

Compounds Included in the Fentanyl Analog Screening Kit



The Fentanyl Analog Screening (FAS) Kit and its expansion pack (Emergent Panel Version 1) contain analog reference materials that will enable labs to screen for 150 synthetic opioid compounds, including more than 140 fentanyl analogs.

No.	Compound Name
1	<i>para</i> -methoxy Butyryl fentanyl (hydrochloride)
2	Acetyl norfentanyl (hydrochloride)
3	α -methyl Acetyl fentanyl (hydrochloride)
4	Crotonyl fentanyl
5	Fentanyl
6	Acrylfentanyl (hydrochloride)
7	Cyclopropyl fentanyl (hydrochloride)
8	Isobutyryl fentanyl (hydrochloride)
9	Methoxyacetyl fentanyl (hydrochloride)
10	Valeryl fentanyl (hydrochloride)
11	(\pm)- <i>cis</i> -3-methyl Fentanyl (hydrochloride)
12	2'-fluoro <i>ortho</i> -Fluorofentanyl (hydrochloride)
13	<i>ortho</i> -Fluorobutyryl fentanyl (hydrochloride)
14	<i>ortho</i> -Fluorofentanyl (hydrochloride)
15	<i>ortho</i> -isopropyl Furanyl fentanyl
16	<i>ortho</i> -methyl Furanyl fentanyl (hydrochloride)
17	<i>ortho</i> -methyl Furanyl fentanyl
18	<i>meta</i> -Fluorobutyryl fentanyl (hydrochloride)
19	(\pm)- <i>cis</i> -3-methyl Butyryl fentanyl (hydrochloride)
20	4'-methyl Acetyl fentanyl (hydrochloride)
21	<i>para</i> -chloro Furanyl fentanyl
22	<i>para</i> -Methoxyfentanyl (hydrochloride)
23	<i>para</i> -methoxy Furanyl fentanyl (hydrochloride)
24	α -methyl Butyryl fentanyl (hydrochloride)
25	α -methyl Thiofentanyl (hydrochloride)
26	Benzodioxole fentanyl
27	Phenyl fentanyl (hydrochloride)
28	Benzyl fentanyl (hydrochloride)
29	N-benzyl Furanyl norfentanyl (hydrochloride)
30	Cyclohexyl fentanyl (hydrochloride)
31	Cyclopentyl fentanyl (hydrochloride)
32	<i>meta</i> -Methylfentanyl (hydrochloride)
33	N-methyl Norcarfentanil (hydrochloride)

No.	Compound Name
34	Norcarfentanil (hydrochloride)
35	<i>ortho</i> -methyl Acrylfentanyl (hydrochloride)
36	<i>ortho</i> -methyl Methoxyacetyl fentanyl (hydrochloride)
37	<i>para</i> -Chloroisobutyryl fentanyl (hydrochloride)
38	Tetrahydrofuran fentanyl (hydrochloride)
39	2,2,3,3-tetramethyl-Cyclopropyl fentanyl (hydrochloride)
40	Thienyl fentanyl (hydrochloride)
41	Thiophene fentanyl (hydrochloride)
42	Thiofentanyl (hydrochloride)
43	<i>ortho</i> -methyl Acetyl fentanyl (hydrochloride)
44	α' -methoxy Fentanyl (hydrochloride)
45	α' -methyl Butyryl fentanyl (hydrochloride)
46	Cyclobutyl fentanyl (hydrochloride)
47	<i>meta</i> -Fluorofentanyl (hydrochloride)
48	N-methyl Cyclopropyl norfentanyl (hydrochloride)
49	(\pm)- <i>trans</i> -3-methyl Thiofentanyl (hydrochloride)
50	4'-fluoro, <i>para</i> -fluoro (\pm)- <i>trans</i> -3-methyl Fentanyl (hydrochloride)
51	4'-methyl Fentanyl (hydrochloride)
52	Alfentanil (hydrochloride)
53	Benzyl Carfentanil (hydrochloride)
54	Benzyl Acrylfentanyl (hydrochloride)
55	β' -Phenyl fentanyl
56	β -methyl Acetyl fentanyl (hydrochloride)
57	Ethoxyacetyl fentanyl (hydrochloride)
58	Fentanyl Carbamate
59	Furanyl fentanyl 3-furancarboxamide isomer (hydrochloride)
60	Furanylethyl fentanyl (hydrochloride)
61	Isovaleryl fentanyl (hydrochloride)
62	<i>meta</i> -Fluoroisobutyryl fentanyl (hydrochloride)
63	<i>meta</i> -fluoro Methoxyacetyl fentanyl (hydrochloride)
64	<i>meta</i> -methyl Furanyl fentanyl (hydrochloride)
65	<i>meta</i> -methyl Methoxyacetyl fentanyl (hydrochloride)
66	Methacrylfentanyl

