

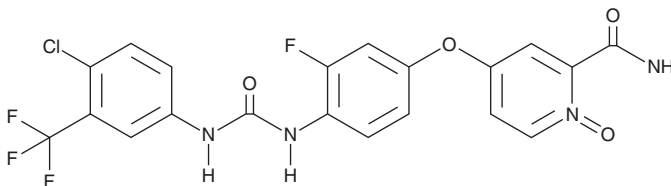
PRODUCT INFORMATION



N-desmethyl Regorafenib N-oxide

Item No. 35102

CAS Registry No.: 835621-12-0
Formal Name: 4-[4-[[[4-chloro-3-(trifluoromethyl)phenyl]amino]carbonyl]amino]-3-fluorophenoxy]-2-pyridinecarboxamide 1-oxide
MF: C₂₀H₁₃ClF₄N₄O₄
FW: 484.8
Purity: ≥90%
UV/Vis.: λ_{max}: 264 nm
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

N-desmethyl Regorafenib N-oxide is supplied as a solid. A stock solution may be made by dissolving the N-desmethyl Regorafenib N-oxide in the solvent of choice, which should be purged with an inert gas. N-desmethyl Regorafenib N-oxide is soluble in DMSO.

Description

N-desmethyl Regorafenib N-oxide is an active metabolite of the multi-kinase inhibitor regorafenib (Item No. 18498).² It is formed from regorafenib by the cytochrome P450 (CYP) isoform CYP3A4. N-desmethyl Regorafenib N-oxide inhibits VEGFR2, Tie2, c-Kit, and B-RAF *in vitro*, as well as inhibits tumor growth in HT-29 and MDA-MB-231 mouse xenograft models when administered at a dose of 1 mg/kg.¹

References

1. Zopf, D., Heinig, R., Thierauch, K.-H., *et al.* Abstract 1666: Regorafenib (BAY 73-4506): Preclinical pharmacology and clinical identification and quantification of its major metabolites. *Cancer Res.* **70(8 Suppl.)**, (2010).
2. Li, Y.-H., Lin, Q.-M., Pang, N.-H., *et al.* Functional characterization of 27 CYP3A4 protein variants to metabolize regorafenib *in vitro*. *Basic Clin. Pharmacol. Toxicol.* **125(4)**, 337-344 (2019).

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, 12/09/2022

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