

# PRODUCT INFORMATION

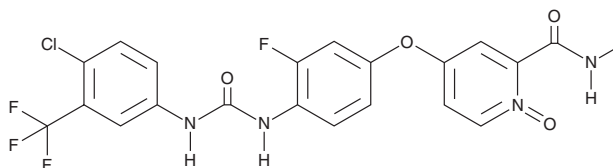


## Regorafenib N-oxide

Item No. 35101

CAS Registry No.: 835621-11-9  
Formal Name: 4-[4-[[[4-chloro-3-(trifluoromethyl)phenyl]amino]carbonyl]amino]-3-fluorophenoxy]-N-methyl-2-pyridinecarboxamide 1-oxide

MF: C<sub>21</sub>H<sub>15</sub>ClF<sub>4</sub>N<sub>4</sub>O<sub>4</sub>  
FW: 498.8  
Purity: ≥95%  
UV/Vis.: λ<sub>max</sub>: 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

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

### Description

Regorafenib N-oxide is an active metabolite of the multi-kinase inhibitor regorafenib (Item No. 18498).<sup>2</sup> It is formed from regorafenib by the cytochrome P450 (CYP) isoform CYP3A4. 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.<sup>1</sup>

### 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, 11/28/2022

#### 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