

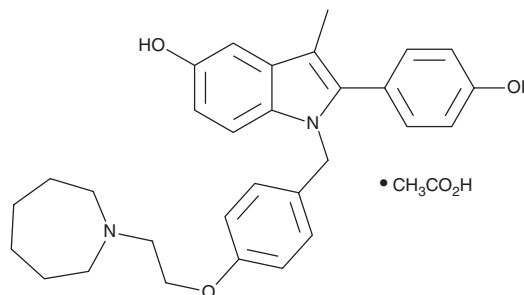
PRODUCT INFORMATION



Bazedoxifene (acetate)

Item No. 15005

CAS Registry No.: 198481-33-3
Formal Name: 1-[[4-[2-(hexahydro-1H-azepin-1-yl)ethoxy]phenyl]methyl]-2-(4-hydroxyphenyl)-3-methyl-1H-indol-5-ol, acetate
Synonym: TSE 424
MF: C₃₀H₃₄N₂O₃ • C₂H₄O₂
FW: 530.7
Purity: ≥98%
UV/Vis.: λ_{max}: 227, 299 nm
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

Bazedoxifene (acetate) is supplied as a crystalline solid. A stock solution may be made by dissolving the bazedoxifene (acetate) in the solvent of choice, which should be purged with an inert gas. Bazedoxifene (acetate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of bazedoxifene (acetate) in these solvents is approximately 5, 14, and 50 mg/ml, respectively.

Bazedoxifene (acetate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, bazedoxifene (acetate) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Bazedoxifene (acetate) has a solubility of approximately 0.04 mg/ml in a 1:20 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Bazedoxifene (acetate) is a third generation selective estrogen receptor modulator (SERM). It is an indole-based ER ligand that binds to both ERα (IC₅₀ = 26 nM) and ERβ (IC₅₀ = 99 nM).¹ Bazedoxifene (acetate) antagonizes 17β-estradiol-dependent MCF-7 and T47D breast cancer cell proliferation *in vitro* as well as hormone-independent growth of MCF-7:⁵C cells that are resistant to long-term estrogen deprivation (80% reduction with 10 nM).² It has been shown to arrest cell cycling by downregulating cyclin D1 and ERα.²

References

1. Komm, B.S., Kharode, Y.P., Bodine, P.V.N., *et al.* Bazedoxifene acetate: A selective estrogen receptor modulator with improved selectivity. *Endocrinology* **146**(9), 3999-4008 (2005).
2. Lewis-Wambi, J.S., Kim, H., Curpan, R., *et al.* The selective estrogen receptor modulator bazedoxifene inhibits hormone-independent breast cancer cell growth and down-regulates estrogen receptor α and cyclin D1. *Mol. Pharmacol.* **80**(4), 610-620 (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, 10/06/2022

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