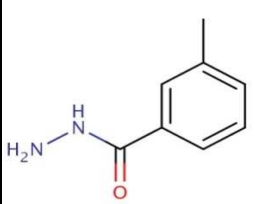
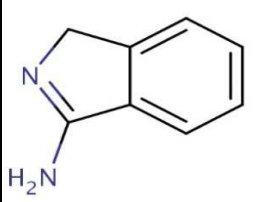

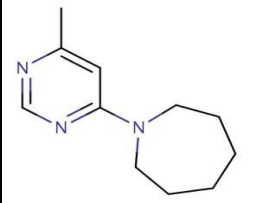
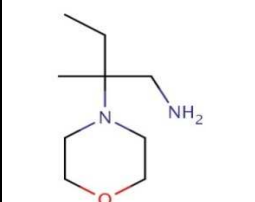
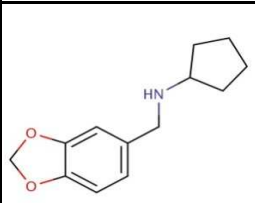
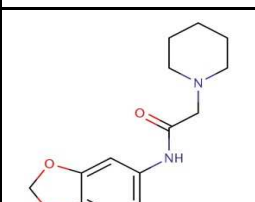
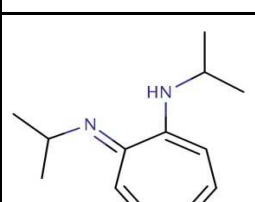


# Frag Xtal Screen

Fragment Screen for Crystallographic Screening

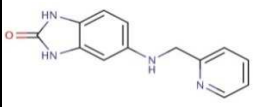
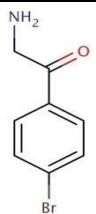
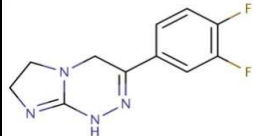
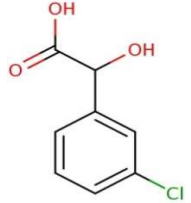
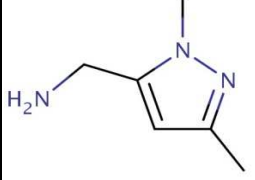

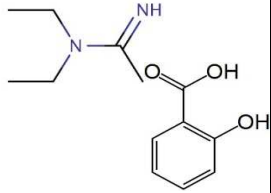
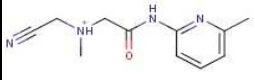
Cat.# X-FS-101

Plate ID	Fragment	Scifinder Name	CAS ID	MW
A1		Benzoic acid, 3-methyl-, hydrazide	13050-47-0	150,18
A2		1H-Isoindol-3-amine	76644-74-1	168,62
A3		Thiourea, N-(4-bromo-2-chlorophenyl)-N'-methyl-	838813-44-8	279,58
A4		1H-Azepine, hexahydro-1-(6-methyl-4-pyrimidinyl)-	1055196-50-3	227,74
A5		4-Morpholineethanamine, $\beta$ -ethyl- $\beta$ -methyl-	914206-26-1	172,30
A6		1,3-Benzodioxole-5-methanamine, N-cyclopentyl-	114413-77-3	219,30
A7		1-Piperidineacetamide, N-(1,3-benzodioxol-5-yl)-	487006-82-6	262,30
A8		1,3,5-Cycloheptatrien-1-amine, N-(1-methylethyl)-7-[(1-methylethyl)imino]-	3337-06-2	204,32

