

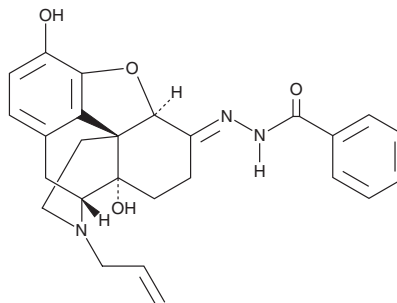
PRODUCT INFORMATION



Naloxone benzoylhydrazone

Item No. 39827

CAS Registry No.: 119630-94-3
Formal Name: 2-[(5 α)-4,5-epoxy-3,14-dihydroxy-17-(2-propenyl)ylidene]hydrazide, benzoic acid
MF: C₂₆H₂₇N₃O₄
FW: 445.5
Purity: $\geq$ 98%
Supplied as: A solid
Storage: -20°C
Stability: $\geq$ 4 years



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

Laboratory Procedures

Naloxone benzoylhydrazone is supplied as a solid. A stock solution may be made by dissolving the naloxone benzoylhydrazone in the solvent of choice, which should be purged with an inert gas. Naloxone benzoylhydrazone is soluble in DMSO.

Description

Naloxone benzoylhydrazone is an antagonist of μ -, κ_1 -, and δ -opioid receptors (pA_{2s} = 8.82, 7.45, and 7.76, respectively, for the mouse receptors).¹ It is also an agonist of the κ_3 -opioid receptor (IC_{50} = 0.7 nM for the guinea pig receptor). Naloxone benzoylhydrazone (1 μ g/animal, i.c.v.) inhibits increases in the latency to tail withdrawal induced by the μ -opioid receptor agonist DAMGO (Item No. 21553) in a hot plate assay in mice.² It increases food intake in rats when intracerebrally administered at doses of 0.1, 1, and 20 μ g/animal.³

References

1. Berzetei-Gurske, I.P., White, A., Polgar, W., *et al.* The in vitro pharmacological characterization of naloxone benzoylhydrazone. *Eur. J. Pharmacol.* **277**(2-3), 257-263 (1995)
2. Paul, D., Levison, J.A., Howard, D.H., *et al.* Naloxone benzoylhydrazone (NalBzoH) analgesia. *J. Pharmacol. Exp. Ther.* **255**(2), 769-774 (1990).
3. Koch, J.E., Pasternak, G.W., Arjune, D., *et al.* Naloxone benzoylhydrazone, a κ_3 opioid agonist, stimulates food intake in rats. *Brain. Res.* **581**(2), 311-314 (1992).

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, 10/09/2023

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