

PRODUCT INFORMATION



Cathinone Analytical Standards Panel

Item No. 31616

Storage: -20°C
Stability: ≥3 years

Laboratory Procedures

Uncap each vial to be used. Add 500 µl of methanol to each vial. This will provide a 200 µg/ml standard solution for each analyte. Re-cap the vials and place plate on a plate mixer or vortexer. Mix for a minimum of 15 minutes to ensure full reconstitution.

The plate included in this panel contains a residual amount of glycerol to aid in the reconstitution of the analyte in methanol. Recovery in methanol has been validated for all analytes on the plate. Recovery from other solvents has not been verified.

Description

The Cathinone Analytical Standards Panel contains 239 analytical reference materials and standards categorized as cathinones and cathinone metabolites. These compounds are supplied at 100 µg/well in a 96-well Matrix™ tube rack format for rapid screening or building in-house spectral libraries. The plate included in this panel contains a residual amount of glycerol to aid in the reconstitution of the analyte in methanol. Recovery in methanol has been validated for all analytes on the plate. Recovery from other solvents has not been verified.

Please review the product insert for a full list of targets. The Cathinone Analytical Standards Panel contains compounds regulated as Schedule I compounds in the United States and is regulated as a Schedule I item. This product is intended for research and forensic applications.

Panel Contents

Plate	Well	Contents	Item Number	Molecular Formula	CAS Number
1	A1	<i>Unused</i>			
1	A2	(-)-(S)-Cathinone (hydrochloride)	13869	C ₉ H ₁₁ NO • HCl	72739-14-1
1	A3	4-methyl Pentedrone (hydrochloride)	9002180	C ₁₃ H ₁₉ NO • HCl	
1	A4	Methcathinone (hydrochloride)	11709	C ₁₀ H ₁₃ NO • HCl	49656-78-2
1	A5	2-chloro-N,N-Dimethylcathinone (hydrochloride)	24846	C ₁₁ H ₁₄ ClNO • HCl	856609-26-2
1	A6	2-Fluoroisocathinone (hydrochloride)	11607	C ₉ H ₁₀ FNO • HCl	1909305-71-0
1	A7	4-chloro Buphedrone (hydrochloride)	25298	C ₁₁ H ₁₄ ClNO • HCl	
1	A8	2-Thiothinone (hydrochloride)	9002101	C ₈ H ₁₁ NOS • HCl	54817-67-3
1	A9	Mephedrone (hydrochloride)	10801	C ₁₁ H ₁₅ NO • HCl	1189726-22-4
1	A10	3-methyl-α-Pyrrolidinopropiophenone (hydrochloride)	11485	C ₁₄ H ₁₉ NO • HCl	2749302-38-1
1	A11	3-Methylmethcathinone (hydrochloride)	11224	C ₁₁ H ₁₅ NO • HCl	1246816-62-5
1	A12	<i>Unused</i>			
1	B1	<i>Unused</i>			
1	B2	Ethcathinone (hydrochloride)	11241	C ₁₁ H ₁₅ NO • HCl	51553-17-4
1	B3	3,4-dichloro-α-Pyrrolidinoisohexanophenone (hydrochloride)	30249	C ₁₆ H ₂₁ Cl ₂ NO • HCl	
1	B4	N,N-Dimethylcathinone (hydrochloride)	9001144	C ₁₁ H ₁₅ NO • HCl	10105-90-5
1	B5	Mephedrone metabolite (hydrochloride) ((±)-Ephedrine stereochemistry)	14756	C ₁₁ H ₁₇ NO • HCl	117342-78-6
1	B6	4-ethyl Pentedrone (hydrochloride)	22993	C ₁₄ H ₂₁ NO • HCl	2748592-71-2
1	B7	Mephedrone metabolite (hydrochloride) ((±)-Pseudoephedrine stereochemistry)	9001434	C ₁₁ H ₁₇ NO • HCl	
1	B8	α-ethylamino-3,3-Dimethylbutyrophenone (hydrochloride)	25657	C ₁₄ H ₂₁ NO • HCl	2707567-71-1

