

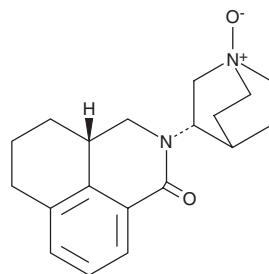
PRODUCT INFORMATION



Palonosetron N-oxide

Item No. 34212

CAS Registry No.: 813425-83-1
Formal Name: (3aS)-2,3,3a,4,5,6-hexahydro-2-[(3S)-1-oxido-1-azabicyclo[2.2.2]oct-3-yl]-1H-benz[de]isoquinolin-1-one
MF: C₁₉H₂₄N₂O₂
FW: 312.4
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Palonosetron N-oxide is supplied as a crystalline solid. A stock solution may be made by dissolving the palonosetron N-oxide in the solvent of choice, which should be purged with an inert gas. Palonosetron N-oxide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of palonosetron N-oxide in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Palonosetron N-oxide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, palonosetron N-oxide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Palonosetron N-oxide has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Palonosetron N-oxide is a metabolite of the serotonin (5-HT) receptor subtype 5-HT₃ antagonist palonosetron (Item No. 16752).¹ It is also a potential impurity in palonosetron preparations.² Palonosetron N-oxide is a degradation product formed by exposure to oxidative stress.

References

1. Stoltz, R.A., Cyong, J.-C., Shah, A., *et al.* Pharmacokinetic and safety evaluation of palonosetron, a 5-hydroxytryptamine-3 receptor antagonist, in U.S. and Japanese healthy subjects. *J. Clin. Pharmacol.* **44(5)**, 520-531 (2004).
2. Vishnu Murthy, M., Srinivas, K., Kumar, R., *et al.* Development and validation of a stability-indicating LC method for determining palonosetron hydrochloride, its related compounds and degradation products using naphthaethyl stationary phase. *J. Pharm. Biomed. Anal.* **56(2)**, 429-435 (2011).

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, 11/20/2024

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