# Frag Xtal Screen

Fragment Screen for Crystallographic Screening

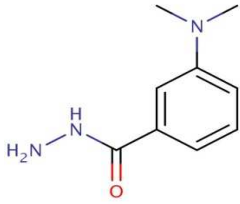
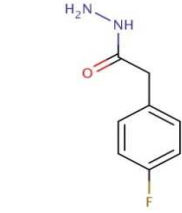
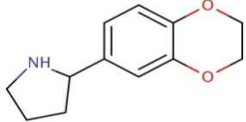
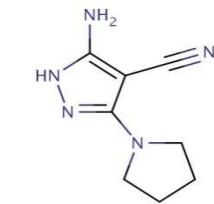
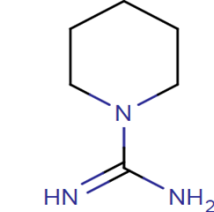
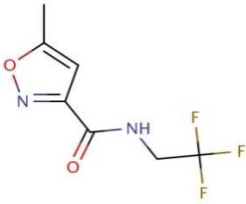
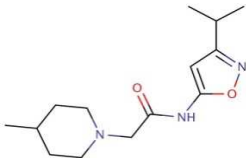
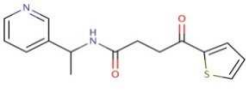
Cat.# X-FS-101

Plate ID	Fragment	Scifinder Name	CAS ID	MW
A9		2H -Benzimidazol-2-one, 1,3-dihydro-5-[(2-pyridinylmethyl)amino]-	876898-01-0	240,26
A10		Ethanone, 2-amino-1-(4-bromophenyl)-, hydrochloride (1:1)	5467-72-1	250,52
A11		Imidazo[2,1-c][1,2,4]triazine, 3-(3,4-difluorophenyl)-1,4,6,7-tetrahydro-	219929-40-5	236,22
A12		Benzeneacetic acid, 3-chloro-α-hydroxy-	61008-98-8	186,59
B1		1H -Pyrazole-5-methanamine, 1,3-dimethyl-	499770-63-7	125,20
B2		2-Furanmethanamine, N-methyl-5-(3-pyridinyloxy)-	859850-77-4	204,23
B3		Benzoic acid, 2-hydroxy-, compd. with N,N-diethylethanamide (1:1)	301843-27-6	252,30
B4		Acetamide, 2-[(cyanomethyl)methylamino]-N-(6-methyl-2-pyridinyl)-	1311649-76-9	218,25

# Frag Xtal Screen

Fragment Screen for Crystallographic Screening

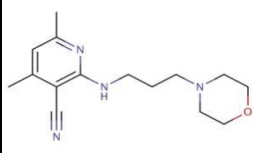
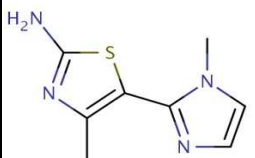

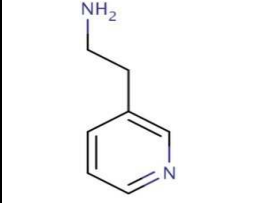

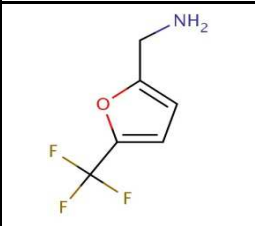
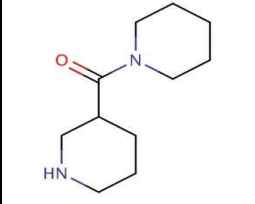
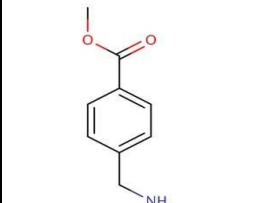
Cat.# X-FS-101

Plate ID	Fragment	Scifinder Name	CAS ID	MW
B5		Benzoic acid, 3-(dimethylamino)-, hydrazide	205927-64-6	179,22
B6		Benzeneacetic acid, 4-fluoro-, hydrazide	34547-28-9	168,17
B7		Pyrrolidine, 2-(2,3-dihydro-1,4-benzodioxin-6-yl)-	524674-08-6	205,25
B8		1H -Pyrazole-4-carbonitrile, 5-amino-3-(1-pyrrolidinyl)-	1119391-48-8	177,20
B9		1-Piperidinecarboximidamide, hydriodide (1:1)	102392-91-6	255,10
B10		3-Isoxazolecarboxamide, 5-methyl-N-(2,2,2-trifluoroethyl)-	878595-52-9	208,13
B11		1-Piperidineacetamide, 4-methyl-N-[3-(1-methylethyl)-5-isoxazolyl]-	923192-35-2	265,35
B12		2-Thiophenebutanamide, $\gamma$ -oxo-N-[1-(3-pyridinyl)ethyl]-	923810-83-7	288,36

