

PRODUCT INFORMATION



Synthetic Opioid Analytical Standards Panel 1

Item No. 44436

Supplied as: 100 µg/well
Storage: -20°C
Stability: ≥4 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 plates included in this panel contain 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; however, recovery from other solvents has not been verified.

Description

Synthetic Opioid Standards Panel 1 contains reference standards for 250 synthetic opioid compounds, including more than 210 fentanyl analogs. 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. Please review the product insert for a full list of targets. Synthetic Opioid Analytical Standards Panel 1 contains compounds regulated as Schedule I compounds in the United States; therefore, this panel is regulated as a Schedule I product in the United States. This product is intended for research and forensic applications.

Panel Contents

Plate	Well	Contents	Item Number	Molecular Formula	CAS Number
1	A1	Unused			
1	A2	α-methyl Acetyl fentanyl (hydrochloride)	20816	C ₂₂ H ₂₈ N ₂ O • HCl	2306822-12-6
1	A3	Acetyl norfentanyl (hydrochloride)	ISO00130	C ₁₃ H ₁₈ N ₂ O • HCl	22352-82-5
1	A4	ortho-methoxy Furanyl fentanyl	22805	C ₂₅ H ₂₈ N ₂ O ₃	101343-50-4
1	A5	ortho-methyl Furanyl fentanyl	22806	C ₂₅ H ₂₈ N ₂ O ₂	2309383-07-9
1	A6	4'-methyl Acetyl fentanyl (hydrochloride)	9002271	C ₂₂ H ₂₈ N ₂ O • HCl	1071703-95-1
1	A7	Thiophene fentanyl (hydrochloride)	22802	C ₂₄ H ₂₆ N ₂ OS • HCl	2306823-39-0
1	A8	para-chloro Cyclobutyl fentanyl (hydrochloride)	23339	C ₂₄ H ₂₉ ClN ₂ O • HCl	2748290-00-6
1	A9	N-methyl Cyclopropyl norfentanyl (hydrochloride)	24278	C ₁₆ H ₂₂ N ₂ O • HCl	2703474-67-1
1	A10	β-methyl Acetyl fentanyl (hydrochloride)	21736	C ₂₂ H ₂₈ N ₂ O • HCl	2749910-75-4
1	A11	4-ANPP	18810	C ₁₉ H ₂₄ N ₂	21409-26-7
1	A12	Unused			
1	B1	Unused			
1	B2	para-methyl Acrylfentanyl (hydrochloride)	23038	C ₂₃ H ₂₈ N ₂ O • HCl	2748409-55-2
1	B3	ortho-methyl Cyclopropyl fentanyl (hydrochloride)	23043	C ₂₄ H ₃₀ N ₂ O • HCl	2749910-65-2
1	B4	Fentanyl	23580	C ₂₂ H ₂₈ N ₂ O	437-38-7
1	B5	Norfentanyl	15899	C ₁₄ H ₂₀ N ₂ O	1609-66-1
1	B6	Cyclopropyl fentanyl (hydrochloride)	21739	C ₂₃ H ₂₈ N ₂ O • HCl	2306825-44-3
1	B7	Butyryl norfentanyl (hydrochloride)	19428	C ₁₅ H ₂₂ N ₂ O • HCl	2701720-35-4
1	B8	ortho-methyl Acetyl fentanyl (hydrochloride)	23548	C ₂₂ H ₂₈ N ₂ O • HCl	2748409-58-5
1	B9	α'-methyl Butyryl fentanyl (hydrochloride)	22994	C ₂₄ H ₃₂ N ₂ O • HCl	2748591-35-5

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, 03/10/2026

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	B10	Methacrylfentanyl	22884	C ₂₃ H ₂₈ N ₂ O	2361296-62-8
1	B11	α-methyl Butyryl fentanyl (hydrochloride)	20818	C ₂₄ H ₃₂ N ₂ O • HCl	2306825-91-0
1	B12	<i>Unused</i>			
1	C1	<i>Unused</i>			
1	C2	<i>para</i> -fluoro Crotonyl fentanyl	23544	C ₂₃ H ₂₇ FN ₂ O	
1	C3	α-methyl Fentanyl (hydrochloride)	18280	C ₂₃ H ₃₀ N ₂ O • HCl	1443-44-3
1	C4	Crotonyl fentanyl	22801	C ₂₃ H ₂₈ N ₂ O	760930-59-4
1	C5	<i>para</i> -methyl Cyclopropyl fentanyl (hydrochloride)	23045	C ₂₄ H ₃₀ N ₂ O • HCl	2749910-43-6
