

PRODUCT INFORMATION

A1480LS

Item No. 45340

Formal Name: (S)-(2-benzyl-4-(6-(trifluoromethyl)pyrazin-2-yl)piperazin-1-yl)(5-phenyl-2H-indazol-2-yl)methanone

MF: C₃₀H₂₅F₃N₆O

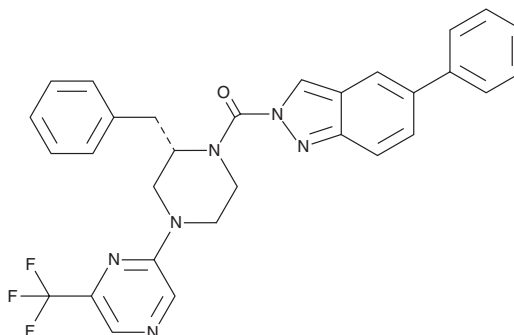
FW: 542.6

Purity: ≥98%

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

A1480LS is supplied as a solid. A stock solution may be made by dissolving the A1480LS in the solvent of choice, which should be purged with an inert gas. A1480LS is soluble (≥10 mg/ml) in DMSO.

Description

A1480LS is a peripherally biased inhibitor of diacylglycerol lipase α (DAGLα) and DAGLβ (IC₅₀s = 6 and 4 nM, respectively, in human brain membrane homogenates).¹ It reduces levels of the DAGL-derived endocannabinoid 2-arachidonoyl glycerol (2-AG) in peripheral tissue but not in the brain in mice when administered at 20 mg/kg. A1480LS (20 mg/kg, p.o.) reduces mechanical allodynia in rat models of inflammatory pain and chemotherapy-induced neuropathy without inducing CNS adverse effects in the Irwin primary observation test.

Reference

1. Tufail, Y.Z., Guijas, C., Kummer, D.A., et al. Suppression of pain transmission and behavior by inhibition of peripheral diacylglycerol metabolism. *Cell Chem. Bio.* **33**(1), 74-90.e19 (2025).

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, 06/09/2026

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