Compounds Included in the Fentanyl Analog Screening Kit *Continued*

No.	Compound Name
67	<i>ortho</i> -fluoro Acrylfentanyl (hydrochloride)
68	<i>ortho</i> -fluoro Furanyl fentanyl (hydrochloride)
69	<i>ortho</i> -Fluoroisobutyryl fentanyl (hydrochloride)
70	<i>ortho</i> -methoxy Butyryl fentanyl (hydrochloride)
71	<i>ortho</i> -methyl Cyclopropyl fentanyl (hydrochloride)
72	<i>ortho</i> -Methylfentanyl (hydrochloride)
73	<i>ortho</i> -methyl Phenyl fentanyl (hydrochloride)
74	<i>para</i> -Chlorofentanyl (hydrochloride)
75	<i>para</i> -chloro Acrylfentanyl (hydrochloride)
76	<i>para</i> -Chlorobutyryl fentanyl (hydrochloride)
77	<i>para</i> -chloro Cyclobutyl fentanyl (hydrochloride)
78	<i>para</i> -chloro Cyclopentyl fentanyl (hydrochloride)
79	<i>para</i> -chloro Methoxyacetyl fentanyl (hydrochloride)
80	<i>para</i> -chloro Valeryl fentanyl (hydrochloride)
81	<i>para</i> -chloro Cyclopropyl fentanyl (hydrochloride)
82	<i>para</i> -fluoro Acrylfentanyl
83	<i>para</i> -fluoro Crotonyl fentanyl
84	<i>para</i> -fluoro Cyclopentyl fentanyl (hydrochloride)
85	<i>para</i> -fluoro Cyclopropyl fentanyl (hydrochloride)
86	<i>para</i> -fluoro Furanyl fentanyl 3-furancarboxamide (hydrochloride)
87	<i>para</i> -fluoro Furanyl fentanyl (hydrochloride)
88	<i>para</i> -fluoro Methoxyacetyl fentanyl (hydrochloride)
89	<i>para</i> -fluoro Tetrahydrofuran fentanyl (hydrochloride)
90	<i>para</i> -fluoro Valeryl fentanyl (hydrochloride)
91	<i>para</i> -Fluoroacetyl fentanyl (hydrochloride)
92	<i>para</i> -methoxy Acrylfentanyl (hydrochloride)
93	<i>para</i> -methoxy Tetrahydrofuran fentanyl (hydrochloride)
94	<i>para</i> -methoxy Valeryl fentanyl (hydrochloride)
95	<i>para</i> -Methylfentanyl (hydrochloride)
96	<i>para</i> -methyl Acetyl fentanyl (hydrochloride)
97	<i>para</i> -methyl Acrylfentanyl (hydrochloride)
98	<i>para</i> -methyl Cyclopropyl fentanyl (hydrochloride)
99	<i>para</i> -methyl Furanyl fentanyl (hydrochloride)
100	<i>para</i> -methyl Isobutyryl fentanyl (hydrochloride)
101	<i>para</i> -methyl Methoxyacetyl fentanyl (hydrochloride)
102	<i>para</i> -methyl Tetrahydrofuran fentanyl (hydrochloride)
103	Phenylacetyl fentanyl (hydrochloride)

No.	Compound Name
104	Pivaloyl fentanyl (hydrochloride)
105	Remifentanil (hydrochloride)
106	Sufentanil
107	Tetrahydrofuran fentanyl 3-tetrahydrofurancarboxamide (hydrochloride)
108	(±)- <i>cis</i> -3-methyl Thiofentanyl (hydrochloride)
109	β-Hydroxythiofentanyl (hydrochloride)
110	Norfentanyl
111	4-ANPP
112	β-methyl Fentanyl (hydrochloride)
113	(±)- <i>trans</i> -3-methyl Fentanyl (hydrochloride)
114	<i>para</i> -Fluorobutyryl fentanyl (hydrochloride)
115	FIBF (hydrochloride)
116	α-methyl Fentanyl (hydrochloride)
117	Ocfentanil (hydrochloride)
118	<i>para</i> -Fluorofentanyl (hydrochloride)
119	Butyryl norfentanyl (hydrochloride)
120	Butyryl fentanyl (hydrochloride)
121	Furanyl fentanyl (hydrochloride)
122	Acetyl fentanyl (hydrochloride)
123	4'-Fluorofentanyl (hydrochloride)
124	4-Phenyl fentanyl (hydrochloride)
125	Cyclopentenyl fentanyl (hydrochloride)
126	<i>para</i> -methyl Cyclopentyl fentanyl (hydrochloride)
127	<i>meta</i> -methyl Cyclopropyl fentanyl (hydrochloride)
128	<i>para</i> -methyl Butyryl fentanyl (hydrochloride)
129	<i>para</i> -methoxy Acetyl fentanyl (hydrochloride)
130	<i>para</i> -methoxy Methoxyacetyl fentanyl (hydrochloride)
131	Tetrahydrothiophene fentanyl
132	N,N-Dimethylamido-despropionyl fentanyl
133	Hexanoyl fentanyl (hydrochloride)
134	Heptanoyl fentanyl (hydrochloride)
135	β-Hydroxythioacetylfentanyl
136	2,3-seco-Fentanyl (hydrochloride)
137	Senecioylfentanyl
138	Phenoxyacetyl fentanyl (hydrochloride)
139	Fentanyl Methyl Carbamate
140	β-hydroxy Fentanyl (hydrochloride)

Compounds Included in the Fentanyl Analog Screening Kit *Continued*

No.	Compound Name
141	N-(3-ethylindole) Norfentanyl
142	U-47700
143	U-49900
144	U-48800 (hydrochloride)
145	Isopropyl U-47700
146	AH 7921
147	MT-45 (hydrochloride)
148	2-fluoro MT-45 (hydrochloride)
149	Furanyl norfentanyl (hydrochloride)
150	Norsufentanil