

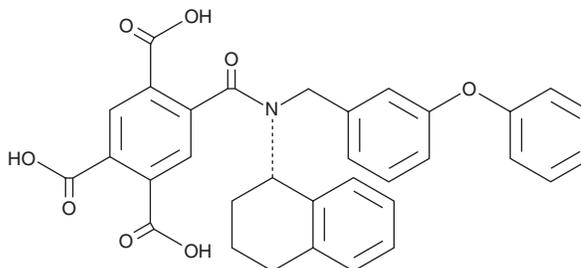
PRODUCT INFORMATION



A-317491

Item No. 19256

CAS Registry No.: 475205-49-3
Formal Name: 5-[[[(3-phenoxyphenyl)methyl] [(1S)-1,2,3,4-tetrahydro-1-naphthalenyl]amino]carbonyl]-1,2,4-benzenetricarboxylic acid
MF: C₃₃H₂₇NO₈
FW: 565.6
Purity: ≥98%
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

A-317491 is supplied as a crystalline solid. A stock solution may be made by dissolving the A-317491 in the solvent of choice. A-317491 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of A-317491 in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of A-317491 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of A-317491 in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

A-317491 is a non-nucleotide antagonist of the purinergic receptors P2X₃ and P2X_{2/3} (K_is = 22-92 nM in a calcium influx assay) that is selective over other purinergic and neurotransmitter receptors, ion channels, and enzymes (IC₅₀s = >10 μM).¹ It reduces thermal hyperalgesia induced by complete Freund's adjuvant in rats (ED₅₀ = 30 μmol/kg, s.c.) and nociceptive responses to chemically-induced pain. Chronic administration of A-317491 (30 μmol/kg, twice daily) reduces early, but not late, phase pain in a mouse model of bone cancer.² A-317491 does not readily cross the blood-brain barrier and reduces pain responses when applied topically in a rat model of inflammatory pain, indicating that it acts primarily at peripheral purinergic receptors.³

References

1. Jarvis, M.F., Burgard, E.C., McGaraughty, S., *et al.* A-317491, a novel potent and selective non-nucleotide antagonist of P2X₃ and P2X_{2/3} receptors, reduces chronic inflammatory and neuropathic pain in the rat. *Proc. Natl. Acad. Sci. U.S.A.* **99**(26), 17179-17184 (2002).
2. Hansen, R.R., Nasser, A., Falk, S., *et al.* Chronic administration of the selective P2X₃, P2X_{2/3} receptor antagonist, A-317491, transiently attenuates cancer-induced bone pain in mice. *Eur. J. Pharmacol.* **688**(1-3), 27-34 (2012).
3. Wu, G., Whiteside, G.T., Lee, G., *et al.* A-317491, a selective P2X₃/P2X_{2/3} receptor antagonist, reverses inflammatory mechanical hyperalgesia through action at peripheral receptors in rats. *Eur. J. Pharmacol.* **504**(1-2), 45-53 (2004).

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, 10/21/2022

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