

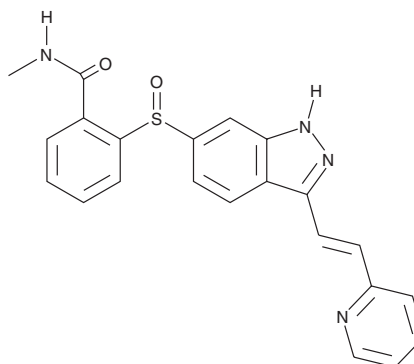
PRODUCT INFORMATION



Axitinib Sulfoxide

Item No. 28042

CAS Registry No.: 1347304-18-0
Formal Name: N-methyl-2-[[3-[(1E)-2-(2-pyridinyl)ethenyl]-1H-indazol-6-yl]sulfinyl]-benzamide
MF: C₂₂H₁₈N₄O₂S
FW: 402.5
Purity: ≥98%
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

Axitinib sulfoxide is supplied as a solid. A stock solution may be made by dissolving the axitinib sulfoxide in the solvent of choice, which should be purged with an inert gas. Axitinib sulfoxide is soluble in the organic solvent DMSO.

Description

Axitinib sulfoxide is a major inactive metabolite of the tyrosine kinase inhibitor axitinib (Item No. 13813).^{1,2} It is formed from axitinib primarily by the cytochrome P450 (CYP) isoform CYP3A4 with minor contributions from CYP3A5 and CYP2C19.¹

References

1. Zientek, M.A., Goosen, T.C., Tseng, E., *et al.* In vitro kinetic characterization of axitinib metabolism. *Drug Metab. Dispos.* **44(1)**, 102-114 (2016).
2. Smith, B.J., Pithavala, Y., Bu, H.Z., *et al.* Pharmacokinetics, metabolism, and excretion of [¹⁴C]axitinib, a vascular endothelial growth factor receptor tyrosine kinase inhibitor, in humans. *Drug Metab. Dispos.* **42(5)**, 918-931 (2014).

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