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/23/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Plate	Well	Contents	Item Number	Molecular Formula	CAS Number
1	B9	3-Fluoromethcathinone (hydrochloride)	10730	C ₁₀ H ₁₂ FNO • HCl	1346600-40-5
1	B10	α-Ethylaminopentiophenone (hydrochloride)	14280	C ₁₃ H ₁₉ NO • HCl	18268-16-1
1	B11	2-Fluoromethcathinone (hydrochloride)	9001135	C ₁₀ H ₁₂ FNO • HCl	1346599-37-8
1	B12	<i>Unused</i>			
1	C1	<i>Unused</i>			
1	C2	α-Dimethylaminopentiophenone (hydrochloride)	9001579	C ₁₃ H ₁₉ NO • HCl	
1	C3	4-Fluoromethcathinone metabolite (hydrochloride) ((±)-Ephedrine stereochemistry)	14757	C ₁₀ H ₁₄ FNO • HCl	7799-72-6
1	C4	α-Methylaminohexanophenone (hydrochloride)	9002181	C ₁₃ H ₁₉ NO • HCl	
1	C5	MTTA (hydrochloride)	17121	C ₁₂ H ₁₅ NO • HCl	2749302-69-8
1	C6	4'-fluoro-α-Pyrrolidinopropiophenone (hydrochloride)	14448	C ₁₃ H ₁₆ FNO • HCl	2748590-18-1
1	C7	2-Methylethcathinone (hydrochloride)	11221	C ₁₂ H ₁₇ NO • HCl	2705355-07-1
1	C8	3,4-Dimethoxymethcathinone (hydrochloride)	17447	C ₁₂ H ₁₇ NO ₃ • HCl	22930-82-1
1	C9	2,3-Dimethylmethcathinone (hydrochloride)	11225	C ₁₂ H ₁₇ NO • HCl	1797981-99-7
1	C10	4-chloro Pentedrone (hydrochloride)	19383	C ₁₂ H ₁₆ ClNO • HCl	2749282-78-6
1	C11	4-Methylbuphedrone (hydrochloride)	11486	C ₁₂ H ₁₇ NO • HCl	1336911-98-8
1	C12	<i>Unused</i>			
1	D1	<i>Unused</i>			
1	D2	Isopentredone (hydrochloride)	11563	C ₁₂ H ₁₇ NO • HCl	1429402-13-0
1	D3	bk-EABDI (hydrochloride)	17990	C ₁₅ H ₂₁ NO • HCl	2726084-56-4
1	D4	N-Ethylbuphedrone (hydrochloride)	11665	C ₁₂ H ₁₇ NO • HCl	2731709-73-0
1	D5	4-fluoro Buphedrone (hydrochloride)	15165	C ₁₁ H ₁₄ FNO • HCl	2624137-27-3
1	D6	2,3-Methylenedioxymethcathinone (hydrochloride)	9001133	C ₁₁ H ₁₃ NO ₃ • HCl	1797884-10-6
1	D7	3-Chloromethcathinone (hydrochloride)	17394	C ₁₀ H ₁₂ ClNO • HCl	1607439-32-6
1	D8	N-methoxy Mephedrone (hydrochloride)	20808	C ₁₂ H ₁₇ NO ₂ • HCl	2703493-29-0
1	D9	α-Pyrrolidinopropiophenone (hydrochloride)	10445	C ₁₃ H ₁₇ NO • HCl	92040-10-3
1	D10	4-fluoro Pentedrone (hydrochloride)	9002182	C ₁₂ H ₁₆ FNO • HCl	2469350-88-5
1	D11	3-Ethylethcathinone (hydrochloride)	11198	C ₁₃ H ₁₉ NO • HCl	2446466-61-9
1	D12	<i>Unused</i>			
1	E1	<i>Unused</i>			
1	E2	2-Ethylethcathinone (hydrochloride)	11199	C ₁₃ H ₁₉ NO • HCl	2446466-59-5
1	E3	Methedrone (hydrochloride)	10529	C ₁₁ H ₁₅ NO ₂ • HCl	879665-92-6
1	E4	2,3-Dimethylethcathinone (hydrochloride)	11227	C ₁₃ H ₁₉ NO • HCl	2705531-06-0
1	E5	3-Methoxymethcathinone (hydrochloride)	9001187	C ₁₁ H ₁₅ NO ₂ • HCl	1435933-70-2
1	E6	2,4-Dimethylethcathinone (hydrochloride)	11242	C ₁₃ H ₁₉ NO • HCl	2702352-15-4
1	E7	Pentredone metabolite (hydrochloride) ((±)-Ephedrine stereochemistry)	9001428	C ₁₂ H ₁₉ NO • HCl	1422513-91-4
1	E8	4-methyl-α-Ethylaminobutiophenone (hydrochloride)	11489	C ₁₃ H ₁₉ NO • HCl	18268-19-4
1	E9	Pentredone metabolite (hydrochloride) ((±)-Pseudoephedrine stereochemistry)	9001435	C ₁₂ H ₁₉ NO • HCl	
1	E10	Butylone (hydrochloride)	10393	C ₁₂ H ₁₅ NO ₃ • HCl	17762-90-2
1	E11	3,4-Dimethylmethcathinone metabolite (hydrochloride) (±)-Pseudoephedrine stereochemistry)	9001437	C ₁₂ H ₁₉ NO • HCl	4865-62-7
1	E12	<i>Unused</i>			
1	F1	<i>Unused</i>			

