

PRODUCT INFORMATION



JYL1421

Item No. 20458

CAS Registry No.: 401907-26-4

Formal Name: N-[4-[[[4-(1,1-dimethylethyl)phenyl]methyl]amino]thioxomethyl]amino]methyl]-2-fluorophenyl]-methanesulfonamide

MF: C₂₀H₂₆FN₃O₂S₂

FW: 423.6

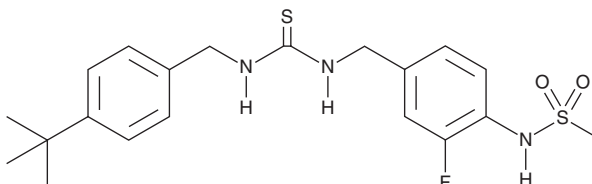
Purity: ≥95%

UV/Vis.: λ_{max}: 225 nm

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

JYL1421 is supplied as a crystalline solid. A stock solution may be made by dissolving the JYL1421 in the solvent of choice. JYL1421 is soluble in the organic solvent DMSO, which should be purged with an inert gas.

Description

JYL1421 is an antagonist of transient receptor potential vanilloid 1 (TRPV1).¹ It inhibits calcium uptake induced by capsaicin (Item Nos. 92350 | 10010743) in CHO cells expressing rat TRPV1 (EC₅₀ = 9.2 nM). JYL1421 inhibits capsaicin-induced release of the neuropeptides somatostatin, substance P (Item No. 24035), and calcitonin gene-related peptide (CGRP) from isolated rat trachea (IC₅₀s = 227-491 nM).² It inhibits capsaicin-induced hypothermia and hypotension in rats when administered at doses of 2 and 0.4 mg/kg, respectively. JYL1421 (2 mg/kg) also reduces the number of wiping movements induced by ocular administration of capsaicin in rats. Unlike several other TRPV1 antagonists, JYL1421 does not induce hyperthermia in rats when administered at doses ranging from 1.02 to 32.77 μmol/kg.³

References

1. Wang, Y., Szabo, T., Welter, J.D., *et al.* High affinity antagonists of the vanilloid receptor. *Mol. Pharmacol.* **62**(4), 947-956 (2002).
2. Jakab, B., Helyes, Z., Varga, A., *et al.* Pharmacological characterization of the TRPV1 receptor antagonist JYL1421 (SC0030) in vitro and in vivo in the rat. *Eur. J. Pharmacol.* **517**(1-2), 35-44 (2005).
3. Garami, A., Shimansky, Y.P., Pakai, E., *et al.* Contributions of different modes of TRPV1 activation to TRPV1 antagonist-induced hyperthermia. *J. Neurosci.* **30**(4), 1435-1440 (2010).

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, 03/27/2025

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