

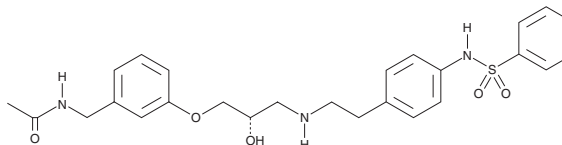
PRODUCT INFORMATION



L-748,337

Item No. 19643

CAS Registry No.: 244192-94-7
Formal Name: N-[[3-[(2S)-2-hydroxy-3-[[2-[4-[(phenylsulfonyl)amino]phenyl]ethyl]amino]propoxy]phenyl]methyl]-acetamide
MF: C₂₆H₃₁N₃O₅S
FW: 497.6
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

L-748,337 is supplied as a crystalline solid. A stock solution may be made by dissolving the L-748,337 in the solvent of choice, which should be purged with an inert gas. L-748,337 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of L-748,337 in ethanol and DMF is approximately 30 mg/ml and approximately 25 mg/ml in DMSO.

L-748,337 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, L-748,337 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. L-748,337 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

L-748,337 is a β_3 -adrenergic receptor antagonist that displays selectivity over β_1 and β_2 receptors (K_i s = 4, 390, and 204 nM, respectively).¹ It has been shown to couple with G_i to activate MAPK signaling.² L-748,337 can inhibit agonist-stimulated lipolysis in rhesus monkey adipocytes, as well as agonist-induced relaxation of smooth muscle.¹

References

1. Candelmore, M.R., Deng, L., Tota, L., *et al.* Potent and selective human β_3 -adrenergic receptor antagonists. *J. Pharmacol. Exp. Ther.* **290**(2), 649-655 (1999).
2. Sato, M., Hutchinson, D.S., Evans, B.A., *et al.* The β_3 -adrenoceptor agonist 4-[[[(Hexylamino)carbonyl]amino]-N-[4-[2-[[[(2S)-2-hydroxy-3-(4-hydroxyphenoxy)propyl]amino]ethyl]-phenyl]-benzenesulfonamide (L755507) and antagonist (S)-N-[4-[2-[[3-[3-(acetamidomethyl)phenoxy]-2-hydroxypropyl]amino]-ethyl]phenyl]benzenesulfonamide (L748337) activate different signaling pathways in Chinese hamster ovary-K1 cells stably expressing the human β_3 -adrenoceptor. *Mol. Pharmacol.* **74**(5), 1417-1428 (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, 12/16/2022

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