# Frag Xtal Screen

Fragment Screen for Crystallographic Screening

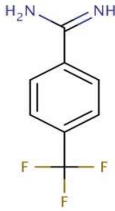
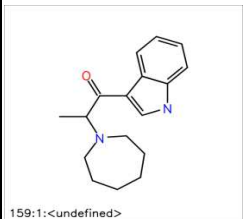
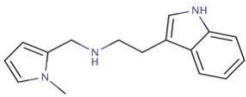
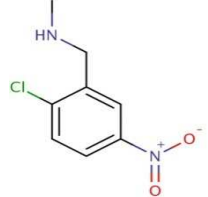
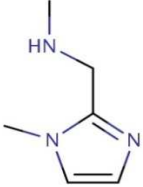
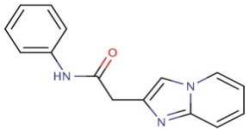
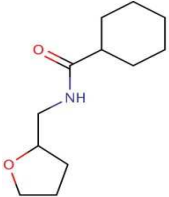
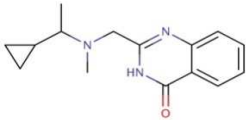
Cat.# X-FS-101

Plate ID	Fragment	Scifinder Name	CAS ID	MW
C1		3-Pyridinecarbonitrile, 4,6-dimethyl-2-[[3-(4-morpholinyl)propyl]amino]-	831183-48-3	274,36
C2		2-Thiazolamine, 4-methyl-5-(1-methyl-1H-imidazol-2-yl)-	929975-79-1	194,26
C3	 <small>337:1:&lt;undefined&gt;</small>	1,3,5-Triazine-2,4-diamine, 6-[1-(hexahydro-1H-azepin-1-yl)ethyl]-N2,N2-dimethyl-	923208-22-4	264,37
C4		3-Pyridineethanamine	20173-24-4	122,17
C5		3-Thiophenecarboximidamide, hydrochloride (1:1)	51528-02-0	162,64
C6		2-Furanmethanamine, 5-(trifluoromethyl)-	868755-68-4	165,11
C7		Methanone, 1-piperidiny-3-piperidiny-	845885-83-8	232,75
C8		Benzoic acid, 4-(aminomethyl)-, methyl ester, hydrochloride (1:1)	6232-11-7	201,65

# Frag Xtal Screen

Fragment Screen for Crystallographic Screening

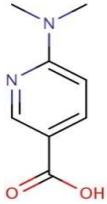
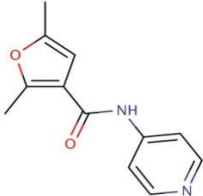
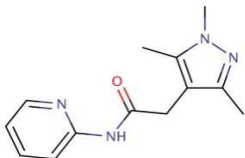
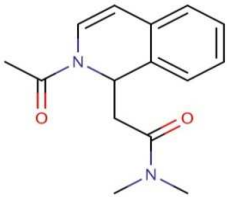
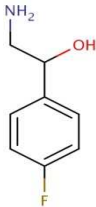
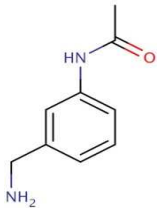
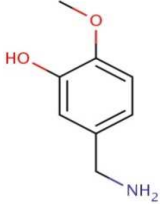

Cat.# X-FS-101

Plate ID	Fragment	Scifinder Name	CAS ID	MW
C9		Benzenecarboximidamide, 4-(trifluoromethyl)-	38980-96-0	224,61
C10	 <small>159:1:&lt;undefined&gt;</small>	Guanidine, N-(4-ethoxy-8-methyl-2-quinazolinyl)-	923803-92-3	270,37
C11		1H-Indole-3-ethanamine, N-[(1-methyl-1H-pyrrol-2-yl)methyl]-	289487-79-2	253,34
C12		Benzenemethanamine, 2-chloro-N-methyl-5-nitro-	1177318-36-3	237,08
D1		1H-Imidazole-2-methanamine, N,1-dimethyl-	1049717-38-5	198,09
D2		Imidazo[1,2-a]pyridine-2-acetamide, N-phenyl-	1090352-33-2	251,28
D3		Cyclohexanecarboxamide, N-[(tetrahydro-2-furanyl)methyl]-	401597-32-8	211,30
D4		4(3H)-Quinazolinone, 2-[[[(1-cyclopropylethyl)methylamino]methyl]-	1090518-66-3	257,33

