

PRODUCT INFORMATION



U-69593

Item No. 13255

Formal Name: N-methyl-2-phenyl-N-(7-(pyrrolidin-1-yl)-1-oxaspiro[4.5]decan-8-yl)acetamide

MF: C₂₂H₃₂N₂O₂

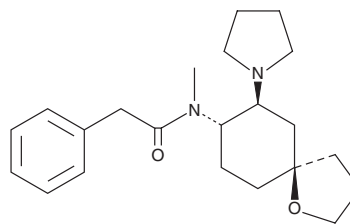
FW: 356.5

Purity: ≥95%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥5 years



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

Description

U-69593 (Item No. 13255) is an analytical reference standard categorized as an opioid.¹ This product is intended for research and forensic applications.

Reference

1. La Regina, A., Petrillo, P., Sbacchi, M., *et al.* Interaction of U-69,593 with mu-, alpha- and kappa-opioid binding sites and its analgesic and intestinal effects in rats. *Life Sci.* **42(3)**, 293-301 (1988).

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, 07/21/2023

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