



Emergent Drug Panel - Internal Standard

Item No. 9004330

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 Dilution 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 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) Kits, Emergent Drug Panel (EDP) Kits, and the EDP expansion pack, Emergent Drug Panel – Internal Standard (EDP-IS) Kit.

NOTE: This expansion pack has been prepared in an alternate format from the previous versions of the TOM Kits®, with the analytes presented as methanolic solutions as opposed to dried residues. This expansion pack provides a panel of 24 stable-labeled emergent drugs of abuse corresponding to a subset of compounds included in the FAS and EDP Kits. This expansion pack and the analytes within can be used as internal standards when optimizing analytical instrumentation and detection methods.

Instructions for Use

The EDP-IS plates are provided in solution for ease of use. Each tube well contains a nominal 100 µl of each component at a 2 mg/ml concentration in a solution of methanol with 0.5% glycerol. The preferred solvent for additional dilution is methanol.

Before opening, warm the EDP-IS plate to room temperature while upright to allow the solution within each tube to settle to the bottom. Use proper Personal Protective Equipment (PPE) when handling the tube wells. The plate is ready to use as is, but if a lower concentration and higher volume is preferred, an example dilution procedure is presented below.

TOM Kits®, **TRACEABLE OPIOID MATERIAL®**, and **TOM Kits®** are registered marks of the U.S. Department of Health and Human Services.

Example Dilution Procedure - Dilute in Tube Well

1. Carefully uncap each tube to be used. Add 300 µl methanol to each tube. This will create 400 µl of a 500 µg/ml (nominal†) solution of each analyte. *NOTE: Tube capacity is 700 µl. No more than 500 µl of solvent should be added.*
2. Recap the tube(s) and place the plate on a plate mixer or vortex mixer. Plates and/or tube wells should be well mixed or vortexed to ensure a homogenous solution within the tube.
3. Carefully remove the septa cap(s) for analysis.

We strongly recommend plates be stored upright at -20°C until use. 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.

†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."

PLATE 1												
	1	2	3	4	5	6	7	8	9	10	11	12
A	Unused	27782 Clonazolam-d ₄	38190 Etizolam-d ₄	29433 Flualprazolam-d ₄	32537 Flubromazolam-d ₄	36918 Bromazolam-d ₅	38353 Xylazine-d ₆	32536 Flubromazepam-d ₄	19597 <i>para</i> - Fluorofentanyl-d ₅ (hydrochloride)	31648 Metonitazene-d ₃ (citrate)	29319 Isotonitazene-d ₇	Unused
B	Unused	31566 Brorphine-d ₇	25562 Carfentanil- ¹³ C ₆	38173 N-Pyrrolidino Etonitazene-d ₅	38174 Etodesnitazene-d ₅ (citrate)	31647 2-methyl AP-237-d ₇ (hydrochloride)	38177 Protonitazene-d ₇ (hydrochloride)	31629 Eutylone-d ₅ (hydrochloride)	38168 α-Pyrrolidinoisoh exanophenone-d ₈ (hydrochloride)	38167 3-hydroxy PCP-d ₁₀ (hydrochloride)	38172 α- Pyrrolidinohexan- ophenone-d ₈ (hydrochloride)	Unused
C	Unused	36754 MDMB-4en- PINACA-d ₄	27487 5-fluoro MDMB- PICA-d ₅	36753 (S)-ADB-BINACA- d ₅	37890 4-fluoro MDMB- BUTINACA-d ₄	Unused	Unused	Unused	Unused	Unused	Unused	Unused
D	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused
E	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused
F	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	Unused	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.

©04/17/2024, Cayman Chemical Company, Ann Arbor, MI, All rights reserved. Printed in U.S.A.

