

Compounds in the FAS - Emergent Drug Panel & Internal Standard Kits



The Emergent Drug Panel (EDP) and EDP - Version 1 (EDP-V1) expand coverage to co-drugs that may be found in fentanyl-containing samples. The EDP and EDP-V1 provide coverage of 80 emergent synthetic opioids, benzodiazepines, stimulants, hallucinogens, synthetic cannabinoids, and xylazine.

The Emergent Drug Panel - Internal Standard (EDP-IS) contains isotopically labeled versions of select compounds included in the EDP and FAS Kits for use as internal standards in mass spectrometry-based analyses.

Emergent Drug Panel (EDP)

- Provides coverage of 60 fentanyl co-drugs
- Released April 2023
- Includes listed compound Nos. 1-60

No.	Compound Name
1	Clonazepam
2	Etizolam
3	Flualprazolam
4	Flubromazolam
5	Desalkylflurazepam
6	Bromazolam
7	Adinazolam
8	Deschloroetizolam
9	(±)-Meclonazepam
10	Flubromazepam
11	Pyrazolam
12	Diclazepam
13	Bromazepam
14	Metizolam
15	8-Aminoclonazepam
16	Phenazepam
17	Fluclotizolam
18	N-Pyrrolidino Etonitazene
19	Etodesnitazene (citrate)
20	Protonitazene (hydrochloride)
21	Dipyanone (hydrochloride)
22	N-Piperidinyl Etonitazene (citrate)

No.	Compound Name
23	Butonitazene
24	Eutylone (hydrochloride)
25	α-Pyrrolidinoisohexanophenone (hydrochloride)
26	Deschloroketamine (hydrochloride)
27	3-hydroxy PCP (hydrochloride)
28	2-fluoro Deschloroketamine (hydrochloride)
29	3-methoxy PCP (hydrochloride)
30	3-chloro PCP (hydrochloride)
31	α-Ethylaminohexanophenone (hydrochloride)
32	Pentylone (hydrochloride)
33	α-Pyrrolidinohexanophenone (hydrochloride)
34	N-ethyl Pentylone (hydrochloride)
35	3,4-Methylenedioxy-N-benzylcathinone (hydrochloride)
36	4-methoxy PCP (hydrochloride)
37	MDMB-4en-PINACA
38	N-butyl Pentylone (hydrochloride)
39	3,4-Methylenedioxy PV8 (hydrochloride)
40	α-Ethylaminopentiophenone (hydrochloride)
41	deschloro-N-ethyl-Ketamine (hydrochloride)
42	4'-chloro-α-Pyrrolidinovalerophenone (hydrochloride)
43	α-Pyrrolidinopentiophenone (hydrochloride)
44	4-hydroxy DIPT (hydrochloride)
45	3-hydroxy PCE (hydrochloride)
46	5-fluoro MDMB-PICA
47	Xylazine
48	4-fluoro MDMB-BUTINACA

Compounds in the FAS - Emergent Drug Panel & Internal Standard Kits

No.	Compound Name
49	ADB-BINACA
50	(S)-5-fluoro ADB
51	4-fluoro MDMB-BUTICA
52	ADB-4en-PINACA
53	ADB-HEXINACA
54	4-fluoro ABUTINACA
55	ADB-BUTINACA
56	ACHMINACA
57	4-cyano CUMYL-BUTINACA
58	MMB-FUBINACA
59	5-fluoro EDMB-PINACA
60	5-fluoro EDMB-PICA

Emergent Drug Panel - Version 1 (EDP-V1)

- Provides coverage of 20 fentanyl co-drugs
- Released April 2024
- Includes listed compound Nos. 61-80

No.	Compound Name
61	4'-chloro Deschloroalprazolam
62	Desalkylgizapem
63	N,N-Dimethylpentylone (hydrochloride)
64	3,4-Methylenedioxy- α -Pyrrolidinohexanophenone (hydrochloride)
65	3,4-Methylenedioxy- α -propylaminobutiophenone (hydrochloride)
66	3,4-Methylenedioxy- α -Cyclohexylaminopropiophenone (hydrochloride)
67	2-Methylmethcathinone (hydrochloride)
68	N-cyclohexyl Butylone (hydrochloride)
69	3-Methylmethcathinone (hydrochloride)
70	Butylone (hydrochloride)
71	3-methoxy PCE (hydrochloride)
72	ADB-FUBIATA
73	4-methoxy PCE (hydrochloride)
74	BZO-POXIZID
75	BZO-CHMOXIZID
76	5-fluoro BZO-POXIZID
77	MDA 19

No.	Compound Name
78	CH-PIATA
79	BZO-4en-POXIZID
80	ADB-5'Br-BUTINACA

Emergent Drug Panel - Internal Standard (EDP-IS)

- Contains isotopically labeled versions of select compounds in the EDP and FAS Kits
- Released April 2024
- Includes listed compound Nos. 81-104

No.	Compound Name
81	Clonazolam-d ₄
82	Etizolam-d ₄
83	Flualprazolam-d ₄
84	Flubromazolam-d ₄
85	Bromazolam-d ₅
86	Xylazine-d ₆
87	Flubromazepam-d ₄
88	<i>para</i> -Fluorofentanyl-d ₅ (hydrochloride)
89	Metonitazene-d ₃ (citrate)
90	Isotonitazene-d ₇
91	Brorphine-d ₇
92	Carfentanil- ¹³ C ₆
93	N-Pyrrolidino Etonitazene-d ₅
94	Etodesnitazene-d ₅ (citrate)
95	2-methyl AP-237-d ₇ (hydrochloride)
96	Protonitazene-d ₇ (hydrochloride)
97	Eutylone-d ₅ (hydrochloride)
98	α -Pyrrolidinoisohexanophenone-d ₈ (hydrochloride)
99	3-hydroxy PCP-d ₁₀ (hydrochloride)
100	α -Pyrrolidinohexanophenone-d ₈ (hydrochloride)
101	MDMB-4en-PINACA-d ₄
102	5-fluoro MDMB-PICA-d ₅
103	(S)-ADB-BINACA-d ₅
104	4-fluoro MDMB-BUTINACA-d ₄