1	C6	<i>para</i> -Methoxyfentanyl (hydrochloride)	20035	C ₂₃ H ₃₀ N ₂ O ₂ • HCl	23609-41-8
1	C7	Cyclobutyl fentanyl (hydrochloride)	22389	C ₂₄ H ₃₀ N ₂ O • HCl	2306827-56-3
1	C8	<i>ortho</i> -methyl Acrylfentanyl (hydrochloride)	23036	C ₂₃ H ₂₈ N ₂ O • HCl	2749299-21-4
1	C9	(±)- <i>cis</i> -3-methyl Fentanyl (hydrochloride)	9002747	C ₂₃ H ₃₀ N ₂ O • HCl	78995-18-3
1	C10	Ocfentanil (hydrochloride)	18583	C ₂₂ H ₂₇ FN ₂ O ₂ • HCl	112964-97-3
1	C11	<i>meta</i> -Methylfentanyl (hydrochloride)	22633	C ₂₃ H ₃₀ N ₂ O • HCl	1465-22-1
1	C12	<i>Unused</i>			
1	D1	<i>Unused</i>			
1	D2	<i>ortho</i> -fluoro Acrylfentanyl (hydrochloride)	21932	C ₂₂ H ₂₅ FN ₂ O • HCl	2309383-10-4
1	D3	<i>para</i> -Chlorofentanyl (hydrochloride)	20034	C ₂₂ H ₂₇ ClN ₂ O • HCl	117994-27-1
1	D4	<i>para</i> -fluoro Cyclopropyl fentanyl (hydrochloride)	22749	C ₂₃ H ₂₇ FN ₂ O • HCl	2306825-53-4
1	D5	Butyryl fentanyl (hydrochloride)	14728	C ₂₃ H ₃₀ N ₂ O • HCl	1443-52-3
1	D6	Fentanyl Carbamate	23749	C ₂₂ H ₂₈ N ₂ O ₂	1465-20-9
1	D7	Isobutyryl fentanyl (hydrochloride)	18584	C ₂₃ H ₃₀ N ₂ O • HCl	117332-90-8
1	D8	<i>ortho</i> -methyl Methoxyacetyl fentanyl (hydrochloride)	22977	C ₂₃ H ₃₀ N ₂ O ₂ • HCl	2748485-23-4
1	D9	Benzyl Carfentanil (hydrochloride)	20350	C ₂₃ H ₂₈ N ₂ O ₃ • HCl	2749504-78-5
1	D10	α-methyl Thiofentanyl (hydrochloride)	20821	C ₂₁ H ₂₈ N ₂ OS • HCl	117332-94-2
1	D11	<i>para</i> -methoxy-Butyryl fentanyl (hydrochloride)	18089	C ₂₄ H ₃₂ N ₂ O ₂ • HCl	2306825-45-4
1	D12	<i>Unused</i>			
1	E1	<i>Unused</i>			
1	E2	<i>para</i> -chloro Cyclopropyl fentanyl (hydrochloride)	23342	C ₂₃ H ₂₇ ClN ₂ O • HCl	2748591-61-7
1	E3	<i>ortho</i> -Fluorobutyryl fentanyl (hydrochloride)	19567	C ₂₃ H ₂₉ FN ₂ O • HCl	2306823-44-7
1	E4	<i>para</i> -fluoro Acrylfentanyl	21116	C ₂₂ H ₂₅ FN ₂ O	2306823-27-6
1	E5	Cyclopentyl fentanyl (hydrochloride)	20439	C ₂₅ H ₃₂ N ₂ O • HCl	2306824-94-0
1	E6	Phenyl fentanyl (hydrochloride)	22551	C ₂₆ H ₂₈ N ₂ O • HCl	2309383-16-0
1	E7	Tetrahydrofuran fentanyl (hydrochloride)	20859	C ₂₄ H ₃₀ N ₂ O ₂ • HCl	2306822-94-4
1	E8	Methoxyacetyl fentanyl (hydrochloride)	20782	C ₂₂ H ₂₈ N ₂ O ₂ • HCl	101365-54-2
1	E9	<i>ortho</i> -Fluoroisobutyryl fentanyl (hydrochloride)	20207	C ₂₃ H ₂₉ FN ₂ O • HCl	2748591-21-9
1	E10	<i>para</i> -fluoro Furanyl fentanyl 3-furancarboxamide (hydrochloride)	22751	C ₂₄ H ₂₅ FN ₂ O ₂ • HCl	2749433-11-0
1	E11	<i>para</i> -Fluorobutyryl fentanyl (hydrochloride)	17049	C ₂₃ H ₂₉ FN ₂ O • HCl	2306823-23-2
1	E12	<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	F1	<i>Unused</i>			
1	F2	<i>para</i> -methyl Tetrahydrofuran fentanyl (hydrochloride)	23054	C ₂₅ H ₃₂ N ₂ O ₂ • HCl	2749433-21-2
1	F3	Phenylacetyl fentanyl (hydrochloride)	22743	C ₂₇ H ₃₀ N ₂ O • HCl	2749618-58-2
1	F4	<i>para</i> -Chloroisobutyryl fentanyl (hydrochloride)	20018	C ₂₃ H ₂₉ ClN ₂ O • HCl	2306827-85-8
1	F5	<i>meta</i> -Fluorobutyryl fentanyl (hydrochloride)	19566	C ₂₃ H ₂₉ FN ₂ O • HCl	2306827-52-9
1	F6	<i>ortho</i> -fluoro Furanyl fentanyl (hydrochloride)	23628	C ₂₄ H ₂₅ FN ₂ O ₂ • HCl	2749619-19-8
1	F7	<i>meta</i> -Fluoroisobutyryl fentanyl (hydrochloride)	20206	C ₂₃ H ₂₉ FN ₂ O • HCl	2306827-89-2
