

# PRODUCT INFORMATION

## GC-MS Drug Standard Mixture 1

Item No. 9002770

**Supplied as:** A solution in acetonitrile

**Fill Volume:** >1 ml

**Storage:** -20°C

**Stability:** ≥7 years

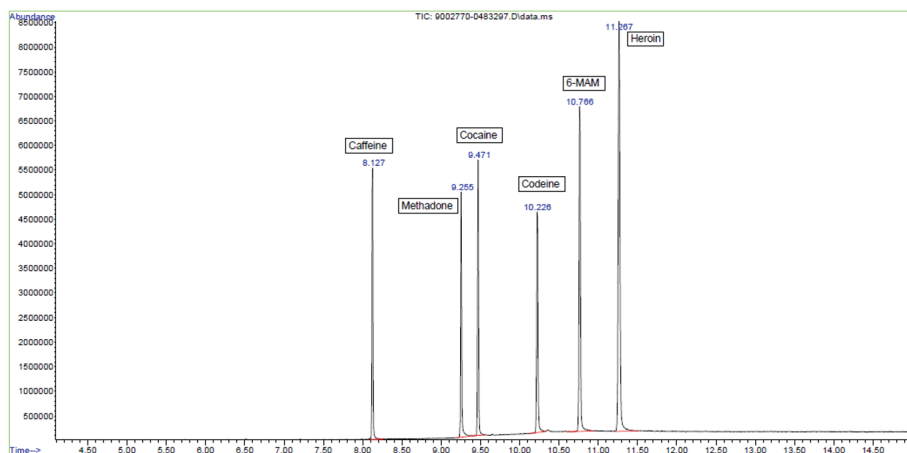
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Description and Contents

GC-MS Drug Standard Mixture 1 (Item No. 9002770) is a mixture of common alkaloids and opioids, as well as a prominent metabolite. The mixture is supplied in an amber ampule in which the headspace has been purged with argon. It is designed to provide authentic standards for analyses using GC-MS or LC-MS. This product is intended for forensic and research applications.

This mixture contains compounds that are regulated as Schedule I, II, and IV compounds in the United States. Included are:

| Item Number | Item Name                           | Concentration |
|-------------|-------------------------------------|---------------|
| 14118       | Caffeine                            | 40 µg/ml      |
| ISO60145    | (±)-Methadone (hydrochloride) (CRM) | 30 µg/ml      |
| ISO60176    | Cocaine (CRM)                       | 30 µg/ml      |
| ISO60140    | Codeine (CRM)                       | 50 µg/ml      |
| 19418       | 6-Acetylmorphine (6-MAM) (CRM)      | 75 µg/ml      |
| ISO60187    | Heroin (CRM)                        | 75 µg/ml      |



#### 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, 07/02/2024

#### 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