



Fentanyl Analog Screening Kit

Item No. 9003237

www.caymanchem.com

Customer Service 800.364.9897

Technical Support 888.526.5351

1180 E. Ellsworth Rd • Ann Arbor, MI • USA

TABLE OF CONTENTS

GENERAL INFORMATION	3	Safety Data
	3	If You Have Problems
INSTRUCTIONS	4	About this Kit
	4	Instructions for Use
	4	Example Reconstitution Procedure
	5	Electronic Resources
	5	Acknowledgement of FAS Kit Use
PLATE LAYOUT	6	Plate 1
	8	Plate 2
RESOURCES	10	Notes
	11	Warranty and Limitation of Remedy

GENERAL INFORMATION

Kit will arrive packaged at -20°C.

If any of the items listed are damaged or missing, please contact our Customer Service department at (800) 364-9897 or (734) 971-3335. We cannot accept any returns without prior authorization.



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.

If You Have Problems

Technical Service Contact Information

Phone: 888-526-5351 (USA and Canada only) or 734-975-3888
Fax: 734-971-3640
Email: faskit@caymanchem.com

In order for our staff to assist you quickly and efficiently, please be ready to supply the lot number of the kit (found on the outside of the box).

INSTRUCTIONS

About this Kit

The Centers for Disease Control and Prevention (CDC) has developed Traceable Opioid Material* Kits (TOM Kits*) to support laboratory detection of emerging opioids in the US. The CDC has contracted Cayman Chemical to manufacture and distribute the Fentanyl Analog Screening (FAS) Kit. This kit enables labs to screen for 120 synthetic fentanyl compounds.

Instructions for Use

The dried plates included in this kit contain a small amount of glycerol in each well to aid with compound recovery. Recovery in methanol has been validated for all analytes on the dried plates and is the recommended reconstitution solvent for this product. Recovery from other solvents has not been verified.

Example Reconstitution Procedure

1. Uncap each vial to be used. Add 500 μ l methanol to each vial. This will provide 400 μ g/ml (nominal†) of each analyte. *NOTE: Vial capacity is 700 μ l. No more than 500 μ l of solvent should be added.*
2. Recap the vials and place the plate on a plate mixer or vortexer. Plates should be mixed at 400 rpm for 1 hour or vortexed at medium speed for a minimum of 15 minutes to ensure full recovery.
3. Carefully remove the caps for analysis.

* TRACEABLE OPIOID MATERIAL, TOM KITS, and the TOM KITS Logo are marks of the U.S. Department of Health and Human Services

†Analyte mass included per vial is nominal and is not adjusted for salt content or organic purity.

Testing of the kits determined them to be stable for 1 week after reconstitution. Longer periods have not been tested. We recommend plates be stored dry until use. Dry plates are expected to be stable for a minimum of 18 months when stored unopened at the recommended storage conditions. Testing is ongoing to ensure long-term stability of this product.

Electronic Resources

To obtain the Certificate of Analysis, Safety Data Sheets, searchable compound database, and other electronic documents for this kit, please visit www.caymanchem.com/faskitdocuments.

Acknowledgement of FAS Kit Use

Acknowledge the use of Traceable Opioid Material* Kits (TOM Kits*) in presentations, publications, reports, and other communications by using the following citation: "Laboratory findings were made possible, in part, by the Centers for Disease Control and Prevention's design and support of Traceable Opioid Material* Kits."

* TRACEABLE OPIOID MATERIAL, TOM KITS, and the TOM KITS Logo are marks of the U.S. Department of Health and Human Services

