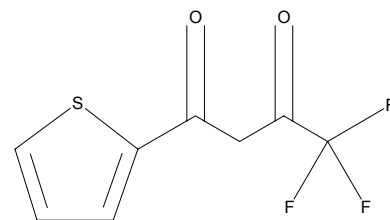
Brand	Catalog Number	MW
Aldrich	636657	221.25

**Sigma - Aldrich Building Blocks**

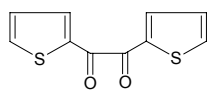
Brand	Catalog Number	MW
Aldrich	131873	221.92

**Sigma - Aldrich Building Blocks**

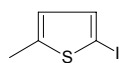
Brand	Catalog Number	MW
Aldrich	T27006	222.19

**Sigma - Aldrich Building Blocks**

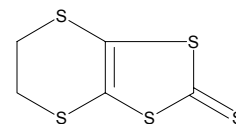
Brand	Catalog Number	MW
Aldrich	454389	222.29

**Sigma - Aldrich Building Blocks**

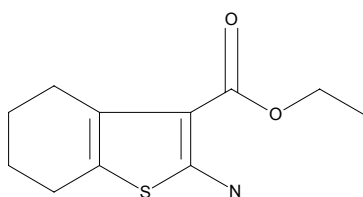
Brand	Catalog Number	MW
Aldrich	460532	224.06

**Sigma - Aldrich Building Blocks**

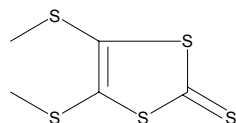
Brand	Catalog Number	MW
Aldrich	533459	224.41

**Sigma - Aldrich Building Blocks**

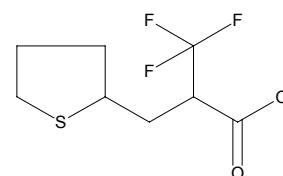
Brand	Catalog Number	MW
Aldrich	159182	225.31

**Sigma - Aldrich Building Blocks**

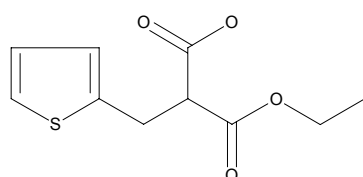
Brand	Catalog Number	MW
Aldrich	533432	226.42

**Sigma - Aldrich Building Blocks**

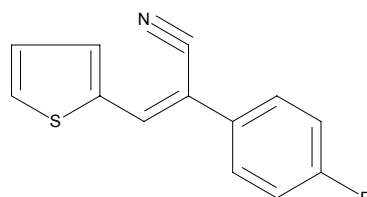
Brand	Catalog Number	MW
Aldrich	559628	228.23

**Sigma - Aldrich Building Blocks**

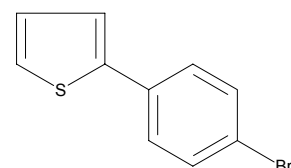
Brand	Catalog Number	MW
Aldrich	553573	228.27

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	596345	229.28

**Sigma - Aldrich Building Blocks**

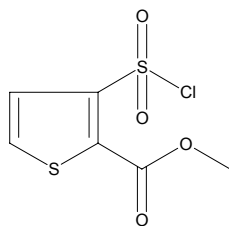
Brand	Catalog Number	MW
Aldrich	596663	239.14





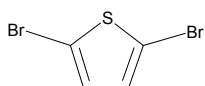
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	540501	240.69



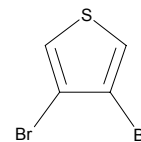
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	108472	241.93



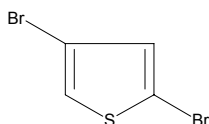
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	247154	241.93



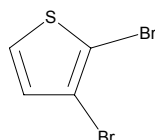
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	540749	241.93



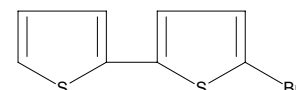
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	D43905	241.93



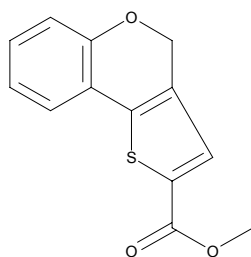
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	522856	245.16



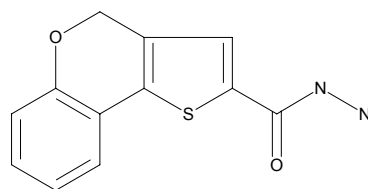
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	560898	246.29



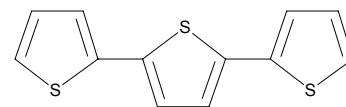
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	555746	246.29



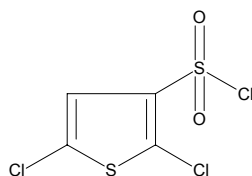
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	311073	248.39



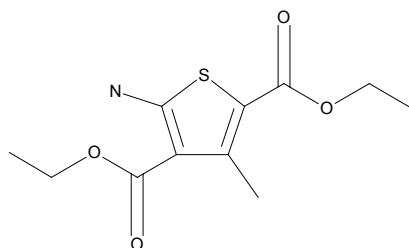
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	544280	251.54