# Frag Xtal Screen

Fragment Screen for Crystallographic Screening

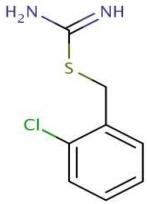
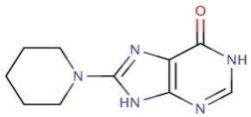
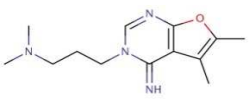
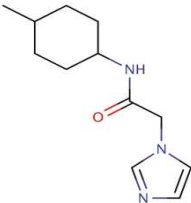
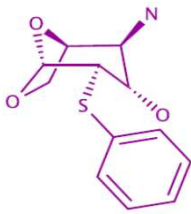
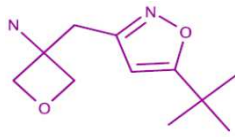
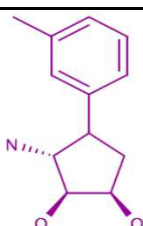
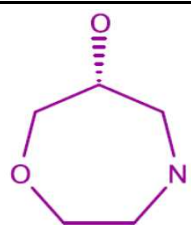
Cat.# X-FS-101

Plate ID	Fragment	Scifinder Name	CAS ID	MW
D5		3-Pyridinecarboxylic acid, 6-(dimethylamino)-	82846-28-4	166,18
D6		3-Furancarboxamide, 2,5-dimethyl-N-4-pyridinyl-	501442-73-5	216,24
D7		1H-Pyrazole-4-acetamide, 1,3,5-trimethyl-N-2-pyridinyl-	1171575-61-3	244,29
D8		1-Isoquinolineacetamide, 2-acetyl-1,2-dihydro-N,N-dimethyl-	1090496-96-0	258,32
D9		Benzenemethanol, $\alpha$ -(aminomethyl)-4-fluoro-	456-05-3	155,17
D10		Acetamide, N-[3-(aminomethyl)phenyl]-	96783-68-5	164,20
D11		Phenol, 5-(aminomethyl)-2-methoxy-, hydrochloride (1:1)	89702-89-6	153,18
D12		4-Pyridinemethanamine, N-(2,3-dihydro-1,4-benzodioxin-6-yl)-	774553-49-0	242,27

# Frag Xtal Screen

Fragment Screen for Crystallographic Screening

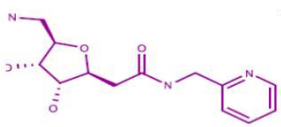
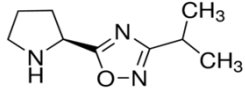
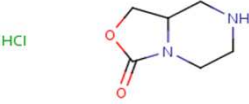
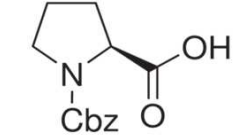
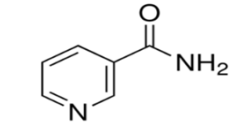
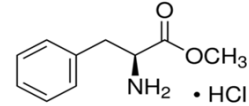
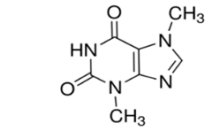
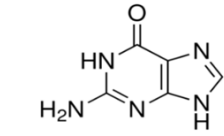
Cat.# X-FS-101

Plate ID	Fragment	Scifinder Name	CAS ID	MW
E1		Carbamimidothioic acid, (2-chlorophenyl)methyl ester	14122-38-4	200,69
E2		6H-Purin-6-one, 1,9-dihydro-8-(1-piperidinyl)-	312923-57-2	219,25
E3		Furo[2,3-d]pyrimidine-3(4H)-propanamine, 4-imino-N,N,5,6-tetramethyl-	929868-45-1	248,33
E4		1H-Imidazole-1-acetamide, N-(4-methylcyclohexyl)-	1118862-09-1	221,30
E5		INDEX NAME NOT YET ASSIGNED	1212574-75-8	253,32
E6		3-Oxetanamine, 3-[[5-(1,1-dimethylethyl)-3-isoxazolyl]methyl]-	1371533-72-0	210,28
E7		1,2-Cyclopentanediol, 3-amino-4-(3-methylphenyl)-, (1R,2S,3R)-	1421058-66-3	243,74
E8		1,4-Oxazepin-6-ol, hexahydro-, (6R)-	1022915-33-8	153,65

