



## Fentanyl Analog Screening Kit - Emergent Panel Version 4

---

Item No. 9003381

[www.caymanchem.com](http://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

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 and its expansion pack, Emergent Panel Version 4. This expansion pack enables labs to screen for an additional 38 synthetic opioid 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 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 18 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](http://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."

**TOM**Kits®, 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 | 20126<br>Despropionyl<br>ortho-<br>Fluorofentanyl | 21961<br>4-Piperidone<br>(hydrochloride<br>hydrate)               | 28069<br>2',4'-dimethoxy<br>Fentanyl<br>(hydrochloride) | 13976<br>O-Desmethyl- <i>cis</i> -<br>tramadol<br>(hydrochloride) | 28072<br>3',4'-dimethoxy<br>Fentanyl<br>(hydrochloride)                    | 26485<br>2-methyl AP-237<br>(hydrochloride)                                   | 25950<br>Tianeptine<br>(sodium salt)              | 26335<br>Piperidylthiambutene<br>(hydrochloride)                     | 27255<br>Isotonitazene                        | 23084<br><i>meta</i> -fluoro<br>Valeryl fentanyl<br>(hydrochloride)    | Unused |
| B | Unused | 21931<br><i>meta</i> -fluoro<br>Acrylfentanyl     | 31303<br>Metonitazene<br>(citrate)                                | 20086<br>4-Anilino-1-<br>benzylpiperidine               | 29119<br>4-Anilino-1-Boc-<br>piperidine                           | 22843<br><i>para</i> -chloro<br>Furanyl fentanyl<br>3-<br>furancarboxamide | 23009<br>Thiophene fentanyl<br>3-thiophenecarbox-<br>amide<br>(hydrochloride) | 25884<br>3'-Fluorofentanyl<br>(hydrochloride)     | 23083<br><i>ortho</i> -fluoro<br>Valeryl fentanyl<br>(hydrochloride) | 21732<br>4-methyl Fentanyl<br>(hydrochloride) | 26259<br>Cyclopropaneacetyl<br>fentanyl<br>(hydrochloride)             | Unused |
| C | Unused | 21824<br>Etonitazene                              | 23016<br><i>para</i> -Chloroacetyl<br>fentanyl<br>(hydrochloride) | 27233<br>2-fluoro Viminol                               | 26053<br><i>para</i> -hydroxy<br>Butyryl fentanyl                 | 25883<br>2'-Fluorofentanyl<br>(hydrochloride)                              | 26484<br>AP-237<br>(hydrochloride)                                            | 26050<br><i>meta</i> -methoxy<br>Furanyl fentanyl | 27816<br>Borophine<br>(hydrochloride)                                | 26258<br>2,3-Benzodioxole<br>fentanyl         | 23971<br>3'-fluoro <i>ortho</i> -<br>Fluorofentanyl<br>(hydrochloride) | Unused |
| D | Unused | 26261<br>Tigloyl fentanyl                         | 28068<br>2',3'-dimethoxy<br>Fentanyl<br>(hydrochloride)           | 21962<br>N-Benzyl-4-<br>piperidone                      | 28071<br>2',6'-dimethoxy<br>Fentanyl<br>(hydrochloride)           | 20528<br>NPP                                                               | 28073<br>3',5'-dimethoxy<br>Fentanyl<br>(hydrochloride)                       | 31128<br>AP-238<br>(hydrochloride)                | 20085<br>4-<br>Anilinopiperidine<br>(hydrochloride)                  | 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 |

## 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.

©04/13/2021, Cayman Chemical Company, Ann Arbor, MI, All rights reserved.  
Printed in U.S.A.

