

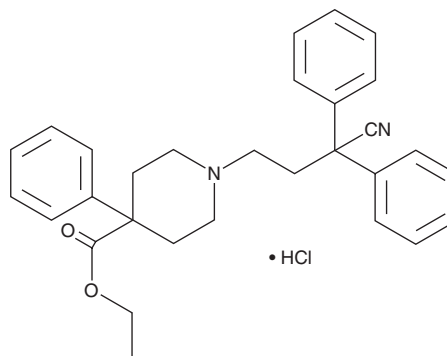
PRODUCT INFORMATION



Diphenoxylate (hydrochloride)

Item No. 20374

CAS Registry No.: 3810-80-8
Formal Name: 1-(3-cyano-3,3-diphenylpropyl)-4-phenyl-4-piperidinecarboxylic acid, ethyl ester, monohydrochloride
Synonym: R 1132
MF: $C_{30}H_{32}N_2O_2 \cdot HCl$
FW: 489.1
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Description

Diphenoxylate (hydrochloride) (Item No. 20374) is an analytical reference standard categorized as an opioid.¹ Diphenoxylate has been detected in products marketed as traditional herbal medicine and has been associated with abuse.^{2,3} Formulations containing diphenoxylate have been used in the treatment of diarrhea. Diphenoxylate is regulated as a Schedule II compound in the United States. This product is intended for research and forensic applications.

References

1. Bonifazi, A., Battiti, F.O., Sanchez, J., *et al.* Novel dual-target μ -opioid receptor and dopamine D_3 receptor ligands as potential nonaddictive pharmacotherapeutics for pain management. *J. Med. Chem.* **64**(11), 7778-7808 (2021).
2. Foroughi, M.H., Akhgari, M., Jokar, F., *et al.* Identification of undeclared active pharmaceutical ingredients in counterfeit herbal medicines used as opioid substitution therapy. *Aus. J. Forensic Sci.* **49**(6), 720-729 (2016).
3. Mehra, A., Sarkar, S., and Basu, D. Lomotil (diphenoxylate) dependence in India. *Indian J. Psychol. Med.* **35**(3), 248-250 (2013).

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, 03/24/2025

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