

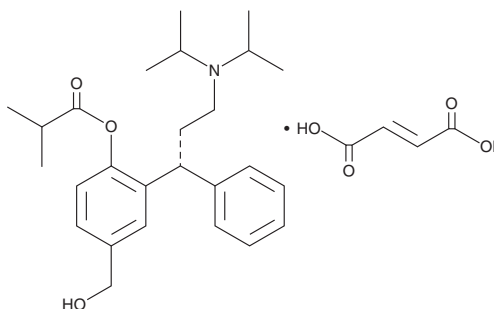
PRODUCT INFORMATION



Fesoterodine (fumarate)

Item No. 23777

CAS Registry No.: 286930-03-8
Formal Name: 2-methyl-propanoic acid, 2-[(1R)-3-[bis(1-methylethyl)amino]-1-phenylpropyl]-4-(hydroxymethyl) phenyl ester, (2E)-2-butenedioate
Synonyms: SMP 8272, SPM 907
MF: C₂₆H₃₇NO₃ • C₄H₄O₄
FW: 527.7
Purity: ≥98%
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

Fesoterodine (fumarate) is supplied as a solid. A stock solution may be made by dissolving the fesoterodine (fumarate) in the solvent of choice, which should be purged with an inert gas. Fesoterodine (fumarate) is slightly soluble in methanol. Fesoterodine (fumarate) is also slightly soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Fesoterodine is an antagonist of muscarinic acetylcholine receptors (K_is = 15.8, 11.2, 12.6, 22.4, and 25.1 nM for human M₁₋₅, respectively).¹ *In vivo*, fesoterodine (0.01 mg/kg, i.v.) reduces micturition pressure and increases bladder capacity and intercontraction interval (ICI) in conscious female rats. Fesoterodine also reduces intermicturition pressure and prevents bladder overactivity in a rat model of spinal cord injury-induced overactive bladder.²

References

1. Ney, P., Pandita, R.K., Newgreen, D.T., *et al.* Pharmacological characterization of a novel investigational antimuscarinic drug, fesoterodine, *in vitro* and *in vivo*. *BJU Int.* **101(8)**, 1036-1042 (2008).
2. Biardeau, X., Przydacz, M., Aharony, S., *et al.* Early fesoterodine fumarate administration prevents neurogenic detrusor overactivity in a spinal cord transected rat model. *PLoS One* **12(1)**, e0169694 (2017).

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