

PRODUCT INFORMATION

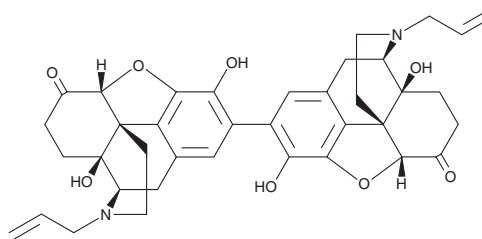


2,2'-Bisnaloxone

Item No. 36678

CAS Registry No.: 211738-08-8
Formal Name: (5 α)-(5' α)-4,5:4',5'-diepoxy-3,3',14,14'-tetrahydroxy-17,17'-bis(2-propen-1-yl)-[2,2'-bimorphinan]-6,6'-dione

MF: C₃₈H₄₀N₂O₈
FW: 625.7
Purity: \geq 95%
UV/Vis.: λ_{max} : 229, 275 nm
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

2,2'-Bisnaloxone is supplied as a solid. A stock solution may be made by dissolving the 2,2'-bisnaloxone in the solvent of choice, which should be purged with an inert gas. 2,2'-Bisnaloxone is slightly soluble in chloroform, DMSO and methanol.

Description

2,2'-Bisnaloxone is a potential impurity in commercial preparations of the opioid antagonist naloxone.¹

Reference

1. Hsu, H.-J., Yang, Y., Pavuluri, V., *et al.* Effect of formulation variables on the nasal permeability and stability of naloxone intranasal formulations. *AAPS PharmSciTech* **20(6)**, 232 (2019).

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, 06/20/2023

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