

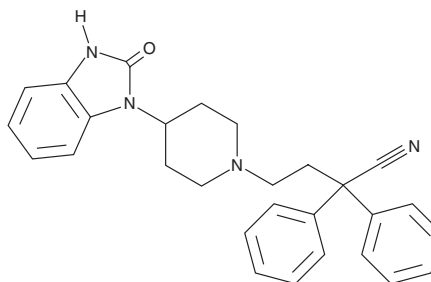
PRODUCT INFORMATION



Despropionyl Bezitramide

Item No. 43337

CAS Registry No.: 83898-28-6
Formal Name: 4-(2,3-dihydro-2-oxo-1H-benzimidazol-1-yl)- α,α -diphenyl-1-piperidinebutanenitrile
Synonym: R-4618
MF: $C_{28}H_{28}N_4O$
FW: 436.6
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Description

Despropionyl bezitramide (Item No. 43337) is an analytical reference standard categorized as an opioid.¹ Despropionyl bezitramide is a metabolite of bezitramide (Item No. 42345). This product is intended for research and forensic applications.

Reference

1. De Baere, S.M., Lambert, W.E., Esmans, E.L., *et al.* Identification and quantitation of despropionyl-bezitramide in postmortem samples by liquid chromatography coupled to electrospray ionization tandem mass spectrometry. *Anal. Chem.* **71**(14), 2908-2914 (1999).

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, 09/10/2025

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