

PRODUCT INFORMATION

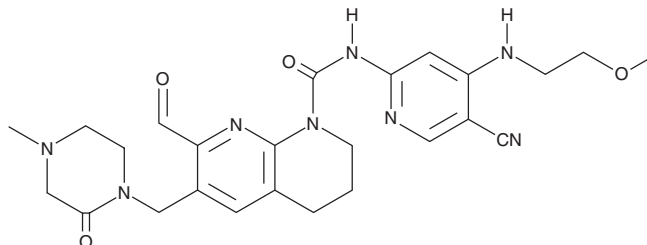


FGF401

Item No. 23029

CAS Registry No.: 1708971-55-4
Formal Name: N-[5-cyano-4-[(2-methoxyethyl)amino]-2-pyridinyl]-7-formyl-3,4-dihydro-6-[(4-methyl-2-oxo-1-piperazinyl)methyl]-1,8-naphthyridine-1(2H)-carboxamide

Synonym: Roblitinib
MF: C₂₅H₃₀N₈O₄
FW: 506.6
Purity: ≥95%
UV/Vis.: λ_{max}: 227, 248, 284 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

FGF401 is supplied as a crystalline solid. A stock solution may be made by dissolving the FGF401 in the solvent of choice. FGF401 is soluble in organic solvents such as DMSO (gently warm and vortex) and chloroform which should be purged with an inert gas. The solubility of FGF401 in these solvents is approximately 5 and 30 mg/ml, respectively.

Description

FGF401 is an inhibitor of FGF receptor 4 (FGFR4; IC₅₀ = 1.9 nM).¹ It inhibits proliferation of Huh7 carcinoma cells with an IC₅₀ value of 12 nM. FGF401 induces bile acid synthesis in dogs, increasing plasma and fecal levels of bile acids when administered at a dose of 45 mg/kg per day for four weeks.² It also decreases serum cholesterol when administered at doses ranging from 5 to 45 mg/kg per day.

References

1. Buschmann, N., Fairhurst, R.A., Furet, P., *et al.* Ring-fused bicyclic pyridyl derivatives as fgfr4 inhibitors. *Novartis*. **WO2015059668A1** (2015).
2. Schadt, H.S., Wolf, A., Mahl, J.A., *et al.* Bile acid sequestration by cholestyramine mitigates FGFR4 inhibition-induced ALT elevation. *Toxicol. Sci.* **163(1)**, 265-278 (2018).

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, 12/02/2022

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