# Frag Xtal Screen

Fragment Screen for Crystallographic Screening

Cat.# X-FS-101

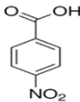
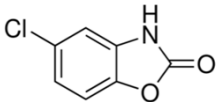
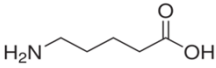
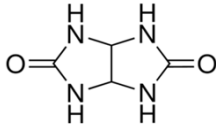
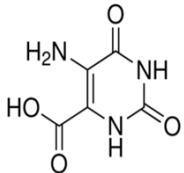
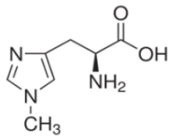
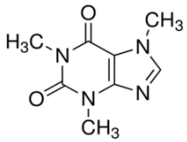
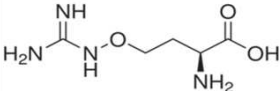
Plate ID	Fragment	Scifinder Name	CAS ID	MW
E9		NO NAME ASSIGNED IN SCIFINDER	NO CAS AVAILABLE	354,24
E10		1,2,4-Oxadiazole, 3-(1-methylethyl)-5-(2-pyrrolidinyl)-	1185170-55-1	126,16
E11	HCl 	3H -Oxazolo[3,4-a]pyrazin-3-one, hexahydro-	1020349-31-8	178,61
E12		1,2-Pyrrolidinedicarboxylic acid, 1-(phenylmethyl) ester, (2S)-	1148-11-4	249,26
F1		3-Pyridinecarboxamide	98-92-0	122,12
F2		L-Phenylalanine, methyl ester, hydrochloride (1:1)	7524-50-7	179,22
F3		1H-Purine-2,6-dione, 3,7-dihydro-3,7-dimethyl-	83-67-0	180,16
F4		6H-Purin-6-one, 2-amino-1,9-dihydro-	73-40-5	151,13



# Frag Xtal Screen

Fragment Screen for Crystallographic Screening

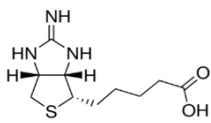
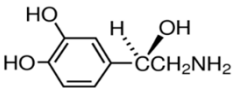
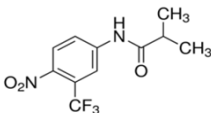
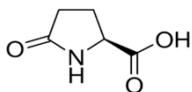
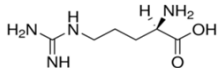
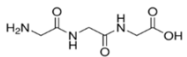
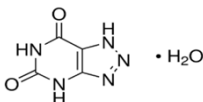
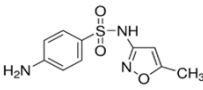
Cat.# X-FS-101

Plate ID	Fragment	Scifinder Name	CAS ID	MW
F5		Benzoic acid, 4-nitro-	62-23-7	167,12
F6		2(3H)-Benzoxazolone, 5-chloro-	95-25-0	169,57
F7		Pentanoic acid, 5-amino-	660-88-8	117,15
F8		Imidazo[4,5-d]imidazole-2,5(1H,3H)-dione, tetrahydro-	496-46-8	142,12
F9		4-Pyrimidinecarboxylic acid, 5-amino-1,2,3,6-tetrahydro-2,6-dioxo-	7164-43-4	171,11
F10		L-Histidine, 1-methyl-	332-80-9	169,18
F11		1H-Purine-2,6-dione, 3,7-dihydro-1,3,7-trimethyl-	58-08-2	194,19
F12		L-Homoserine, O-[(aminoiminomethyl)amino]-	543-38-4	176,17

