

PRODUCT INFORMATION

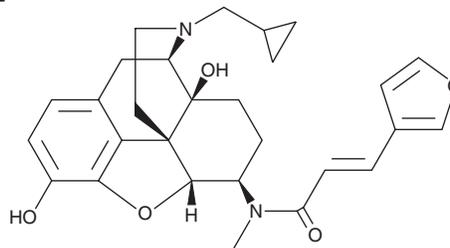


Nalfurafine

Item No. 23186

CAS Registry No.: 152657-84-6
Formal Name: (2E)-N-[(5 α ,6 β)-17-(cyclopropylmethyl)-4,5-epoxy-3,14-dihydroxymorphinan-6-yl]-3-(3-furanyl)-N-methyl-2-propenamamide

Synonym: TRK-820
MF: C₂₈H₃₂N₂O₅
FW: 476.6
Purity: \geq 95%
UV/Vis.: λ_{max} : 210, 278 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



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

Description

Nalfurafine (Item No. 23186) is an analytical reference standard categorized as an opioid.¹ Nalfurafine decreases ethanol consumption in male mice and increases ethanol consumption in female mice.² It is regulated as a Schedule II compound in the United States. This product is intended for research and forensic applications.

References

1. Watanabe, Y., Kitazawa, S., Fujii, H., *et al.* Design, synthesis, and structure-activity relationship of novel opioid κ receptor selective agonists: α -Iminoamide derivatives with an azabicyclo[2.2.2]octene skeleton. *Bioorg. Med. Chem. Lett* **24(21)**, 4980-4983 (2014).
2. French, A.R., Gutridge, A.M., Yuan, J., *et al.* Sex- and β -arrestin-dependent effects of kappa opioid receptor-mediated ethanol consumption. *Pharmacol. Biochem. Behav.* **216**, 173377

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, 07/15/2022

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