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Plate	Well	Contents	Item Number	Molecular Formula	CAS Number
1	F2	2-Fluoroethcathinone (hydrochloride)	11229	C ₁₁ H ₁₄ FNO • HCl	2446466-63-1
1	F3	3,4-Dimethylmethcathinone (hydrochloride)	9001098	C ₁₂ H ₁₇ NO • HCl	1081772-06-6
1	F4	4-Fluoroethcathinone (hydrochloride)	11231	C ₁₁ H ₁₄ FNO • HCl	2705473-25-0
1	F5	3-Methylbuphedrone (hydrochloride)	12001	C ₁₂ H ₁₇ NO • HCl	
1	F6	4-Chloromethcathinone (hydrochloride)	16436	C ₁₀ H ₁₂ ClNO • HCl	2319878-22-1
1	F7	Propylcathinone (hydrochloride)	30941	C ₁₂ H ₁₇ NO • HCl	879667-46-6
1	F8	2-Chloromethcathinone (hydrochloride)	17744	C ₁₀ H ₁₂ ClNO • HCl	90869-66-2
1	F9	4-Ethylmethcathinone (hydrochloride)	9001078	C ₁₂ H ₁₇ NO • HCl	1391053-87-4
1	F10	4-Ethylethcathinone (hydrochloride)	11197	C ₁₃ H ₁₉ NO • HCl	2446466-62-0
1	F11	3-Ethylmethcathinone (hydrochloride)	9001082	C ₁₂ H ₁₇ NO • HCl	2446466-60-8
1	F12	Unused			
1	G1	Unused			
1	G2	4-Chlorocathinone (hydrochloride)	30379	C ₉ H ₁₀ ClNO • HCl	23184-97-6
1	G3	4-ethyl-N,N-Dimethylcathinone (hydrochloride)	11207	C ₁₃ H ₁₉ NO • HCl	2702382-98-5
1	G4	3,4-Methylenedioxy PV9 (hydrochloride)	16359	C ₁₉ H ₂₇ NO ₃ • HCl	24646-40-0
1	G5	3,4-Dimethylethcathinone (hydrochloride)	11228	C ₁₃ H ₁₉ NO • HCl	2748561-69-3
1	G6	4-Methylethcathinone metabolite (hydrochloride) ((±)-Ephedrine stereochemistry)	14758	C ₁₂ H ₁₉ NO • HCl	
1	G7	Diethylcathinone (hydrochloride)	11333	C ₁₃ H ₁₉ NO • HCl	134-80-5
1	G8	5-methoxy N,N-diethyl Hexylone (hydrochloride)	28234	C ₁₈ H ₂₇ NO ₄ • HCl	
1	G9	Mexedrone (hydrochloride)	9002622	C ₁₂ H ₁₇ NO ₂ • HCl	2749282-69-5
1	G10	4-Methylethcathinone metabolite (hydrochloride) ((±)-Pseudoephedrine stereochemistry)	9001436	C ₁₂ H ₁₉ NO • HCl	
1	G11	4-Chloroethcathinone (hydrochloride)	18912	C ₁₁ H ₁₄ ClNO • HCl	22198-75-0
1	G12	Unused			
1	H1	Unused			
1	H2	3-Fluoroethcathinone (hydrochloride)	11230	C ₁₁ H ₁₄ FNO • HCl	
1	H3	3-Chloroethcathinone (hydrochloride)	21261	C ₁₁ H ₁₄ ClNO • HCl	2748225-41-2
1	H4	2,3-Methylenedioxy-α-ethylaminopropiophenone (hydrochloride)	9002142	C ₁₂ H ₁₅ NO ₃ • HCl	
1	H5	Methylone (hydrochloride)	10986	C ₁₁ H ₁₃ NO ₃ • HCl	186028-80-8
1	H6	4-methyl-N,N-Dimethylcathinone (hydrochloride)	17393	C ₁₂ H ₁₇ NO • HCl	1448845-14-4
1	H7	4-methoxy-N,N-Dimethylcathinone (hydrochloride)	11666	C ₁₂ H ₁₇ NO ₂ • HCl	1089307-23-2
1	H8	4-Methylethcathinone (hydrochloride)	9001069	C ₁₂ H ₁₇ NO • HCl	1266688-86-1
1	H9	4-methyl-N-Methylbuphedrone (hydrochloride)	11667	C ₁₃ H ₁₉ NO • HCl	2748589-93-5
1	H10	2-Ethylmethcathinone (hydrochloride)	9001081	C ₁₂ H ₁₇ NO • HCl	
1	H11	α-Methylaminoisohexanophenone (hydrochloride)	30240	C ₁₃ H ₁₉ NO • HCl	
1	H12	Unused			
2	A1	Unused			
2	A2	2-Chloroethcathinone (hydrochloride)	21260	C ₁₁ H ₁₄ ClNO • HCl	879660-23-8
2	A3	DL-Cathinone (hydrochloride)	14138	C ₉ H ₁₁ NO • HCl	16735-19-6
2	A4	4-chloro-N,N-Dimethylcathinone (hydrochloride)	24432	C ₁₁ H ₁₄ ClNO • HCl	2493977-47-0
2	A5	nor-Mephedrone (hydrochloride)	9000940	C ₁₀ H ₁₃ NO • HCl	6941-17-9

