

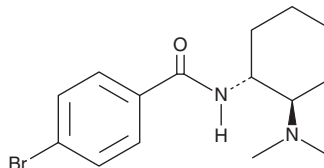
PRODUCT INFORMATION



U-47931E

Item No. 20530

CAS Registry No.: 67579-24-2
Formal Name: *trans*-4-bromo-N-2-(dimethylamino)cyclohexyl]-benzamide
Synonym: Bromadoline
MF: C₁₅H₂₁BrN₂O
FW: 325.3
Purity: ≥98%
UV/Vis.: λ_{max}: 239 nm
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-47931E (Item No. 20530) is an analytical reference standard categorized as a utopioid.¹ U-47931E has antinociceptive activity in rodents.² This product is intended for research and forensic applications.

References

- Otte, L., Wilde, M., Auwärter, V., *et al.* Investigation of the μ- and κ-opioid receptor activation by eight new synthetic opioids using the [³⁵S]-GTPγS assay: U-47700, isopropyl U-47700, U-49900, U-47931E, N-methyl U-47931E, U-51754, U-48520, and U-48800. *Drug Test. Anal.* **14**(7), 1187-1199 (2022).
- Hayes, A.G., Sheehan, M.J., and Tyers, M.B. Differential sensitivity of models of antinociception in the rat, mouse, and guinea-pig to μ- and κ-opioid receptor agonists. *Br. J. Pharmacol.* **91**(4), 823-832 (1987).

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, 06/21/2023

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