# Frag Xtal Screen

Fragment Screen for Crystallographic Screening

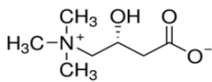
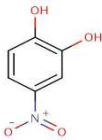
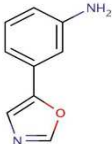
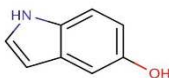
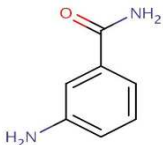
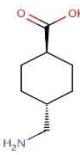
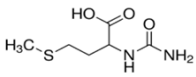
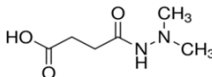
Cat.# X-FS-101

Plate ID	Fragment	Scifinder Name	CAS ID	MW
G1		1H-Thieno[3,4-d]imidazole-6-pentanoic acid, 2-amino-3a,4,6,6a-tetrahydro-, (3aR,6S,6aS)-	13395-35-2	243,33
G2		1,2-Benzenediol, 4-[(1R)-2-amino-1-hydroxyethyl]-	51-41-2	169,18
G3		Propanamide, 2-methyl-N-[4-nitro-3-(trifluoromethyl)phenyl]-	13311-84-7	276,21
G4		L-Proline, 5-oxo-	98-79-3	129,11
G5		D-Arginine	157-06-2	174,20
G6		Glycine, glycylglycyl-	556-33-2	189,17
G7		3H-1,2,3-Triazolo[4,5-d]pyrimidine-5,7(4H,6H)-dione	1468-26-4	171,11
G8		Benzenesulfonamide, 4-amino-N-(5-methyl-3-isoxazolyl)-	723-46-6	253,28

# Frag Xtal Screen

Fragment Screen for Crystallographic Screening

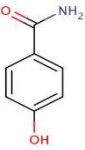
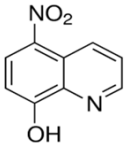
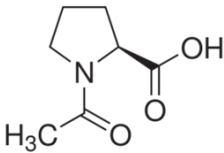
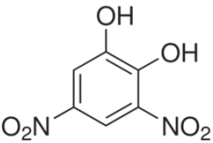
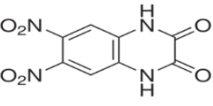
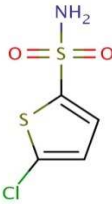
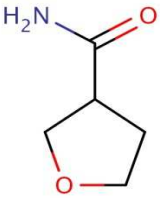
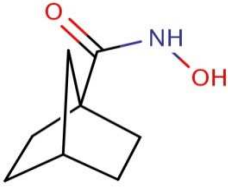
Cat.# X-FS-101

Plate ID	Fragment	Scifinder Name	CAS ID	MW
G9		1-Propanaminium, 3-carboxy-2-hydroxy-N,N,N-trimethyl-, inner salt, (2R)-	541-15-1	161,20
G10		1,2-Benzenediol, 4-nitro-	3316-09-4	155,11
G11		Benzenamine, 3-(5-oxazolyl)-	157837-31-5	160,17
G12		1H-Indol-5-ol	1953-54-4	133,15
H1		Benzamide, 3-amino-	3544-24-9	136,15
H2		Cyclohexanecarboxylic acid, 4-(aminomethyl)-, trans-	1197-18-8	157,21
H3		Methionine, N-(aminocarbonyl)-	30411-84-8	192,24
H4		Butanedioic acid, 1-(2,2-dimethylhydrazide)	1596-84-5	160,17

# Frag Xtal Screen

Fragment Screen for Crystallographic Screening

Cat.# X-FS-101

Plate ID	Fragment	Scifinder Name	CAS ID	MW
H5		Benzamide, 4-hydroxy-	619-57-8	137,14
H6		8-Quinolinol, 5-nitro-	4008-48-4	190,16
H7		L-Proline, 1-acetyl-	68-95-1	157,17
H8		1,2-Benzenediol, 3,5-dinitro-	7659-29-2	200,11
H9		2,3-Quinoxalinedione, 1,4-dihydro-6,7-dinitro-	2379-57-9	252,14
H10		2-Thiophenesulfonamide, 5-chloro-	53595-66-7	197,66
H11		3-Furancarboxamide, tetrahydro-	871677-92-8	115,13
H12		Bicyclo[2.2.1]heptane-1-carboxamide, N-hydroxy-	256954-29-7	155,20