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Plate	Well	Contents	Item Number	Molecular Formula	CAS Number
2	A6	3-chloro-N,N-Dimethylcathinone (hydrochloride)	24847	C ₁₁ H ₁₄ CINO • HCl	514168-22-0
2	A7	4-Fluoroisocathinone (hydrochloride)	9001146	C ₉ H ₁₀ FNO • HCl	2093282-02-9
2	A8	α-Piperidinobutiophenone (hydrochloride)	9001513	C ₁₅ H ₂₁ NO • HCl	92728-82-0
2	A9	2-methyl-α-Pyrrolidinopropiophenone (hydrochloride)	11484	C ₁₄ H ₁₉ NO • HCl	2749897-10-5
2	A10	2-Methylmethcathinone (hydrochloride)	11223	C ₁₁ H ₁₅ NO • HCl	1246815-51-9
2	A11	α-Pyrrolidinobutiophenone (hydrochloride)	9001195	C ₁₄ H ₁₉ NO • HCl	13415-54-8
2	A12	Unused			
2	B1	Unused			
2	B2	4-methyl-α-Ethylaminopentiophenone (hydrochloride)	15609	C ₁₄ H ₂₁ NO • HCl	18297-05-7
2	B3	Buphedrone (hydrochloride)	11283	C ₁₁ H ₁₅ NO • HCl	166593-10-8
2	B4	4'-methyl-α-Pyrrolidinopropiophenone (hydrochloride)	10446	C ₁₄ H ₁₉ NO • HCl	1313393-58-6
2	B5	α-Ethylaminohexanophenone (hydrochloride)	20661	C ₁₄ H ₂₁ NO • HCl	18410-62-3
2	B6	Buphedrone metabolite (hydrochloride) ((±)-Ephedrine stereochemistry)	9001432	C ₁₁ H ₁₇ NO • HCl	63199-70-2
2	B7	4-Methyldiethcathinone (hydrochloride)	24799	C ₁₄ H ₂₁ NO • HCl	2749619-00-7
2	B8	Buphedrone metabolite (hydrochloride) ((±)-Pseudoephedrine stereochemistry)	9001438	C ₁₁ H ₁₇ NO • HCl	1246817-96-8
2	B9	4'-Methyl-N-methylhexanophenone (hydrochloride)	9001515	C ₁₄ H ₂₁ NO • HCl	2749391-59-9
2	B10	4-Fluoromethcathinone (hydrochloride)	10859	C ₁₀ H ₁₂ FNO • HCl	7589-35-7
2	B11	3,4-EDMC (hydrochloride)	15167	C ₁₂ H ₁₅ NO ₃ • HCl	30253-44-2
2	B12	Unused			
2	C1	Unused			
2	C2	bk-MDEA (hydrochloride)	9001123	C ₁₂ H ₁₅ NO ₃ • HCl	1454266-19-3
2	C3	3',4'-dimethyl-α-Ethylaminovalerophenone (hydrochloride)	19210	C ₁₅ H ₂₃ NO • HCl	2748485-09-6
2	C4	4-Fluoromethcathinone metabolite (hydrochloride) ((±)-Pseudoephedrine stereochemistry)	9001433	C ₁₀ H ₁₄ FNO • HCl	7799-73-7
2	C5	3'-fluoro-α-Pyrrolidinopropiophenone (hydrochloride)	14447	C ₁₃ H ₁₆ FNO • HCl	1214940-15-4
2	C6	Pentadrone (hydrochloride)	11011	C ₁₂ H ₁₇ NO • HCl	879669-95-1
2	C7	α-Pyrrolidinobuthiophenone (hydrochloride)	15358	C ₁₂ H ₁₇ NOS • HCl	2518265-88-6
2	C8	3-Methylethcathinone (hydrochloride)	11222	C ₁₂ H ₁₇ NO • HCl	2493976-59-1
2	C9	4-fluoro-α-Ethylaminovalerophenone (hydrochloride)	19211	C ₁₃ H ₁₈ FNO • HCl	
2	C10	2,4-Dimethylmethcathinone (hydrochloride)	11243	C ₁₂ H ₁₇ NO • HCl	2412098-39-4
2	C11	4-chloro-N-Isopropylcathinone (hydrochloride)	19392	C ₁₂ H ₁₆ CINO • HCl	2749504-71-8
2	C12	Unused			
2	D1	Unused			
2	D2	3,4-Dichloromethcathinone (hydrochloride)	25376	C ₁₀ H ₁₁ Cl ₂ NO • HCl	15861-85-5
2	D3	N-ethyl-N-Methylcathinone (hydrochloride)	11604	C ₁₂ H ₁₇ NO • HCl	1157739-24-6
2	D4	α-Pyrrolidinopentiophenone (hydrochloride)	9001083	C ₁₅ H ₂₁ NO • HCl	5485-65-4
2	D5	bk-DMBDB (hydrochloride)	9001125	C ₁₃ H ₁₇ NO ₃ • HCl	17763-12-1
2	D6	3-methyl-α-Pyrrolidinobutiophenone (hydrochloride)	9001189	C ₁₅ H ₂₁ NO • HCl	
2	D7	4'-methoxy-α-Ethylaminovalerophenone (hydrochloride)	19275	C ₁₄ H ₂₁ NO ₂ • HCl	17837-89-7
2	D8	α-Isopropylaminopentiophenone (hydrochloride)	20021	C ₁₄ H ₂₁ NO • HCl	18268-14-9
2	D9	6-methoxy Methylone (hydrochloride)	18029	C ₁₂ H ₁₅ NO ₄ • HCl	2749504-36-5
2	D10	4-methoxy-α-Pyrrolidinopropiophenone (hydrochloride)	10394	C ₁₄ H ₁₉ NO ₂ • HCl	1794760-01-2