PLATE 1												
	1	2	3	4	5	6	7	8	9	10	11	12
A	Unused	20816 <i>α</i> -methyl Acetyl fentanyl (hydrochloride)	ISO00130 Acetyl norfentanyl (hydrochloride)	22805 <i>ortho</i> -methoxy Furanyl fentanyl	22806 <i>ortho</i> -methyl Furanyl fentanyl	9002271 4'-methyl Acetyl fentanyl (hydrochloride)	22802 Thiophene fentanyl (hydrochloride)	23339 <i>para</i> -chloro Cyclobutyl fentanyl (hydrochloride)	24278 N-methyl Cyclopropyl norfentanyl (hydrochloride)	21736 <i>β</i> -methyl Acetyl fentanyl (hydrochloride)	18810 4-ANPP	Unused
B	Unused	23038 <i>para</i> -methyl Acrylfentanyl (hydrochloride)	23043 <i>ortho</i> -methyl Cyclopropyl fentanyl (hydrochloride)	23580 Fentanyl	15899 Norfentanyl	21739 Cyclopropyl fentanyl (hydrochloride)	19428 Butyryl norfentanyl (hydrochloride)	23548 <i>ortho</i> -methyl Acetyl fentanyl (hydrochloride)	22994 <i>α</i> '-methyl Butyryl fentanyl (hydrochloride)	22884 Methacrylfentanyl	20818 <i>α</i> -methyl Butyryl fentanyl (hydrochloride)	Unused
C	Unused	23544 <i>para</i> -fluoro Crotonyl fentanyl	18280 <i>α</i> -methyl Fentanyl (hydrochloride)	22801 Crotonyl fentanyl	23045 <i>para</i> -methyl Cyclopropyl fentanyl (hydrochloride)	20035 <i>para</i> - Methoxyfentanyl (hydrochloride)	22389 Cyclobutyl fentanyl (hydrochloride)	23036 <i>ortho</i> -methyl Acrylfentanyl (hydrochloride)	9002747 (±)- <i>cis</i> -3-methyl Fentanyl (hydrochloride)	18583 Ocfentanil (hydrochloride)	22633 <i>meta</i> - Methylfentanyl (hydrochloride)	Unused
D	Unused	21932 <i>ortho</i> -fluoro Acrylfentanyl (hydrochloride)	20034 <i>para</i> - Chlorofentanyl (hydrochloride)	22749 <i>para</i> -fluoro Cyclopropyl fentanyl (hydrochloride)	14728 Butyryl fentanyl (hydrochloride)	23749 Fentanyl Carbamate	18584 Isobutyryl fentanyl (hydrochloride)	22977 <i>ortho</i> -methyl Methoxyacetyl fentanyl (hydrochloride)	20350 Benzyl Carfentanil (hydrochloride)	20821 <i>α</i> -methyl Thiofentanyl (hydrochloride)	18089 <i>para</i> -methoxy Butyryl fentanyl (hydrochloride)	Unused
E	Unused	23342 <i>para</i> -chloro Cyclopropyl fentanyl (hydrochloride)	19567 <i>ortho</i> - Fluorobutyryl fentanyl (hydrochloride)	21116 <i>para</i> -fluoro Acrylfentanyl	20439 Cyclopentyl fentanyl (hydrochloride)	22551 Phenyl fentanyl (hydrochloride)	20859 Tetrahydrofuran fentanyl (hydrochloride)	20782 Methoxyacetyl fentanyl (hydrochloride)	20207 <i>ortho</i> - Fluoroisobutyryl fentanyl (hydrochloride)	22751 <i>para</i> -fluoro Furanyl fentanyl 3-furancarboxamide (hydrochloride)	17049 <i>para</i> - Fluorobutyryl fentanyl (hydrochloride)	Unused
F	Unused	23054 <i>para</i> -methyl Tetrahydrofuran fentanyl (hydrochloride)	22743 Phenylacetyl fentanyl (hydrochloride)	20018 <i>para</i> - Chloroisobutyryl fentanyl (hydrochloride)	19566 <i>meta</i> - Fluorobutyryl fentanyl (hydrochloride)	23628 <i>ortho</i> -fluoro Furanyl fentanyl (hydrochloride)	20206 <i>meta</i> - Fluoroisobutyryl fentanyl (hydrochloride)	23343 <i>para</i> -chloro Methoxyacetyl fentanyl (hydrochloride)	19292 Alfentanil (hydrochloride)	23858 <i>β</i> '-Phenyl fentanyl	20858 Benzodioxole fentanyl	Unused
G	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused
H	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused

