

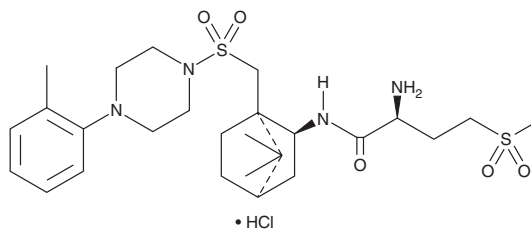
PRODUCT INFORMATION



L-368,899 (hydrochloride)

Item No. 29868

CAS Registry No.: 160312-62-9
Formal Name: [1S-[1 α ,2 α (R*),4 β]]-2-amino-N-[7,7-dimethyl-1-[[[4-(2-methylphenyl)-1-piperazinyl]sulfonyl]methyl]bicyclo[2.2.1]hept-2-yl]-4-(methylsulfonyl)butanamide, monohydrochloride
MF: C₂₆H₄₂N₄O₅S₂ • HCl
FW: 591.2
Purity: ≥98%
UV/Vis.: λ_{\max} : 241 nm
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

L-368,899 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the L-368,899 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. L-368,899 (hydrochloride) is soluble in an organic solvent such as DMSO. It is also soluble in water. The solubility of L-368,899 (hydrochloride) in DMSO and water is approximately 100 mM. We do not recommend storing the aqueous solution for more than one day.

Description

L-368,899 is a non-peptide oxytocin receptor antagonist (IC_{50} s = 8.9 and 26 nM in radioligand binding assays using isolated human and rat uterus, respectively).¹ It is selective for oxytocin receptors over vasopressin V_{1a} and V₂ receptors (IC_{50} s = 370 and 570 nM, respectively). L-368,899 reduces oxytocin-induced contraction of isolated rat uterus (pA_2 = 8.9) and *in situ* rat uterus (ED_{50} = 0.35 mg/kg, i.v.).

Reference

- Williams, P.D., Anderson, P.S., Ball, R.G., *et al.* 1-((7,7-Dimethyl-2(S)-(2(S)-amino-4-(methylsulfonyl)butyramido)bicyclo [2.2.1]-heptan-1(S)-yl)methyl)sulfonyl)-4-(2-methylphenyl)piperazine (L-368,899): An orally bioavailable, non-peptide oxytocin antagonist with potential utility for managing preterm labor. *J. Med. Chem.* **37**(5), 565-571 (1994).

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, 12/12/2022

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