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Plate	Well	Contents	Item Number	Molecular Formula	CAS Number
2	D11	α -Pyrrolidinopentiothiophenone (hydrochloride)	14182	$C_{13}H_{19}NOS \cdot HCl$	2748622-52-6
2	D12	<i>Unused</i>			
2	E1	<i>Unused</i>			
2	E2	α -Pyrrolidinopentiothiophenone metabolite 1 (trifluoroacetate salt)	14093	$C_{15}H_{23}NO \cdot CF_3COOH$	1797986-63-0
2	E3	bk-MDDMA (hydrochloride)	9001124	$C_{12}H_{15}NO_3 \cdot HCl$	109367-07-9
2	E4	N-ethyl Heptedrone (hydrochloride)	27327	$C_{15}H_{23}NO \cdot HCl$	
2	E5	3-Bromomethcathinone (hydrochloride)	14035	$C_{10}H_{12}BrNO \cdot HCl$	676487-42-6
2	E6	NRG-3 (hydrochloride)	14115	$C_{16}H_{19}NO \cdot HCl$	2701928-05-2
2	E7	3,4-Methylenedioxy-5-methylethcathinone (hydrochloride)	14078	$C_{13}H_{17}NO_3 \cdot HCl$	2748289-02-1
2	E8	3,4-dichloro-N,N-Dimethcathinone (hydrochloride)	25377	$C_{11}H_{13}Cl_2NO \cdot HCl$	
2	E9	3,4-Methylenedioxy-N-propylcathinone (hydrochloride)	19371	$C_{13}H_{17}NO_3 \cdot HCl$	2705469-52-7
2	E10	3',4'-trimethylene- α -Ethylaminovalerophenone (hydrochloride)	17991	$C_{16}H_{23}NO \cdot HCl$	2748590-67-0
2	E11	3,4-methylenedioxy- α -methylamino-Isovalerophenone (hydrochloride)	23711	$C_{13}H_{17}NO_3 \cdot HCl$	2702751-83-3
2	E12	<i>Unused</i>			
2	F1	<i>Unused</i>			
2	F2	2,3-Methylenedioxy- α -ethylaminobutiophenone (hydrochloride)	30595	$C_{13}H_{17}NO_3 \cdot HCl$	
2	F3	3,4-Methylenedioxy- α -Pyrrolidinopropiophenone (hydrochloride)	10439	$C_{14}H_{17}NO_3 \cdot HCl$	24698-57-5
2	F4	Eutylone (hydrochloride)	9001103	$C_{13}H_{17}NO_3 \cdot HCl$	17764-18-0
2	F5	2-methyl- α -Pyrrolidinobutiophenone (hydrochloride)	9001188	$C_{15}H_{21}NO \cdot HCl$	
2	F6	α -Butylaminohexanophenone (hydrochloride)	27728	$C_{16}H_{25}NO \cdot HCl$	18296-66-7
2	F7	4-methyl- α -Pyrrolidinobutiophenone (hydrochloride)	9001190	$C_{15}H_{21}NO \cdot HCl$	1214-15-9
2	F8	4'-chloro- α -Pyrrolidinopropiophenone (hydrochloride)	9002179	$C_{13}H_{16}ClNO \cdot HCl$	93307-24-5
2	F9	3',4'-trimethylene- α -methylamino-Valerophenone (hydrochloride)	9001514	$C_{15}H_{21}NO \cdot HCl$	
2	F10	3,4-Methylenedioxy-N,N-diethylcathinone (hydrochloride)	18586	$C_{14}H_{19}NO_3 \cdot HCl$	2749303-12-4
2	F11	4-methoxy- α -Pyrrolidinopropiophenone (tosylate)	10449	$C_{14}H_{19}NO_2 \cdot C_7H_8O_3S$	2748561-60-4
2	F12	<i>Unused</i>			
2	G1	<i>Unused</i>			
2	G2	4-fluoro IPV (hydrochloride)	9002414	$C_{14}H_{20}FNO \cdot HCl$	
2	G3	4-chloro-N-Butylcathinone (hydrochloride)	19972	$C_{13}H_{18}ClNO \cdot HCl$	2748591-11-7
2	G4	4-Bromomethcathinone (hydrochloride)	12089	$C_{10}H_{12}BrNO \cdot HCl$	135333-27-6
2	G5	4-methyl-N-Ethylhexanophenone (hydrochloride)	28365	$C_{15}H_{23}NO \cdot HCl$	
2	G6	2,3-Pentylone isomer (hydrochloride)	11463	$C_{13}H_{17}NO_3 \cdot HCl$	2749428-57-5
2	G7	3,4-Dichloroethcathinone (hydrochloride)	25375	$C_{11}H_{13}Cl_2NO \cdot HCl$	
2	G8	3,4-Methylenedioxy-N-isopropylcathinone (hydrochloride)	19370	$C_{13}H_{17}NO_3 \cdot HCl$	2749298-40-4
2	G9	Pyrovalerone (hydrochloride)	10836	$C_{16}H_{23}NO \cdot HCl$	1147-62-2
2	G10	N-Methylethylone (hydrochloride)	23710	$C_{13}H_{17}NO_3 \cdot HCl$	2703646-17-5
2	G11	α -Pyrrolidinoisohexanophenone (hydrochloride)	21682	$C_{16}H_{23}NO \cdot HCl$	2705245-60-7
2	G12	<i>Unused</i>			
2	H1	<i>Unused</i>			
2	H2	α -Pyrrolidinoisohexanophenone (hydrochloride)	9001934	$C_{16}H_{23}NO \cdot HCl$	13415-59-3
2	H3	Pentylone (hydrochloride)	9000746	$C_{13}H_{17}NO_3 \cdot HCl$	17763-01-8
2	H4	4-methoxy- α -Pyrrolidinobutiophenone (hydrochloride)	15354	$C_{15}H_{21}NO_2 \cdot HCl$	2749897-35-4
