

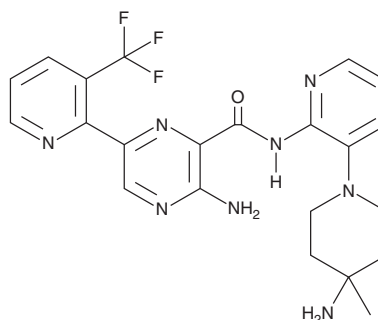
PRODUCT INFORMATION



LXS-196

Item No. 31455

CAS Registry No.: 1874276-76-2
Formal Name: 3-amino-N-[3-(4-amino-4-methyl-1-piperidiny)-2-pyridinyl]-6-[3-(trifluoromethyl)-2-pyridinyl]-2-pyrazinecarboxamide
MF: C₂₂H₂₃F₃N₈O
FW: 472.5
Purity: ≥95%
UV/Vis.: λ_{max}: 288, 375 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

LXS-196 is supplied as a crystalline solid. A stock solution may be made by dissolving the LXS-196 in the solvent of choice, which should be purged with an inert gas. LXS-196 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of LXS-196 in these solvents is approximately 1 mg/ml.

LXS-196 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, LXS-196 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. LXS-196 has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

LXS-196 is a PKC inhibitor (IC₅₀s = 1.9 and 0.4 nM for PKCα and PKCθ, respectively).¹ It is selective for PKCα and PKCθ over GSK3β (IC₅₀ = 3,100 nM). It inhibits proliferation of TMD8 B cell lymphoma and 92.1 uveal melanoma cells (IC₅₀s = 900 and 184 nM, respectively) but not SK-MEL-28 skin melanoma cells (IC₅₀ = >10,000 nM).

Reference

1. Luzzio, M.J., Papillon, J., and Visser, M.S. Protein kinase C inhibitors and methods of their use. *Novartis AG. WO 2016/020864 A1* (2016).

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, 10/25/2022

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