

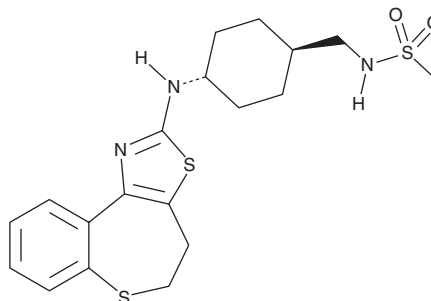
PRODUCT INFORMATION



Lu AA33810

Item No. 32913

CAS Registry No.: 304008-29-5
Formal Name: N-[[*trans*-4-[(4,5-dihydro[1]benzothiepine[5,4-d]thiazol-2-yl)amino]cyclohexyl]methyl]-methanesulfonamide
MF: C₁₉H₂₅N₃O₂S₃
FW: 423.6
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 244, 255, 363 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

Lu AA33810 is supplied as a solid. A stock solution may be made by dissolving the Lu AA33810 in the solvent of choice, which should be purged with an inert gas. Lu AA33810 is soluble in the organic solvent DMSO at a concentration of approximately 100 mM.

Description

Lu AA33810 is an antagonist of neuropeptide Y (NPY) receptor Y₅ (K_i = 1.5 nM for the human receptor).¹ It is selective for Y₅ over Y₁, Y₂, and Y₄ (K_is = >10 μM for all in HEK293 membranes expressing rat receptors). Lu AA33810 also binds the serotonin (5-HT) receptor subtypes 5-HT_{1A} and 5-HT_{2B} (K_is = 478 and 247 nM, respectively). It inhibits NPY-induced calcium mobilization and cAMP accumulation in COS-7 and HEK293 cells expressing the rat Y₅ receptor, respectively, in a concentration-dependent manner. Lu AA33810 (10 mg/kg) increases social interaction time in a social exploration test, as well as decreases immobility time in the forced swim test, in rats, indicating anxiolytic- and antidepressant-like activities, respectively.

Reference

1. Walker, M.W., Wolinsky, T.D., Jubian, V., *et al.* The novel neuropeptide Y Y₅ receptor antagonist Lu AA33810 [N-[[*trans*-4-[(4,5-dihydro[1]benzothiepine[5,4-d]thiazol-2-yl)amino]cyclohexyl]methyl]-methanesulfonamide] exerts anxiolytic- and antidepressant-like effects in rat models of stress sensitivity. *J. Pharmacol. Exp. Ther.* **328**(3), 900-911 (2009).

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, 11/10/2022

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