

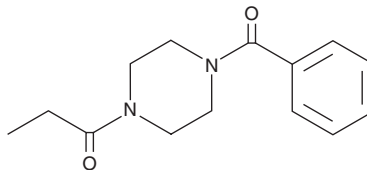
PRODUCT INFORMATION



Sunifiram

Item No. 35018

CAS Registry No.: 314728-85-3
Formal Name: 1-(4-benzoyl-1-piperazinyl)-1-propanone
Synonym: DM235
MF: $C_{14}H_{18}N_2O_2$
FW: 246.3
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 5 years



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

Description

Sunifiram (Item No. 35018) is an analytical reference standard categorized as a nootropic.¹ Sunifiram prevents scopolamine-induced memory deficits in mice. It has been found in seized illicit drug samples.² This product is intended for research and forensic applications.

References

1. Romanelli, M.N., Galeotti, N., Ghelardini, C., *et al.* Pharmacological characterization of DM232 (unifiram) and DM235 (sunifiram), new potent cognition enhancers. *CNS Drug Rev.* **12**(1), 39-52 (2006).
2. Piorunski-Sedlak, K. and Stypulkowska, K. Strategy for identification of new psychoactive substances in illicit samples using attenuated total reflectance infrared spectroscopy. *Forensic Sci. Int.* **312**, 110262 (2020).

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, 07/06/2023

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