

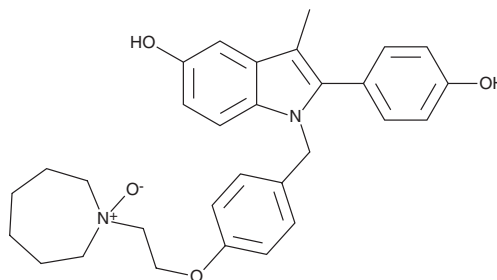
PRODUCT INFORMATION



Bazedoxifene N-oxide

Item No. 35072

CAS Registry No.: 1174289-22-5
Formal Name: 1-[[4-[2-(hexahydro-1-oxido-1H-azepin-1-yl)ethoxy]phenyl]methyl]-2-(4-hydroxyphenyl)-3-methyl-1H-indol-5-ol
MF: C₃₀H₃₄N₂O₄
FW: 486.6
Purity: ≥90%
UV/Vis.: λ_{max}: 225, 297 nm
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

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

Description

Bazedoxifene N-oxide is an oxidative degradation product of the selective estrogen receptor modulator (SERM) bazedoxifene (Item No. 15005).¹

Reference

1. Chandrasekaran, A., McKeand, W.E., Sullivan, P., *et al.* Metabolic disposition of [¹⁴C]bazedoxifene in healthy postmenopausal women. *Drug Metab. Dispos.* **37**(6), 1219-1225 (2009).

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, 11/18/2024

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