1	F8	<i>para</i> -chloro Methoxyacetyl fentanyl (hydrochloride)	23343	C ₂₂ H ₂₇ ClN ₂ O ₂ • HCl	101345-75-9
1	F9	Alfentanil (hydrochloride)	19292	C ₂₁ H ₃₂ N ₆ O ₃ • HCl	69049-06-5
1	F10	β ¹ -Phenyl fentanyl	23858	C ₂₈ H ₃₂ N ₂ O	79279-02-0
1	F11	Benzodioxole fentanyl	20858	C ₂₇ H ₂₈ N ₂ O ₃	2306823-01-6
1	F12	<i>Unused</i>			
1	G1	<i>Unused</i>			
1	G2	<i>para</i> -methyl Acetyl fentanyl (hydrochloride)	23191	C ₂₂ H ₂₈ N ₂ O • HCl	2748496-88-8
1	G3	<i>para</i> -methyl Isobutyryl fentanyl (hydrochloride)	23057	C ₂₄ H ₃₂ N ₂ O • HCl	2748623-68-7
1	G4	4'-methyl Fentanyl (hydrochloride)	21914	C ₂₃ H ₃₀ N ₂ O • HCl	1071703-97-3
1	G5	<i>para</i> -Fluoroacetyl fentanyl (hydrochloride)	23017	C ₂₁ H ₂₅ FN ₂ O • HCl	104818-37-3
1	G6	β-Hydroxythiofentanyl (hydrochloride)	17421	C ₂₀ H ₂₆ N ₂ O ₂ S • HCl	2320475-89-4
1	G7	Pivaloyl fentanyl (hydrochloride)	22991	C ₂₄ H ₃₂ N ₂ O • HCl	2748409-42-7
1	G8	<i>para</i> -Methylfentanyl (hydrochloride)	20038	C ₂₃ H ₃₀ N ₂ O • HCl	1807-12-1
1	G9	Furanyl fentanyl 3-furancarboxamide isomer (hydrochloride)	21213	C ₂₄ H ₂₆ N ₂ O ₂ • HCl	2306823-47-0
1	G10	<i>para</i> -chloro Acrylfentanyl (hydrochloride)	23338	C ₂₂ H ₂₅ ClN ₂ O • HCl	2306823-58-3
1	G11	Valeryl fentanyl (hydrochloride)	18934	C ₂₄ H ₃₂ N ₂ O • HCl	117332-91-9
1	G12	<i>Unused</i>			
1	H1	<i>Unused</i>			
1	H2	(±)- <i>trans</i> -3-methyl Fentanyl (hydrochloride)	9002482	C ₂₃ H ₃₀ N ₂ O • HCl	78995-09-2
1	H3	<i>meta</i> -Fluorofentanyl (hydrochloride)	19424	C ₂₂ H ₂₇ FN ₂ O • HCl	2309383-05-7
1	H4	<i>para</i> -fluoro Methoxyacetyl fentanyl (hydrochloride)	21741	C ₂₂ H ₂₇ FN ₂ O ₂ • HCl	2306823-76-5
1	H5	Isovaleryl fentanyl (hydrochloride)	22990	C ₂₄ H ₃₂ N ₂ O • HCl	2583550-98-3
1	H6	<i>ortho</i> -Methylfentanyl (hydrochloride)	22634	C ₂₃ H ₃₀ N ₂ O • HCl	1443-53-4
1	H7	<i>para</i> -Fluorofentanyl (hydrochloride)	15496	C ₂₂ H ₂₇ FN ₂ O • HCl	117332-92-0
1	H8	Tetrahydrofuran fentanyl 3-tetrahydrofurancarboxamide (hydrochloride)	22664	C ₂₄ H ₃₀ N ₂ O ₂ • HCl	2306827-13-2
1	H9	(±)- <i>cis</i> -3-methyl Butyryl fentanyl (hydrochloride)	20817	C ₂₄ H ₃₂ N ₂ O • HCl	88641-20-7
1	H10	β-methyl Fentanyl (hydrochloride)	9002860	C ₂₃ H ₃₀ N ₂ O • HCl	1443-43-2
1	H11	(±)- <i>cis</i> -3-methyl Thiofentanyl (hydrochloride)	22240	C ₂₁ H ₂₈ N ₂ OS • HCl	131690-45-4
1	H12	<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
2	A1	<i>Unused</i>			
2	A2	<i>para</i> -methoxy Acrylfentanyl (hydrochloride)	23192	C ₂₃ H ₂₈ N ₂ O ₂ • HCl	2748409-46-1
2	A3	<i>ortho</i> -methoxy Butyryl fentanyl (hydrochloride)	22035	C ₂₄ H ₃₂ N ₂ O ₂ • HCl	2749433-35-8
2	A4	<i>meta</i> -methyl Methoxyacetyl fentanyl (hydrochloride)	22978	C ₂₃ H ₃₀ N ₂ O ₂ • HCl	2748304-70-1
2	A5	<i>ortho</i> -Fluorofentanyl (hydrochloride)	19425	C ₂₂ H ₂₇ FN ₂ O • HCl	2306825-26-1
2	A6	Remifentanil (hydrochloride)	19291	C ₂₀ H ₂₈ N ₂ O ₅ • HCl	132539-07-2
2	A7	<i>para</i> -Chlorobutyryl fentanyl (hydrochloride)	23015	C ₂₃ H ₂₉ ClN ₂ O • HCl	2306823-72-1
2	A8	Ethoxyacetyl fentanyl (hydrochloride)	22992	C ₂₃ H ₃₀ N ₂ O ₂ • HCl	101365-55-3
2	A9	(±)- <i>trans</i> -3-methyl Thiofentanyl (hydrochloride)	22241	C ₂₁ H ₂₈ N ₂ OS • HCl	2748623-40-5