2	H5	α -Isobutylaminoisohexanophenone (hydrochloride)	26944	$C_{16}H_{25}NO \cdot HCl$	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Plate	Well	Contents	Item Number	Molecular Formula	CAS Number
2	H6	4-fluoro- α -Pyrrolidinobutiophenone (hydrochloride)	15516	C ₁₄ H ₁₈ FNO • HCl	2749302-50-7
2	H7	α -Diethylaminohexanophenone (hydrochloride)	27729	C ₁₆ H ₂₅ NO • HCl	
2	H8	N-acetyl-3,4-Methylenedioxy-methcathinone	9001492	C ₁₃ H ₁₅ NO ₄	1227293-15-3
2	H9	N-ethyl Pentylone (hydrochloride)	16037	C ₁₄ H ₁₉ NO ₃ • HCl	17763-02-9
2	H10	5-methoxy Methylone (hydrochloride)	9002195	C ₁₂ H ₁₅ NO ₄ • HCl	2702151-27-5
2	H11	3',4'-Methylenedioxy-N-tert-butylcathinone (hydrochloride)	18587	C ₁₄ H ₁₉ NO ₃ • HCl	2469270-98-0
2	H12	Unused			
3	A1	Unused			
3	A2	3,4-Methylenedioxy- α -methylaminohexanophenone (hydrochloride)	29753	C ₁₄ H ₁₉ NO ₃ • HCl	27912-40-9
3	A3	4-fluoro- α -Pyrrolidinoisohexanophenone (hydrochloride)	29388	C ₁₆ H ₂₂ FNO • HCl	
3	A4	N,N-Dimethylpentylone (hydrochloride)	9001933	C ₁₄ H ₁₉ NO ₃ • HCl	17763-13-2
3	A5	4-fluoro-3-methyl- α -Pyrrolidinovalerophenone (hydrochloride)	31050	C ₁₆ H ₂₂ FNO • HCl	
3	A6	3,4-methylenedioxy- α -ethylamino-Isovalerophenone (hydrochloride)	9002600	C ₁₄ H ₁₉ NO ₃ • HCl	
3	A7	DL-4662 (hydrochloride)	9002077	C ₁₅ H ₂₃ NO ₃ • HCl	
3	A8	3-fluoro- α -Pyrrolidinopentiophenone (hydrochloride)	24496	C ₁₅ H ₂₀ FNO • HCl	2748623-76-7
3	A9	Benzedrone (hydrochloride)	11915	C ₁₇ H ₁₉ NO • HCl	1797979-43-1
3	A10	4-Bromoethcathinone (hydrochloride)	17395	C ₁₁ H ₁₄ BrNO • HCl	135333-26-5
3	A11	N-sec-butyl Pentylone (hydrochloride)	27084	C ₁₆ H ₂₃ NO ₃ • HCl	
3	A12	Unused			
3	B1	Unused			
3	B2	4'-methyl- α -Pyrrolidinoisohexanophenone (hydrochloride)	10448	C ₁₇ H ₂₅ NO • HCl	1391052-36-0
3	B3	α -Pyrrolidinocyclohexanophenone (hydrochloride)	30241	C ₁₈ H ₂₅ NO • HCl	1803168-16-2
3	B4	3',4'-dimethyl- α -Pyrrolidinovalerophenone (hydrochloride)	19212	C ₁₇ H ₂₅ NO • HCl	2748485-15-4
3	B5	3-desoxy-3,4-Methylenedioxy Pyrovalerone (hydrochloride)	16437	C ₁₇ H ₂₃ NO ₂ • HCl	2117405-33-9
3	B6	4-ethyl- α -Pyrrolidinopentiophenone (hydrochloride)	30226	C ₁₇ H ₂₅ NO • HCl	
3	B7	4-methyl PV8 (hydrochloride)	16356	C ₁₈ H ₂₇ NO • HCl	
3	B8	4-methoxy- α -Pyrrolidinopentiophenone (hydrochloride)	14097	C ₁₆ H ₂₃ NO ₂ • HCl	5537-19-9
3	B9	3-ethoxy-4-methoxy N,N-Diethylhexedrone (hydrochloride)	28235	C ₁₉ H ₃₁ NO ₃ • HCl	
3	B10	N-ethyl Hexylone (hydrochloride)	25315	C ₁₅ H ₂₁ NO ₃ • HCl	
3	B11	2,3-Methylenedioxy Pyrovalerone (hydrochloride)	9001051	C ₁₆ H ₂₁ NO ₃ • HCl	
3	B12	Unused			
3	C1	Unused			
3	C2	N-methyl-N-isopropyl Pentylone (hydrochloride)	27083	C ₁₆ H ₂₃ NO ₃ • HCl	
3	C3	3,4-Methylenedioxy- α -methylaminoisohexanophenone (hydrochloride)	30705	C ₁₄ H ₁₉ NO ₃ • HCl	
3	C4	N,N-dimethyl Heptylone (hydrochloride)	27115	C ₁₆ H ₂₃ NO ₃ • HCl	
3	C5	3',4'-methylenedioxy- α -dimethylamino-Isovalerophenone (hydrochloride)	9002599	C ₁₄ H ₁₉ NO ₃ • HCl	
3	C6	4'-chloro- α -Pyrrolidinovalerophenone (hydrochloride)	18326	C ₁₅ H ₂₀ ClNO • HCl	5537-17-7
3	C7	4-fluoro- α -Pyrrolidinopentiophenone (hydrochloride)	15166	C ₁₅ H ₂₀ FNO • HCl	850352-31-7
3	C8	4-chloro-N,N-Diethylpentedrone (hydrochloride)	27397	C ₁₅ H ₂₂ ClNO • HCl	
3	C9	N-methyl Benzedrone (hydrochloride)	21214	C ₁₈ H ₂₁ NO • HCl	2749433-00-7
3	C10	4-bromo- α -Methylaminovalerophenone (hydrochloride)	23309	C ₁₂ H ₁₆ BrNO • HCl	
3	C11	3',4'-tetramethylene- α -Pyrrolidinobutiophenone (hydrochloride)	17956	C ₁₈ H ₂₅ NO • HCl	2749504-65-0
3	C12	Unused			

