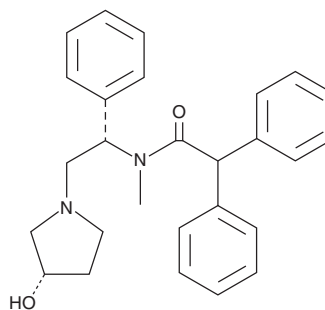


PRODUCT INFORMATION



Asimadoline Item No. 26411

CAS Registry No.: 153205-46-0
Formal Name: N-[(1S)-2-[(3S)-3-hydroxy-1-phenylethyl]-1-pyrrolidinyl]-N-methyl- α -phenyl-benzeneacetamide
Synonym: EMD 61753
MF: C₂₇H₃₀N₂O₂
FW: 414.5
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Asimadoline is supplied as a crystalline solid. A stock solution may be made by dissolving the asimadoline in the solvent of choice, which should be purged with an inert gas. Asimadoline is soluble in DMSO at a concentration of approximately 103 mg/ml.

Description

Asimadoline is a potent κ -opioid receptor (KOR) agonist (IC₅₀s = 5.6 and 1.2 nM for guinea pig and human receptors, respectively).¹ It is 501- and 498-fold selective for κ -opioid over μ - and δ -opioid receptors, respectively. Asimadoline is spasmolytic in isolated rat duodenum (IC₅₀ = 4.2 μ M) and inhibits spontaneous contractions of isolated rat uterus (IC₅₀ = 12.7 μ M). *In vivo*, asimadoline reduces joint damage in a rat model of arthritis induced by complete Freund's adjuvant (CFA). Asimadoline (25 mg/kg) also reduces the abdominal withdrawal reflex in a model of visceral pain induced by colonic distension in wild-type, but not KOR^{-/-}, mice.²

References

1. Camilleri, M. Asimadoline, a κ -opioid agonist, and visceral sensation. *Neurogastroenterol. Motil.* **20(9)**, 971-979 (2008).
2. Larsson, M.H., Bayati, A., Lindström, E., *et al.* Involvement of kappa-opioid receptors in visceral nociception in mice. *Neurogastroenterol. Motil.* **20(10)**, 1157-1164 (2008).

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, 10/12/2022

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