

PRODUCT INFORMATION

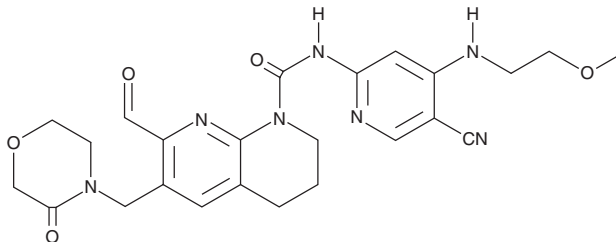


FGFR4-IN-1

Item No. 21342

CAS Registry No.: 1708971-72-5
Formal Name: N-[5-cyano-4-[(2-methoxyethyl)amino]-2-pyridinyl]-7-formyl-3,4-dihydro-6-[[3-oxo-4-morpholinyl)methyl]-1,8-naphthyridine-1(2H)-carboxamide

MF: C₂₄H₂₇N₇O₅
FW: 493.5
Purity: ≥98%
UV/Vis.: λ_{max}: 255, 274 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

FGFR4-IN-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the FGFR4-IN-1 in the solvent of choice, which should be purged with an inert gas. FGFR4-IN-1 is soluble in organic solvents such as DMSO (sonicated) and chloroform. The solubility of FGFR4-IN-1 in these solvents is approximately 1 and 30 mg/ml, respectively.

Description

FGFR-IN-1 is an inhibitor of FGF receptor 4 (FGFR4; IC₅₀ = 1.3 nM).¹ It is selective for FGFR4 over FGFR1 and a panel of 36 kinases (IC₅₀s = >10 μM for all). It inhibits proliferation of Hep3B and Huh7 hepatic cancer cells with IC₅₀ values of 1.1 and 2.5 nM, respectively. FGFR-IN-1 reduces tumor growth by 62.7 and 70.8% in a Huh7 mouse xenograft model when administered at doses of 30 and 100 mg/kg per day, respectively.

Reference

1. Buschmann, N., Fairhurst, R.A., Furet, P., *et al.* Ring-fused bicyclic pyridyl derivatives as FGFR4 inhibitors. *Novartis*. **WO2015059668A1** (2015).

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, 10/19/2022

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