

PRODUCT INFORMATION



XL147 analog 2

Item No. 16299

CAS Registry No.: 1349796-36-6
Formal Name: N-[4-[[[3-[(3,5-dimethoxyphenyl)amino]-2-quinoxaliny]amino]sulfonyl]phenyl]-3-methoxy-4-methyl-benzamide

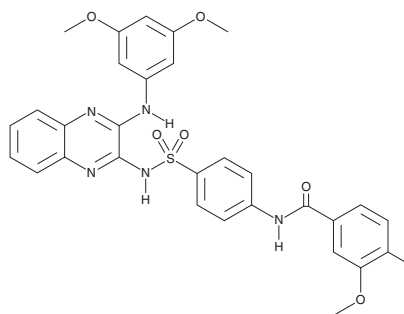
MF: C₃₁H₂₉N₅O₆S
FW: 599.7

Purity: ≥95%

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

XL147 analog 2 is supplied as a crystalline solid. A stock solution may be made by dissolving the XL147 analog 2 in the solvent of choice, which should be purged with an inert gas. XL147 analog 2 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of XL147 analog 2 in these solvents is approximately 10 mg/ml.

XL147 analog 2 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, XL147 analog 2 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. XL147 analog 2 has a solubility of approximately 0.1 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

XL147 analog 2 is an analog of XL147 (Item No. 18011).

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