PLATE 2												
	1	2	3	4	5	6	7	8	9	10	11	12
A	Unused	23191 <i>para</i> -methyl Acetyl fentanyl (hydrochloride)	23057 <i>para</i> -methyl Isobutyl fentanyl (hydrochloride)	21914 4'-methyl Fentanyl (hydrochloride)	23017 <i>para</i> -Fluoroacetyl fentanyl (hydrochloride)	17421 β- Hydroxythiofentanyl (hydrochloride)	22991 Pivaloyl fentanyl (hydrochloride)	20038 <i>para</i> - Methylfentanyl (hydrochloride)	21213 Furanyl fentanyl 3-furancarboxamide isomer (hydrochloride)	23338 <i>para</i> -chloro Acrylfentanyl (hydrochloride)	18934 Valeryl fentanyl (hydrochloride)	Unused
B	Unused	9002482 (±)- <i>trans</i> -3-methyl Fentanyl (hydrochloride)	19424 <i>meta</i> - Fluorofentanyl (hydrochloride)	21741 <i>para</i> -fluoro Methoxyacetyl fentanyl (hydrochloride)	22990 Isovaleryl fentanyl (hydrochloride)	22634 <i>ortho</i> - Methylfentanyl (hydrochloride)	15496 <i>para</i> - Fluorofentanyl (hydrochloride)	22664 Tetrahydrofuran fentanyl 3- tetrahydrofurancarbox- amide (hydrochloride)	20817 (±)- <i>cis</i> -3-methyl Butyryl fentanyl (hydrochloride)	9002860 β-methyl Fentanyl (hydrochloride)	22240 (±)- <i>cis</i> -3-methyl Thiofentanyl (hydrochloride)	Unused
C	Unused	23192 <i>para</i> -methoxy Acrylfentanyl (hydrochloride)	22035 <i>ortho</i> -methoxy Butyryl fentanyl (hydrochloride)	22978 <i>meta</i> -methyl Methoxyacetyl fentanyl (hydrochloride)	19425 <i>ortho</i> - Fluorofentanyl (hydrochloride)	19291 Remifentanyl (hydrochloride)	23015 <i>para</i> - Chlorobutyryl fentanyl (hydrochloride)	22992 Ethoxyacetyl fentanyl (hydrochloride)	22241 (±)- <i>trans</i> -3-methyl Thiofentanyl (hydrochloride)	21740 <i>meta</i> -fluoro Methoxyacetyl fentanyl (hydrochloride)	22750 <i>para</i> -fluoro Furanyl fentanyl (hydrochloride)	Unused
D	Unused	19313 FIBF (hydrochloride)	23114 <i>para</i> -fluoro Cyclopentyl fentanyl (hydrochloride)	23652 <i>ortho</i> -methyl Phenyl fentanyl (hydrochloride)	23085 <i>para</i> -fluoro Valeryl fentanyl (hydrochloride)	22979 <i>para</i> -methyl Methoxyacetyl fentanyl (hydrochloride)	22390 Cyclohexyl fentanyl (hydrochloride)	23471 <i>para</i> -methoxy Valeryl fentanyl (hydrochloride)	15917 Sufentanil	23013 α'-methoxy Fentanyl (hydrochloride)	22807 <i>para</i> -chloro Furanyl fentanyl	Unused
E	Unused	22713 <i>para</i> -methyl Furanyl fentanyl (hydrochloride)	24503 4'-fluoro, <i>para</i> - fluoro (±)- <i>trans</i> - 3-methyl Fentanyl (hydrochloride)	23472 <i>para</i> -methoxy Tetrahydrofuran fentanyl (hydrochloride)	23197 <i>para</i> -methoxy Furanyl fentanyl (hydrochloride)	22844 <i>meta</i> -methyl Furanyl fentanyl (hydrochloride)	23341 <i>para</i> -chloro Cyclopentyl fentanyl (hydrochloride)	19859 Norcarfentanil (hydrochloride)	22752 <i>para</i> -fluoro Tetrahydrofuran fentanyl (hydrochloride)	22744 2,2,3,3-tetramethyl- Cyclopropyl fentanyl (hydrochloride)	20819 Furanylethyl fentanyl (hydrochloride)	Unused
F	Unused	19883 Benzyl fentanyl (hydrochloride)	20785 Thienyl fentanyl (hydrochloride)	19312 Acrylfentanyl (hydrochloride)	22804 <i>ortho</i> -isopropyl Furanyl fentanyl (hydrochloride)	23058 <i>para</i> -chloro Valeryl fentanyl (hydrochloride)	20786 Thiofentanyl (hydrochloride)	25239 N-benzyl Furanyl norfentanyl (hydrochloride)	20784 N-methyl Norcarfentanil (hydrochloride)	22660 Benzyl Acrylfentanyl (hydrochloride)	23400 2'-fluoro <i>ortho</i> - Fluorofentanyl (hydrochloride)	Unused
G	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused
H	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused

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.

This document is copyrighted. All rights are reserved. This document may not, in whole or part, be copied, photocopied, reproduced, translated, or reduced to any electronic medium or machine-readable form without prior consent, in writing, from Cayman Chemical Company.

©08/15/2019, Cayman Chemical Company, Ann Arbor, MI, All rights reserved.
Printed in U.S.A.

