

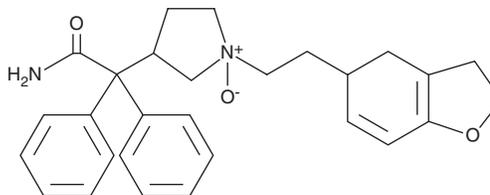
PRODUCT INFORMATION



Darifenacin N-oxide

Item No. 35083

CAS Registry No.: 2416229-88-2
Formal Name: 1-[2-(2,3-dihydro-5-benzofuranyl)ethyl]- α,α -diphenyl-3-pyrrolidineacetamide, 1-oxide
MF: C₂₈H₃₀N₂O₃
FW: 442.5
Purity: $\geq 90\%$
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.

Laboratory Procedures

Darifenacin N-oxide is supplied as a solid. A stock solution may be made by dissolving the darifenacin N-oxide in the solvent of choice, which should be purged with an inert gas. Darifenacin N-oxide is soluble in DMSO.

Description

Darifenacin N-oxide is a potential impurity found in commercial preparations of the muscarinic acetylcholine receptor (mAChR) antagonist darifenacin.¹ It is a degradation product formed under oxidative conditions.

Reference

1. Thomas, S., Paul, S.K., Shandilya, S., *et al.* Identification and structural elucidation of two process impurities and stress degradants in darifenacin hydrobromide active pharmaceutical ingredient by LC-ESI/MSⁿ. *Analyst* **137**(15), 3571-3582 (2012).

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, 12/09/2022

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