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Plate	Well	Contents	Item Number	Molecular Formula	CAS Number
3	D1	<i>Unused</i>			
3	D2	3',4'-trimethylene- α -Pyrrolidinovalerophenone (hydrochloride)	18554	C ₁₈ H ₂₅ NO • HCl	
3	D3	PV8 (hydrochloride)	14762	C ₁₇ H ₂₅ NO • HCl	13415-55-9
3	D4	bk-2C-B (hydrochloride)	15328	C ₁₀ H ₁₂ BrNO ₃ • HCl	
3	D5	2,5-dimethyl- α -Pyrrolidinovalerophenone (hydrochloride)	23457	C ₁₇ H ₂₅ NO • HCl	
3	D6	PV9 (hydrochloride)	15062	C ₁₈ H ₂₇ NO • HCl	2749897-19-4
3	D7	3,4-Methylenedioxy- α -Pyrrolidinobutiophenone (hydrochloride)	10437	C ₁₅ H ₁₉ NO ₃ • HCl	24622-60-4
3	D8	3,4-Methylenedioxy Pyrovalerone (hydrochloride)	10684	C ₁₆ H ₂₁ NO ₃ • HCl	24622-62-6
3	D9	3,4-Methylenedioxy Pyrovalerone metabolite 2 (hydrochloride)	12004	C ₁₅ H ₂₁ NO ₃ • HCl	2748289-38-3
3	D10	3-methoxy-4-ethoxy N,N-Diethylhexedrone (hydrochloride)	28236	C ₁₉ H ₃₁ NO ₃ • HCl	
3	D11	4-bromo- α -Pyrrolidinovalerophenone (hydrochloride)	19329	C ₁₅ H ₂₀ BrNO • HCl	850352-17-9
3	D12	<i>Unused</i>			
3	E1	<i>Unused</i>			
3	E2	3,4-Methylenedioxy Pyrovalerone metabolite 1 (hydrochloride)	10518	C ₁₆ H ₂₃ NO ₃ • HCl	2748590-34-1
3	E3	4-Methylthio-N-benzylcathinone (hydrochloride)	27185	C ₁₇ H ₁₉ NOS • HCl	2706488-88-0
3	E4	N-butyl Pentylone (hydrochloride)	26701	C ₁₆ H ₂₃ NO ₃ • HCl	17763-10-9
3	E5	3',4'-trimethylene- α -Pyrrolidinohexanophenone (hydrochloride)	18030	C ₁₉ H ₂₇ NO • HCl	2748304-64-3
3	E6	N-isopropyl Hexylone (hydrochloride)	27076	C ₁₆ H ₂₃ NO ₃ • HCl	27912-43-2
3	E7	3,4-Methylenedioxy- α -Pyrrolidinohexanophenone (hydrochloride)	16361	C ₁₇ H ₂₃ NO ₃ • HCl	24622-61-5
3	E8	N-propyl Hexylone (hydrochloride)	27078	C ₁₆ H ₂₃ NO ₃ • HCl	27912-42-1
3	E9	3,4-dimethoxy- α -Pyrrolidinopentiophenone (hydrochloride)	14106	C ₁₇ H ₂₅ NO ₃ • HCl	850351-99-4
3	E10	N-tert-butyl Pentylone (hydrochloride)	27080	C ₁₆ H ₂₃ NO ₃ • HCl	
3	E11	4-fluoro PV9 (hydrochloride)	15063	C ₁₈ H ₂₆ FNO • HCl	2117405-40-8
3	E12	<i>Unused</i>			
3	F1	<i>Unused</i>			
3	F2	4'-fluoro- α -Pyrrolidinohexanophenone (hydrochloride)	18818	C ₁₆ H ₂₂ FNO • HCl	2748289-49-6
3	F3	3',4'-trimethylene- α -Pyrrolidinobutiophenone (hydrochloride)	18031	C ₁₇ H ₂₃ NO • HCl	2749618-47-9
3	F4	3-fluoro- α -Pyrrolidinoisohexanophenone (hydrochloride)	29389	C ₁₆ H ₂₂ FNO • HCl	
3	F5	4-fluoro PV8 (hydrochloride)	15356	C ₁₇ H ₂₄ FNO • HCl	2748485-01-8