2	A10	<i>meta</i> -fluoro Methoxyacetyl fentanyl (hydrochloride)	21740	C ₂₂ H ₂₇ FN ₂ O ₂ • HCl	2306825-32-9
2	A11	<i>para</i> -fluoro Furanyl fentanyl (hydrochloride)	22750	C ₂₄ H ₂₅ FN ₂ O ₂ • HCl	2748623-32-5
2	A12	<i>Unused</i>			
2	B1	<i>Unused</i>			
2	B2	FIBF (hydrochloride)	19313	C ₂₃ H ₂₉ FN ₂ O • HCl	2309383-06-8
2	B3	<i>para</i> -fluoro Cyclopentyl fentanyl (hydrochloride)	23114	C ₂₅ H ₃₁ FN ₂ O • HCl	2749394-71-4
2	B4	<i>ortho</i> -methyl Phenyl fentanyl (hydrochloride)	23652	C ₂₇ H ₃₀ N ₂ O • HCl	2748290-11-9
2	B5	<i>para</i> -fluoro Valeryl fentanyl (hydrochloride)	23085	C ₂₄ H ₃₁ FN ₂ O • HCl	2749618-68-4
2	B6	<i>para</i> -methyl Methoxyacetyl fentanyl (hydrochloride)	22979	C ₂₃ H ₃₀ N ₂ O ₂ • HCl	2748409-27-8
2	B7	Cyclohexyl fentanyl (hydrochloride)	22390	C ₂₆ H ₃₄ N ₂ O • HCl	2309383-14-8
2	B8	<i>para</i> -methoxy Valeryl fentanyl (hydrochloride)	23471	C ₂₅ H ₃₄ N ₂ O ₂ • HCl	2749910-53-8
2	B9	Sufentanil	15917	C ₂₂ H ₃₀ N ₂ O ₂ S	56030-54-7
2	B10	α'-methoxy Fentanyl (hydrochloride)	23013	C ₂₃ H ₃₀ N ₂ O ₂ • HCl	101365-58-6
2	B11	<i>para</i> -chloro Furanyl fentanyl	22807	C ₂₄ H ₂₅ ClN ₂ O ₂	2748496-77-5
2	B12	<i>Unused</i>			
2	C1	<i>Unused</i>			
2	C2	<i>para</i> -methyl Furanyl fentanyl (hydrochloride)	22713	C ₂₅ H ₂₈ N ₂ O ₂ • HCl	2306827-94-9
2	C3	4'-fluoro, <i>para</i> -fluoro (±)- <i>trans</i> -3-methyl Fentanyl (hydrochloride)	24503	C ₂₃ H ₂₈ F ₂ N ₂ O • HCl	
2	C4	<i>para</i> -methoxy Tetrahydrofuran fentanyl (hydrochloride)	23472	C ₂₅ H ₃₂ N ₂ O ₃ • HCl	2749299-32-7
2	C5	<i>para</i> -methoxy Furanyl fentanyl (hydrochloride)	23197	C ₂₅ H ₂₈ N ₂ O ₃ • HCl	2748591-41-3
2	C6	<i>meta</i> -methyl Furanyl fentanyl (hydrochloride)	22844	C ₂₅ H ₂₈ N ₂ O ₂ • HCl	2748409-34-7
2	C7	<i>para</i> -chloro Cyclopentyl fentanyl (hydrochloride)	23341	C ₂₅ H ₃₁ ClN ₂ O • HCl	2748485-30-3
2	C8	Norcarfentanil (hydrochloride)	19859	C ₁₆ H ₂₂ N ₂ O ₃ • HCl	61085-87-8
2	C9	<i>para</i> -fluoro Tetrahydrofuran fentanyl (hydrochloride)	22752	C ₂₄ H ₂₉ FN ₂ O ₂ • HCl	2748485-19-8
2	C10	2,2,3,3-tetramethyl-Cyclopropyl fentanyl (hydrochloride)	22744	C ₂₇ H ₃₆ N ₂ O • HCl	2309383-12-6
2	C11	Furanylethyl fentanyl (hydrochloride)	20819	C ₂₀ H ₂₆ N ₂ O ₂ • HCl	1443-49-8
2	C12	<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
2	D1	<i>Unused</i>			
2	D2	Benzyl fentanyl (hydrochloride)	19883	C ₂₁ H ₂₆ N ₂ O • HCl	5156-58-1
2	D3	Thienyl fentanyl (hydrochloride)	20785	C ₁₉ H ₂₄ N ₂ O ₂ • HCl	117332-93-1
2	D4	Acrylfentanyl (hydrochloride)	19312	C ₂₂ H ₂₆ N ₂ O • HCl	79279-03-1
2	D5	<i>ortho</i> -isopropyl Furanyl fentanyl	22804	C ₂₇ H ₃₂ N ₂ O ₂	2748591-50-4
2	D6	<i>para</i> -chloro Valeryl fentanyl (hydrochloride)	23058	C ₂₄ H ₃₁ ClN ₂ O • HCl	2748289-84-9
2	D7	Thiofentanyl (hydrochloride)	20786	C ₂₀ H ₂₆ N ₂ O ₂ • HCl	79278-88-9
2	D8	N-benzyl Furanyl norfentanyl (hydrochloride)	25239	C ₂₃ H ₂₄ N ₂ O ₂ • HCl	2711605-09-1
2	D9	N-methyl Norcarfentanil (hydrochloride)	43943	C ₁₇ H ₂₄ N ₂ O ₃ • HCl	2748623-59-6
2	D10	Benzyl Acrylfentanyl (hydrochloride)	22660	C ₂₁ H ₂₄ N ₂ O • HCl	2306823-77-6
2	D11	2'-fluoro <i>ortho</i> -Fluorofentanyl (hydrochloride)	23400	C ₂₂ H ₂₆ F ₂ N ₂ O • HCl	2748343-88-4
2	D12	<i>Unused</i>			
2	E1	<i>Unused</i>			
2	E2	Acetyl fentanyl (hydrochloride)	ISO00128	C ₂₁ H ₂₆ N ₂ O • HCl	117332-89-5
2	E3	U-47700	18596	C ₁₆ H ₂₂ Cl ₂ N ₂ O	82657-23-6
2	E4	U-49900	21098	C ₁₈ H ₂₆ Cl ₂ N ₂ O	67579-76-4
2	E5	<i>meta</i> -methyl Cyclopropyl fentanyl (hydrochloride)	23044	C ₂₄ H ₃₀ N ₂ O • HCl	2748301-54-2
2	E6	<i>para</i> -methyl Cyclopentyl fentanyl (hydrochloride)	23041	C ₂₆ H ₃₄ N ₂ O • HCl	2748409-68-7
2	E7	<i>para</i> -methoxy Acetyl fentanyl (hydrochloride)	23190	C ₂₂ H ₂₈ N ₂ O ₂ • HCl	2749394-67-8
2	E8	Furanyl norfentanyl (hydrochloride)	19429	C ₁₆ H ₁₈ N ₂ O ₂ • HCl	2748343-79-3
2	E9	Fentanyl Methyl Carbamate	25678	C ₂₁ H ₂₆ N ₂ O ₂	2706537-37-1
2	E10	Cyclopentenyl fentanyl (hydrochloride)	23010	C ₂₅ H ₃₀ N ₂ O • HCl	2749298-75-5
2	E11	U-48800 (hydrochloride)	22278	C ₁₇ H ₂₄ Cl ₂ N ₂ O • HCl	2749298-67-5
2	E12	<i>Unused</i>			
2	F1	<i>Unused</i>			
2	F2	4'-Fluorofentanyl (hydrochloride)	25877	C ₂₂ H ₂₇ FN ₂ O • HCl	2748343-99-7
2	F3	β-hydroxy Fentanyl (hydrochloride)	20789	C ₂₂ H ₂₈ N ₂ O ₂ • HCl	1473-95-6
2	F4	MT-45 (hydrochloride)	14082	C ₂₄ H ₃₂ N ₂ • 2HCl	57314-55-3
2	F5	<i>para</i> -methyl Butyryl fentanyl (hydrochloride)	23048	C ₂₄ H ₃₂ N ₂ O • HCl	
2	F6	Norsufentanil	9001142	C ₁₆ H ₂₄ N ₂ O ₂	61086-18-8
2	F7	β-Hydroxythioacetylfentanyl	24280	C ₁₉ H ₂₄ N ₂ O ₂ S	2748485-33-6
2	F8	Seneciyl fentanyl	25440	C ₂₄ H ₃₀ N ₂ O	2630378-28-6
2	F9	Isopropyl U-47700	24263	C ₁₈ H ₂₆ Cl ₂ N ₂ O	2748319-16-4
2	F10	N,N-Dimethylamido-despropionyl fentanyl	23750	C ₂₂ H ₂₉ N ₃ O	1443-50-1
2	F11	AH 7921	12036	C ₁₆ H ₂₂ Cl ₂ N ₂ O	55154-30-8
2	F12	<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
2	G1	<i>Unused</i>			
2	G2	<i>para</i> -methoxy Methoxyacetyl fentanyl (hydrochloride)	23199	C ₂₃ H ₃₀ N ₂ O ₃ • HCl	2748289-90-7
2	G3	Heptanoyl fentanyl (hydrochloride)	25488	C ₂₆ H ₃₆ N ₂ O • HCl	2749326-85-8
2	G4	Tetrahydrothiophene fentanyl	23399	C ₂₄ H ₃₀ N ₂ OS	2748301-41-7
2	G5	4-Phenyl fentanyl (hydrochloride)	20790	C ₂₈ H ₃₂ N ₂ O • HCl	2748301-34-8
2	G6	Phenoxyacetyl fentanyl (hydrochloride)	25604	C ₂₇ H ₃₀ N ₂ O ₂ • HCl	2749299-63-4
2	G7	2-fluoro MT-45 (hydrochloride)	25288	C ₂₄ H ₃₁ FN ₂ • 2HCl	2748592-60-9
2	G8	Furanyl fentanyl (hydrochloride)	18705	C ₂₄ H ₂₆ N ₂ O ₂ • HCl	101365-56-4
2	G9	2,3-seco-Fentanyl (hydrochloride)	25092	C ₂₂ H ₃₀ N ₂ O • HCl	2748592-46-1
2	G10	N-(3-ethylindole) Norfentanyl	24721	C ₂₄ H ₂₉ N ₃ O	58399-46-5
2	G11	Hexanoyl fentanyl (hydrochloride)	23751	C ₂₅ H ₃₄ N ₂ O • HCl	2583550-67-6
2	G12	<i>Unused</i>			
2	H1	<i>Unused</i>			
2	H2	<i>Unused</i>			
2	H3	<i>Unused</i>			
2	H4	<i>Unused</i>			
2	H5	<i>Unused</i>			
2	H6	<i>Unused</i>			
2	H7	<i>Unused</i>			
2	H8	<i>Unused</i>			
2	H9	<i>Unused</i>			
2	H10	<i>Unused</i>			
2	H11	<i>Unused</i>			
2	H12	<i>Unused</i>			
3	A1	<i>Unused</i>			
3	A2	U-48520	26274	C ₁₆ H ₂₃ ClN ₂ O	67579-11-7
3	A3	3'-methyl Acetyl fentanyl (hydrochloride)	25474	C ₂₂ H ₂₈ N ₂ O • HCl	1071704-02-3
3	A4	4-phenyl U-51754 (hydrochloride)	24980	C ₂₃ H ₃₀ N ₂ O • HCl	67197-94-8
