



Fentanyl Analog Screening Kit - Emergent Drug Panel

Item No. 9004121

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
	5	Example Reconstitution Procedure
	6	Electronic Resources
	7	Acknowledgement of FAS Kit Use
PLATE LAYOUT	8	Plate 1
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

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 and its expansion pack, Emergent Drug Panel. This expansion pack enables labs to screen for an additional 60 emerging drugs of abuse.

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 µl methanol to each vial. This will provide 400 µg/ml (nominal†) of each analyte. *NOTE: Vial capacity is 700 µl. No more than 500 µl 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.

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 36 months when stored unopened at the recommended storage conditions. Testing is ongoing to ensure long-term stability of this product.

TOM Kits®, TRACEABLE OPIOID MATERIAL®, and TOM Kits® are registered 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.

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."

TOMKits®, TRACEABLE OPIOID MATERIAL®, and TOM Kits® are registered 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	18173 Clonazolam	12063 Etizolam	24481 Flualprazolam	16435 Flubromazolam	18484 Desalkylflurazepam	22665 Bromazolam	18205 Adinazolam	23107 Deschloroetizolam	23110 (±)-Meclonazepam	15157 Flubromazepam	Unused
B	Unused	14901 Pyrazolam	15554 Diclazepam	14260 Bromazepam	23111 Metizolam	27247 8-Aminoclonazolam	9000849 Phenazepam	23524 Fluclozizolam	33169 N-Pyrrolidino Etonitazene	29916 Etodesnitazene (citrate)	29381 Protonitazene (hydrochloride)	Unused
C	Unused	33720 Dipyanone (hydrochloride)	33168 N - Piperidinyl Etonitazene (ci- trate)	30278 Butonitazene	9001103 Eutylone (hydrochloride)	21682 α -Pyrrolidinoiso- hexanophenone (hydrochloride)	18683 Deschloroket- amine (hydrochloride)	14421 3-hydroxy PCP (hydrochloride)	19786 2-fluoro Deschloroketamine (hydrochloride)	11137 3-methoxy PCP (hydrochloride)	31475 3-chloro PCP (hydrochloride)	Unused
D	Unused	20661 α -Ethylamino- hexanophenone (hydrochloride)	9000746 Pentylone (hydrochloride)	9001934 α -Pyrrolidino- hexanophenone (hydrochloride)	16037 N-ethyl Pentylone (hydrochloride)	9001330 3,4-Methylene- dioxy-N- benzylcathinone (hydrochloride)	11136 4-methoxy PCP (hydrochloride)	26097 MDMB-4en- PINACA	26701 N-butyl Pentylone (hydrochloride)	16358 3,4- Methylenedioxy PV8 (hydrochloride)	14280 α -Ethylamino- pentiofenone (hydrochloride)	Unused
E	Unused	14105 deschloro-N- ethyl- Ketamine (hydrochloride)	18326 4'-chloro- α - Pyrrolidinovalero- phenone (hydro- chloride)	9001083 α -Pyrrolidino- pentiofenone (hydrochloride)	11312 4-hydroxy DiPT (hydrochloride)	26877 3-hydroxy PCE (hydrochloride)	20803 5-fluoro MDMB-PICA	22641 Xylazine	26645 4-fluoro MDMB- BUTINACA	18757 ADB-BINACA	23631 (S)-5-fluoro ADB	Unused
F	Unused	31075 4-fluoro MDMB-BUTICA	33205 ADB-4en- PINACA	33820 ADB-HEXINACA	30724 4-fluoro ABUTINACA	29350 ADB-BUTINACA	25006 ACHMINACA	20194 4-cyano CUMYL- BUTINACA	9001960 MMB-FUBINACA	23005 5-fluoro EDMB-PINACA	30725 5-fluoro EDMB-PICA	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

NOTES

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.

