

PRODUCT INFORMATION



KDS-0071

Item No. 41996

CAS Registry No.: 1973485-18-5
Formal Name: N-[5-[4-[(1,1-dioxido-4-thiomorpholinyl)methyl]-3-fluorophenyl]imidazo[1,2-a]pyridin-2-yl]-cyclopropanecarboxamide

MF: C₂₂H₂₃FN₄O₃S

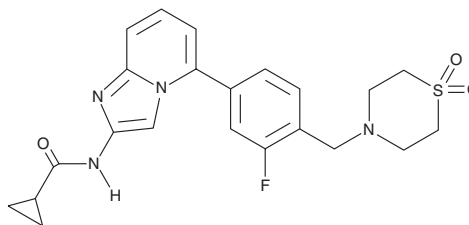
FW: 442.5

Purity: ≥95%

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

KDS-0071 is supplied as a solid. A stock solution may be made by dissolving the KDS-0071 in the solvent of choice, which should be purged with an inert gas. KDS-0071 is soluble (≥10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

KDS-0071 is an imidazole-containing compound.¹

Reference

1. Jin, L., Shung, W., Minmin, Y., *et al.* Substituted imidazo[1,2-a]pyridin-2-ylamine compounds, and pharmaceutical compositions and methods of use thereof. *JN Therapeutics*, **WO2016/119700AI** (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, 10/25/2024

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