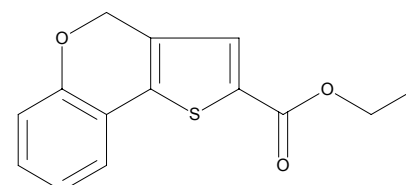
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	159174	257.31



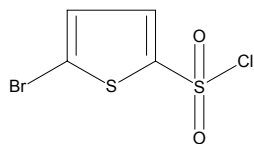
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	555703	260.31

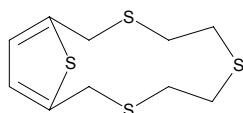


**Sigma - Aldrich Building Blocks**

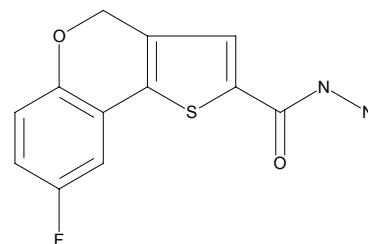
Brand	Catalog Number	MW
Aldrich	636223	261.54

**Sigma - Aldrich Building Blocks**

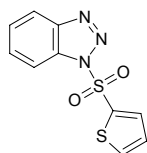
Brand	Catalog Number	MW
Aldrich	361402	262.48

**Sigma - Aldrich Building Blocks**

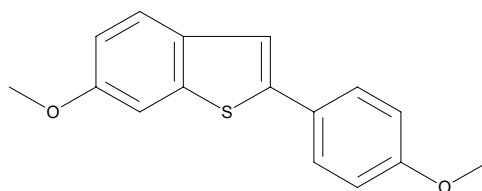
Brand	Catalog Number	MW
Aldrich	555754	264.28

**Sigma - Aldrich Building Blocks**

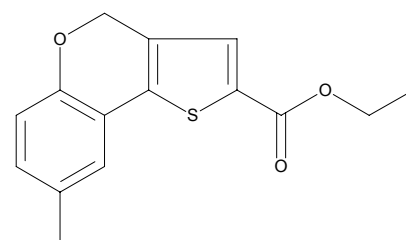
Brand	Catalog Number	MW
Aldrich	647357	265.31

**Sigma - Aldrich Building Blocks**

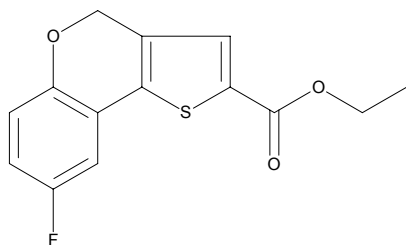
Brand	Catalog Number	MW
Aldrich	561290	270.35

**Sigma - Aldrich Building Blocks**

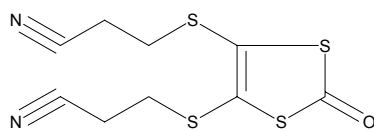
Brand	Catalog Number	MW
Aldrich	555711	274.34

**Sigma - Aldrich Building Blocks**

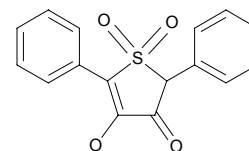
Brand	Catalog Number	MW
Aldrich	555738	278.30

**Sigma - Aldrich Building Blocks**

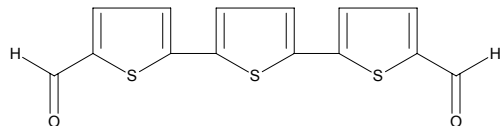
Brand	Catalog Number	MW
Aldrich	533424	288.43

**Sigma - Aldrich Building Blocks**

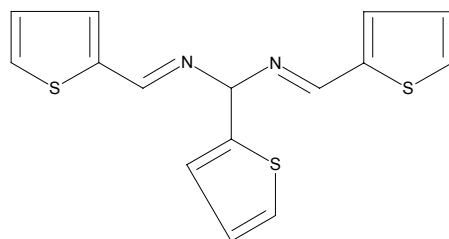
Brand	Catalog Number	MW
Aldrich	343579	300.34

**Sigma - Aldrich Building Blocks**

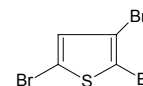
Brand	Catalog Number	MW
Aldrich	499102	304.41

**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	568457	316.47

**Sigma - Aldrich Building Blocks**

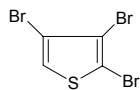
Brand	Catalog Number	MW
Aldrich	172758	320.83





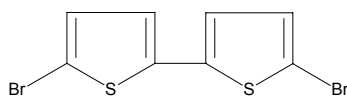
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	454753	320.83



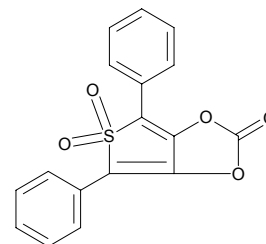
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	515493	324.06



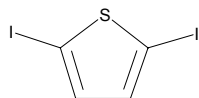
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	345903	326.33



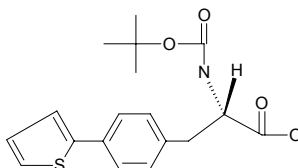
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	331570	335.93



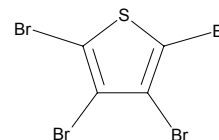
**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	464554	347.44



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	T5400	399.72



**Sigma - Aldrich Building Blocks**

Brand	Catalog Number	MW
Aldrich	447919	430.57

