

Compounds Included in the Fentanyl Analog Screening Kit



The Fentanyl Analog Screening (FAS) Kit and its expansion packs (Emergent Panels Version 1, 2, and 3) contain analog reference materials that enable labs to screen for 212 synthetic opioid compounds, including more than 190 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> -methoxy Furanyl fentanyl
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> -Fluoroisobutyl fentanyl (hydrochloride)
70	<i>ortho</i> -methoxy Butyl 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> -Chlorobutyl 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 Isobutyl 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	Remifentanyl (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> -Fluorobutyl fentanyl (hydrochloride)
115	FIBF (hydrochloride)
116	α-methyl Fentanyl (hydrochloride)
117	Ocfentanyl (hydrochloride)
118	<i>para</i> -Fluorofentanyl (hydrochloride)
119	Butyl norfentanyl (hydrochloride)
120	Butyl fentanyl (hydrochloride)
121	Furanyl fentanyl (hydrochloride)
122	Norsufentanil
123	Acetyl fentanyl (hydrochloride)
124	4'-Fluorofentanyl (hydrochloride)
125	4-Phenyl fentanyl (hydrochloride)
126	Cyclopentenyl fentanyl (hydrochloride)
127	<i>para</i> -methyl Cyclopentyl fentanyl (hydrochloride)
128	<i>meta</i> -methyl Cyclopropyl fentanyl (hydrochloride)
129	<i>para</i> -methyl Butyl fentanyl (hydrochloride)
130	<i>para</i> -methoxy Acetyl fentanyl (hydrochloride)
131	<i>para</i> -methoxy Methoxyacetyl fentanyl (hydrochloride)
132	Tetrahydrothiophene fentanyl
133	N,N-Dimethylamido-despropionyl fentanyl
134	Hexanoyl fentanyl (hydrochloride)
135	Heptanoyl fentanyl (hydrochloride)
136	β-Hydroxythioacetyl fentanyl
137	2,3-seco-Fentanyl (hydrochloride)
138	Seneciylfentanyl
139	Phenoxyacetyl fentanyl (hydrochloride)
140	Fentanyl Methyl Carbamate

Compounds Included in the Fentanyl Analog Screening Kit *Continued*

No.	Compound Name
141	β -hydroxy Fentanyl (hydrochloride)
142	N-(3-ethylindole) Norfentanyl
143	U-47700
144	U-49900
145	U-48800 (hydrochloride)
146	Isopropyl U-47700
147	AH 7921
148	MT-45 (hydrochloride)
149	2-fluoro MT-45 (hydrochloride)
150	Furanyl norfentanyl (hydrochloride)
151	U-48520
152	U-50488 (hydrochloride)
153	U-51754 (hydrochloride)
154	U-47931E
155	U-48753E (maleate)
156	<i>meta</i> -fluoro Furanyl fentanyl (hydrochloride)
157	3,4-Methylenedioxy U-47700 (hydrochloride)
158	3,4-Ethylenedioxy U-51754 (hydrochloride)
159	3,4-Ethylenedioxy U-47700 (hydrochloride)
160	N-methyl U-47931E
161	Propyl U-47700
162	4-phenyl U-51754 (hydrochloride)
163	N-desmethyl U-47700
164	N,N-didesmethyl U-47700 (trifluoroacetate salt)
165	3'-methyl Acetyl fentanyl (hydrochloride)
166	2'-methyl Acetyl fentanyl (hydrochloride)
167	<i>meta</i> -methyl Acetyl fentanyl (hydrochloride)
168	2'-methyl Fentanyl (hydrochloride)
169	3'-methyl Fentanyl (hydrochloride)
170	Remifentanil Acid (trifluoroacetate salt)
171	(\pm)- <i>cis</i> -3-methyl Norfentanyl
172	N-Benzyl phenyl norfentanyl (hydrochloride)
173	<i>para</i> -fluoro 4-ANBP
174	Despropionyl <i>para</i> -Fluorofentanyl
175	Despropionyl <i>meta</i> -Methylfentanyl
176	N-benzyl <i>para</i> -fluoro Norfentanyl (hydrochloride)
177	Despropionyl 2'-fluoro <i>ortho</i> -Fluorofentanyl

No.	Compound Name
178	(\pm)- <i>cis</i> -Isofentanyl (hydrochloride)
179	N-methyl Norfentanyl (hydrochloride)
180	N-benzyl <i>para</i> -fluoro Cyclopropyl norfentanyl (hydrochloride)
181	<i>para</i> -Bromofentanyl
182	<i>para</i> -Toluoyl fentanyl (hydrochloride)
183	N-(MDA) Fentanyl
184	N-(Phentermine) Fentanyl
185	N-(6-APB) Fentanyl
186	N-(6-APDB) Fentanyl
187	N-(2-APB) Fentanyl
188	2',5'-dimethoxy Fentanyl (hydrochloride)
189	N-(2C-I) Fentanyl
190	N-(2C-B) Fentanyl
191	N-(2C-C) Fentanyl
192	N-(2C-N) Fentanyl
193	N-(2C-TFM) Fentanyl
194	N-(2C-D) Fentanyl
195	N-(2C-G) Fentanyl
196	N-(2C-E) Fentanyl
197	N-(2C-iP) Fentanyl
198	N-(2C-P) Fentanyl
199	N-(2C-T) Fentanyl
200	N-(2C-T-2) Fentanyl
201	N-(2C-T-4) Fentanyl
202	N-(2C-T-7) Fentanyl
203	N-(2C-B-fly) Fentanyl
204	N-(3C-B-fly) Fentanyl
205	N-(2,5-DMA) Fentanyl
206	N-(3,4,5-TMA) Fentanyl
207	N-(DOC) Fentanyl
208	N-(DOB) Fentanyl
209	N-(DOI) Fentanyl
210	N-(DOM) Fentanyl (hydrochloride)
211	N-(DOET) Fentanyl
212	N-(DOBU) Fentanyl