

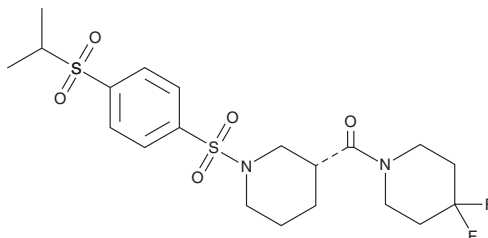
PRODUCT INFORMATION



DX3-213B

Item No. 36783

CAS Registry No.: 2749555-66-4
Formal Name: (R)-(4,4-difluoropiperidin-1-yl) (1-((4-(isopropylsulfonyl)phenyl)sulfonyl)piperidin-3-yl)methanone
MF: C₂₀H₂₈F₂N₂O₅S₂
FW: 478.6
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

DX3-213B is supplied as a crystalline solid. A stock solution may be made by dissolving the DX3-213B in the solvent of choice, which should be purged with an inert gas. DX3-213B is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of DX3-213B in DMSO is approximately 3 mg/ml and approximately 2 mg/ml in ethanol and DMF.

Description

DX3-213B is an inhibitor of mitochondrial complex I, also known as NADH dehydrogenase (IC₅₀ = 3.6 nM).¹ It inhibits the proliferation of MiaPaCa-2 cells cultured in galactose- but not glucose-containing media (IC₅₀s = 9.1 and >3,000 nM, respectively). DX3-213B (10 mg/kg) reduces tumor volume in a Panc02 murine pancreatic ductal adenocarcinoma model.

Reference

1. Xue, D., Xu, Y., Kyani, A., *et al.* Multiparameter optimization of oxidative phosphorylation inhibitors for the treatment of pancreatic cancer. *J. Med. Chem.* **65**(4), 3404-3419 (2022).

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, 11/13/2025

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