3	A5	2'-methyl Acetyl fentanyl (hydrochloride)	25473	C ₂₂ H ₂₈ N ₂ O • HCl	1071703-90-6
3	A6	(±)- <i>cis</i> -3-methyl Norfentanyl	22617	C ₁₅ H ₂₂ N ₂ O	33794-42-2
3	A7	<i>meta</i> -methyl Acetyl fentanyl (hydrochloride)	23997	C ₂₂ H ₂₈ N ₂ O • HCl	1443-51-2
3	A8	U-51754 (hydrochloride)	20285	C ₁₇ H ₂₄ Cl ₂ N ₂ O • HCl	2748595-55-1
3	A9	(±)- <i>cis</i> -Isofentanyl (hydrochloride)	24102	C ₂₂ H ₂₈ N ₂ O • HCl	79278-68-5
3	A10	U-47931E	20530	C ₁₅ H ₂₁ BrN ₂ O	67579-24-2
3	A11	U-48753E (maleate)	20734	C ₁₆ H ₂₂ Cl ₂ N ₂ O • C ₄ H ₄ O ₄	67450-45-7
3	A12	<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
3	B1	<i>Unused</i>			
3	B2	<i>meta</i> -fluoro Furanyl fentanyl (hydrochloride)	26805	C ₂₄ H ₂₅ FN ₂ O ₂ • HCl	
3	B3	3,4-Methylenedioxy U-47700 (hydrochloride)	22753	C ₁₇ H ₂₄ N ₂ O ₃ • HCl	
3	B4	3,4-Ethylenedioxy U-51754 (hydrochloride)	23747	C ₁₉ H ₂₈ N ₂ O ₃ • HCl	2748623-92-7
3	B5	3,4-Ethylenedioxy U-47700 (hydrochloride)	23752	C ₁₈ H ₂₆ N ₂ O ₃ • HCl	2749619-09-6
3	B6	N-methyl U-47931E	24286	C ₁₆ H ₂₃ BrN ₂ O	75570-38-6
3	B7	2'-methyl Fentanyl (hydrochloride)	25475	C ₂₃ H ₃₀ N ₂ O • HCl	2749433-61-0
3	B8	Propyl U-47700	24434	C ₁₈ H ₂₆ Cl ₂ N ₂ O	2749433-76-7
3	B9	N-desmethyl U-47700	20279	C ₁₅ H ₂₀ Cl ₂ N ₂ O	67579-73-1
3	B10	3'-methyl Fentanyl (hydrochloride)	25476	C ₂₃ H ₃₀ N ₂ O • HCl	1082721-49-0
3	B11	N,N-didesmethyl U-47700 (trifluoroacetate salt)	20295	C ₁₄ H ₁₈ Cl ₂ N ₂ O • CF ₃ COOH	2708579-41-1
3	B12	<i>Unused</i>			
3	C1	<i>Unused</i>			
3	C2	Remifentanil Acid (trifluoroacetate salt)	21926	C ₁₉ H ₂₆ N ₂ O ₅ • CF ₃ COOH	132875-69-5
3	C3	N-Benzyl phenyl norfentanyl (hydrochloride)	24249	C ₂₅ H ₂₆ N ₂ O • HCl	2748592-16-5
3	C4	<i>para</i> -fluoro 4-ANBP	25960	C ₁₈ H ₂₁ FN ₂	131587-27-4
3	C5	Despropionyl <i>para</i> -Fluorofentanyl	20104	C ₁₉ H ₂₃ FN ₂	122861-41-0
3	C6	Despropionyl <i>meta</i> -Methylfentanyl	25744	C ₂₀ H ₂₆ N ₂	99918-45-3
3	C7	N-benzyl <i>para</i> -fluoro Norfentanyl (hydrochloride)	25542	C ₂₁ H ₂₅ FN ₂ O • HCl	2748624-34-0
3	C8	Despropionyl 2'-fluoro <i>ortho</i> -Fluorofentanyl	24736	C ₁₉ H ₂₂ F ₂ N ₂	
3	C9	U-50488 (hydrochloride)	19363	C ₁₉ H ₂₆ Cl ₂ N ₂ O • HCl	67197-96-0
3	C10	N-methyl Norfentanyl (hydrochloride)	24446	C ₁₅ H ₂₂ N ₂ O • HCl	24775-71-1
3	C11	N-benzyl <i>para</i> -fluoro Cyclopropyl norfentanyl (hydrochloride)	25238	C ₂₂ H ₂₅ FN ₂ O • HCl	
3	C12	<i>Unused</i>			
3	D1	<i>Unused</i>			
3	D2	Despropionyl <i>ortho</i> -Fluorofentanyl	20126	C ₁₉ H ₂₃ FN ₂	864422-91-3
3	D3	4-Piperidone (hydrochloride hydrate)	21961	C ₅ H ₉ NO • HCl [H ₂ O]	40064-34-4
3	D4	2',4'-dimethoxy Fentanyl (hydrochloride)	28069	C ₂₄ H ₃₂ N ₂ O ₃ • HCl	
3	D5	O-Desmethyl- <i>cis</i> -tramadol (hydrochloride)	13976	C ₁₅ H ₂₃ NO ₂ • HCl	185453-02-5
3	D6	3',4'-dimethoxy Fentanyl (hydrochloride)	28072	C ₂₄ H ₃₂ N ₂ O ₃ • HCl	
3	D7	2-methyl AP-237 (hydrochloride)	26485	C ₁₈ H ₂₆ N ₂ O • HCl	98608-59-4
3	D8	Tianeptine (sodium salt)	25950	C ₂₁ H ₂₄ ClN ₂ O ₄ S • Na	30123-17-2
3	D9	Piperidylthiambutene (hydrochloride)	26335	C ₁₇ H ₂₁ NS ₂ • HCl	64037-51-0
3	D10	<i>para</i> -Bromofentanyl	27536	C ₂₂ H ₂₇ BrN ₂ O	117994-23-7
3	D11	<i>para</i> -Toluoyl fentanyl (hydrochloride)	27537	C ₂₇ H ₃₀ N ₂ O • HCl	
3	D12	<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
3	E1	<i>Unused</i>			
3	E2	<i>meta</i> -fluoro Acrylfentanyl	21931	C ₂₂ H ₂₅ FN ₂ O	2748592-37-0
3	E3	Metonitazene (citrate)	31303	C ₂₁ H ₂₆ N ₄ O ₃ • C ₆ H ₈ O ₇	
3	E4	4-Anilino-1-benzylpiperidine	20086	C ₁₈ H ₂₂ N ₂	1155-56-2
3	E5	4-Anilino-1-Boc-piperidine	29119	C ₁₆ H ₂₄ N ₂ O ₂	125541-22-2
3	E6	<i>para</i> -chloro Furanyl fentanyl 3-furancarboxamide	22843	C ₂₄ H ₂₅ ClN ₂ O ₂	2749618-77-5
3	E7	Thiophene fentanyl 3-thiophenecarboxamide (hydrochloride)	23009	C ₂₄ H ₂₆ N ₂ OS • HCl	2748624-69-1
3	E8	3'-Fluorofentanyl (hydrochloride)	25884	C ₂₂ H ₂₇ FN ₂ O • HCl	2749434-50-0
3	E9	<i>ortho</i> -fluoro Valeryl fentanyl (hydrochloride)	23083	C ₂₄ H ₃₁ FN ₂ O • HCl	2748624-06-6
3	E10	4-methyl Fentanyl (hydrochloride)	21732	C ₂₃ H ₃₀ N ₂ O • HCl	
3	E11	Cyclopropaneacetyl fentanyl (hydrochloride)	26259	C ₂₄ H ₃₀ N ₂ O • HCl	
3	E12	<i>Unused</i>			
3	F1	<i>Unused</i>			
3	F2	Etonitazene	21824	C ₂₂ H ₂₈ N ₄ O ₃	911-65-9
3	F3	<i>para</i> -Chloroacetyl fentanyl (hydrochloride)	23016	C ₂₁ H ₂₅ ClN ₂ O • HCl	
3	F4	2-fluoro Viminol	27233	C ₂₁ H ₃₁ FN ₂ O	63880-43-3
3	F5	<i>para</i> -hydroxy Butyryl fentanyl	26053	C ₂₃ H ₃₀ N ₂ O ₂	2731038-58-5
3	F6	2'-Fluorofentanyl (hydrochloride)	25883	C ₂₂ H ₂₇ FN ₂ O • HCl	2749327-14-6
3	F7	AP-237 (hydrochloride)	26484	C ₁₇ H ₂₄ N ₂ O • HCl	17730-82-4
3	F8	<i>meta</i> -methoxy Furanyl fentanyl	26050	C ₂₅ H ₂₈ N ₂ O ₃	2749434-68-0
3	F9	Brorphine (hydrochloride)	27816	C ₂₀ H ₂₂ BrN ₃ O • HCl	2707204-49-5
3	F10	2,3-Benzodioxole fentanyl	26258	C ₂₇ H ₂₈ N ₂ O ₃	
3	F11	3'-fluoro <i>ortho</i> -Fluorofentanyl (hydrochloride)	23971	C ₂₂ H ₂₆ F ₂ N ₂ O • HCl	2748624-79-3
3	F12	<i>Unused</i>			
3	G1	<i>Unused</i>			
3	G2	Tigloyl fentanyl	26261	C ₂₄ H ₃₀ N ₂ O	
3	G3	2',3'-dimethoxy Fentanyl (hydrochloride)	28068	C ₂₄ H ₃₂ N ₂ O ₃ • HCl	
3	G4	N-Benzyl-4-piperidone	21962	C ₁₂ H ₁₅ NO	3612-20-2
3	G5	2',6'-dimethoxy Fentanyl (hydrochloride)	28071	C ₂₄ H ₃₂ N ₂ O ₃ • HCl	
3	G6	NPP	20528	C ₁₃ H ₁₇ NO	39742-60-4
3	G7	3',5'-dimethoxy Fentanyl (hydrochloride)	28073	C ₂₄ H ₃₂ N ₂ O ₃ • HCl	
3	G8	AP-238 (hydrochloride)	31128	C ₁₈ H ₂₆ N ₂ O • HCl	2749171-05-7
3	G9	4-Anilinopiperidine (hydrochloride)	20085	C ₁₁ H ₁₆ N ₂ • 2HCl	99918-43-1
3	G10	Isotonitazene	27255	C ₂₃ H ₃₀ N ₄ O ₃	14188-81-9
3	G11	<i>meta</i> -fluoro Valeryl fentanyl (hydrochloride)	23084	C ₂₄ H ₃₁ FN ₂ O • HCl	
3	G12	<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
3	H1	Unused			
3	H2	Unused			
3	H3	Unused			
3	H4	Unused			
3	H5	Unused			
3	H6	Unused			
3	H7	Unused			
3	H8	Unused			
3	H9	Unused			
3	H10	Unused			
3	H11	Unused			
3	H12	Unused			
4	A1	Unused			
4	A2	N-(MDA) Fentanyl (hydrochloride)	28994	C ₂₄ H ₃₀ N ₂ O ₃ • HCl	
4	A3	N-(Phentermine) Fentanyl (hydrochloride)	28995	C ₂₄ H ₃₂ N ₂ O • HCl	
4	A4	N-(6-APB) Fentanyl	28968	C ₂₅ H ₃₀ N ₂ O ₂	
4	A5	N-(6-APDB) Fentanyl	28969	C ₂₅ H ₃₂ N ₂ O ₂	
4	A6	N-(2-APB) Fentanyl	28967	C ₂₅ H ₃₀ N ₂ O ₂	
4	A7	2',5'-dimethoxy Fentanyl (hydrochloride)	28070	C ₂₄ H ₃₂ N ₂ O ₃ • HCl	
4	A8	N-(2C-E) Fentanyl (hydrochloride)	28973	C ₂₆ H ₃₆ N ₂ O ₃ • HCl	
4	A9	N-(2C-B) Fentanyl (hydrochloride)	28970	C ₂₄ H ₃₁ BrN ₂ O ₃ • HCl	
4	A10	N-(2C-C) Fentanyl (hydrochloride)	28971	C ₂₄ H ₃₁ ClN ₂ O ₃ • HCl	
4	A11	N-(2C-P) Fentanyl (hydrochloride)	28978	C ₂₇ H ₃₈ N ₂ O ₃ • HCl	
4	A12	Unused			
4	B1	Unused			
4	B2	N-(2C-T-7) Fentanyl (hydrochloride)	28982	C ₂₇ H ₃₈ N ₂ O ₃ S • HCl	
4	B3	N-(2C-D) Fentanyl (hydrochloride)	28972	C ₂₅ H ₃₄ N ₂ O ₃ • HCl	
4	B4	N-(2C-G) Fentanyl (hydrochloride)	28974	C ₂₆ H ₃₆ N ₂ O ₃ • HCl	
4	B5	N-(2C-I) Fentanyl (hydrochloride)	28975	C ₂₄ H ₃₁ IN ₂ O ₃ • HCl	
4	B6	N-(2C-iP) Fentanyl (hydrochloride)	28976	C ₂₇ H ₃₈ N ₂ O ₃ • HCl	
4	B7	N-(2C-N) Fentanyl (hydrochloride)	28977	C ₂₄ H ₃₁ N ₃ O ₅ • HCl	
4	B8	N-(2C-T) Fentanyl	28979	C ₂₅ H ₃₄ N ₂ O ₃ S	
4	B9	N-(2C-T-2) Fentanyl (hydrochloride)	28980	C ₂₆ H ₃₆ N ₂ O ₃ S • HCl	
4	B10	N-(2C-T-4) Fentanyl (hydrochloride)	28981	C ₂₇ H ₃₈ N ₂ O ₃ S • HCl	
4	B11	N-(2C-TFM) Fentanyl (hydrochloride)	28983	C ₂₅ H ₃₁ F ₃ N ₂ O ₃ • HCl	
4	B12	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
4	C1	<i>Unused</i>			
4	C2	N-(2C-B-FLY) Fentanyl (hydrochloride)	28984	$C_{26}H_{31}BrN_2O_3 \cdot HCl$	
4	C3	N-(3C-B-FLY) Fentanyl (hydrochloride)	28985	$C_{27}H_{33}BrN_2O_3 \cdot HCl$	
4	C4	N-(3,4,5-TMA) Fentanyl (hydrochloride)	28997	$C_{26}H_{36}N_2O_4 \cdot HCl$	
4	C5	N-(2,5-DMA) Fentanyl (hydrochloride)	28993	$C_{25}H_{34}N_2O_3 \cdot HCl$	
4	C6	N-(DOC) Fentanyl (hydrochloride)	28988	$C_{25}H_{33}ClN_2O_3 \cdot HCl$	
4	C7	N-(DOB) Fentanyl (hydrochloride)	28986	$C_{25}H_{33}BrN_2O_3 \cdot HCl$	
4	C8	N-(DOI) Fentanyl (hydrochloride)	28990	$C_{25}H_{33}IN_2O_3 \cdot HCl$	
4	C9	N-(DOM) Fentanyl (hydrochloride)	28991	$C_{26}H_{36}N_2O_3 \cdot HCl$	
4	C10	N-(DOET) Fentanyl (hydrochloride)	28989	$C_{27}H_{38}N_2O_3 \cdot HCl$	
4	C11	N-(DOBU) Fentanyl (hydrochloride)	28987	$C_{29}H_{42}N_2O_3 \cdot HCl$	
4	C12	<i>Unused</i>			
4	D1	<i>Unused</i>			
4	D2	<i>Unused</i>			
4	D3	<i>Unused</i>			
4	D4	<i>Unused</i>			
4	D5	<i>Unused</i>			
4	D6	<i>Unused</i>			
4	D7	<i>Unused</i>			
4	D8	<i>Unused</i>			
4	D9	<i>Unused</i>			
4	D10	<i>Unused</i>			
4	D11	<i>Unused</i>			
4	D12	<i>Unused</i>			
4	E1	<i>Unused</i>			
4	E2	<i>Unused</i>			
4	E3	<i>Unused</i>			
4	E4	<i>Unused</i>			
4	E5	<i>Unused</i>			
4	E6	<i>Unused</i>			
4	E7	<i>Unused</i>			
4	E8	<i>Unused</i>			
4	E9	<i>Unused</i>			
4	E10	<i>Unused</i>			
4	E11	<i>Unused</i>			
4	E12	<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
4	F1	Unused			
4	F2	Unused			
4	F3	Unused			
4	F4	Unused			
4	F5	Unused			
4	F6	Unused			
4	F7	Unused			
4	F8	Unused			
4	F9	Unused			
4	F10	Unused			
4	F11	Unused			
4	F12	Unused			
4	G1	Unused			
4	G2	Unused			
4	G3	Unused			
4	G4	Unused			
4	G5	Unused			
4	G6	Unused			
4	G7	Unused			
4	G8	Unused			
4	G9	Unused			
4	G10	Unused			
4	G11	Unused			
4	G12	Unused			
4	H1	Unused			
4	H2	Unused			
4	H3	Unused			
4	H4	Unused			
4	H5	Unused			
4	H6	Unused			
4	H7	Unused			
4	H8	Unused			
4	H9	Unused			
4	H10	Unused			
4	H11	Unused			
4	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