3	F6	α -Phthalimidopropiophenone	17141	C ₁₇ H ₁₃ NO ₃	19437-20-8
3	F7	2-methyl-4'-(methylthio)-2-Morpholinopropiophenone	21651	C ₁₅ H ₂₁ NO ₂ S	71868-10-5
3	F8	6-methoxy-N,N-Diethylpentylone (hydrochloride)	27399	C ₁₇ H ₂₅ NO ₄ • HCl	
3	F9	3-methoxy-4-ethoxy N,N-Diethylpentedrone (hydrochloride)	27626	C ₁₈ H ₂₉ NO ₃ • HCl	
3	F10	Naphyrone 1-naphthyl isomer (hydrochloride)	11240	C ₁₉ H ₂₃ NO • HCl	1349245-31-3
3	F11	4-Dimethylamino-N-benzylcathinone (hydrochloride)	28893	C ₁₈ H ₂₂ N ₂ O • 2HCl	2740524-43-8
3	F12	<i>Unused</i>			
3	G1	<i>Unused</i>			
3	G2	3,4-Methylenedioxy-N-benzylcathinone (hydrochloride)	9001330	C ₁₇ H ₁₇ NO ₃ • HCl	1823274-68-5
3	G3	N,N-Diethylpentylone (hydrochloride)	20801	C ₁₆ H ₂₃ NO ₃ • HCl	17763-15-4
3	G4	3',4'-tetramethylene- α -Pyrrolidinovalerophenone (hydrochloride)	17955	C ₁₉ H ₂₇ NO • HCl	2748590-73-8
3	G5	N-ethyl Heptylone (hydrochloride)	26923	C ₁₆ H ₂₃ NO ₃ • HCl	
3	G6	α -Pyrrolidinononanophenone (hydrochloride)	17138	C ₁₉ H ₂₉ NO • HCl	2748414-29-9
3	G7	3,4-Methylenedioxy- α -ethylmethylaminohexanophenone (hydrochloride)	27077	C ₁₆ H ₂₃ NO ₃ • HCl	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Plate	Well	Contents	Item Number	Molecular Formula	CAS Number
3	G8	4-methoxy PV8 (hydrochloride)	15497	C ₁₈ H ₂₇ NO ₂ • HCl	2749302-60-9
3	G9	N-methyl-N-propyl Pentylone (hydrochloride)	27079	C ₁₆ H ₂₃ NO ₃ • HCl	
3	G10	N,N-diethyl Hexylone (hydrochloride)	28475	C ₁₇ H ₂₅ NO ₃ • HCl	17763-17-6
3	G11	N-isobutyl Pentylone (hydrochloride)	27082	C ₁₆ H ₂₃ NO ₃ • HCl	
3	G12	Unused			
3	H1	Unused			
3	H2	4-fluoro PV8 piperidine analog (hydrochloride)	17519	C ₁₈ H ₂₆ FNO • HCl	2748623-02-9
3	H3	BMDB (hydrochloride)	24062	C ₁₈ H ₁₉ NO ₃ • HCl	1823865-05-9
3	H4	3,4-Methylenedioxy PV8 (hydrochloride)	16358	C ₁₈ H ₂₅ NO ₃ • HCl	24646-39-7
3	H5	4-methoxy PV9 (hydrochloride)	15064	C ₁₉ H ₂₉ NO ₂ • HCl	
3	H6	3,4-dimethoxy- α -Pyrrolidinohexanophenone (hydrochloride)	17137	C ₁₈ H ₂₇ NO ₃ • HCl	2748409-11-0
3	H7	5-methoxy-N,N-Diethylpentylone (hydrochloride)	27398	C ₁₇ H ₂₅ NO ₄ • HCl	
3	H8	4-chloro- α -Pyrrolidinohexanophenone (hydrochloride)	22490	C ₁₆ H ₂₂ ClNO • HCl	2748592-29-0
3	H9	Naphyrone (hydrochloride)	10517	C ₁₉ H ₂₃ NO • HCl	850352-11-3
3	H10	3-ethoxy-4-methoxy N,N-Diethylpentedrone (hydrochloride)	27627	C ₁₈ H ₂₉ NO ₃ • HCl	
3	H11	Unused			
3	H12	Unused			

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM