

PRODUCT INFORMATION

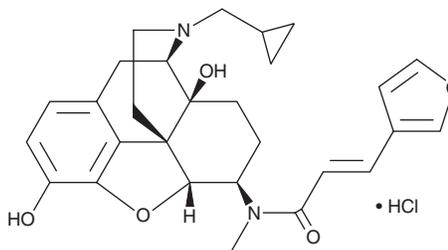


Nalfurafine (hydrochloride)

Item No. 36795

CAS Registry No.: 152658-17-8
Formal Name: N-[(5 α ,6 β)-17-(cyclopropylmethyl)-4,5-epoxy-3,14-dihydroxymorphinan-6-yl]-3-(3-furanyl)-N-methyl-2-propenamide, monohydrochloride

Synonym: TRK-820
MF: C₂₈H₃₂N₂O₅ • HCl
FW: 513.0
Purity: ≥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

Nalfurafine (hydrochloride) (Item No. 36795) 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 (2022).

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, 01/13/2026

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