

PRODUCT INFORMATION

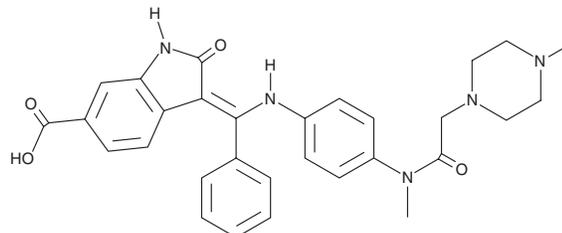


BIBF 1202

Item No. 27176

CAS Registry No.: 894783-71-2
Formal Name: (3Z)-2,3-dihydro-3-[[[4-[methyl[2-(4-methyl-1-piperazinyl)acetyl]amino]phenyl]amino]phenylmethylene]-2-oxo-1H-indole-6-carboxylic acid

MF: C₃₀H₃₁N₅O₄
FW: 525.6
Purity: ≥95%
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

BIBF 1202 is supplied as a solid. A stock solution may be made by dissolving the BIBF 1202 in the solvent of choice, which should be purged with an inert gas. BIBF 1202 is soluble in the organic solvent DMSO.

Description

BIBF 1202 is an active metabolite of the VEGFR, FGFR, and PDGFR inhibitor BIBF 1120 (nintedanib; Item No. 11022).^{1,2} It is formed from BIBF 1120 by intracellular esterases.¹ BIBF 1202 inhibits VEGFR2 with an IC₅₀ value of 62 nM.

References

- Hilberg, F., Roth, G.J., Krssak, M., *et al.* BIBF 1120: Triple angiokinase inhibitor with sustained receptor blockade and good antitumor efficacy. *Cancer Res.* **68**(12), 4774-4782 (2008).
- Lin, D., Qiao, L.M., Zhang, Y.N., *et al.* Simultaneous determination of nintedanib and its metabolite by UPLC-MS/MS in rat plasma and its application to a pharmacokinetic study. *J. Pharm. Biomed. Anal.* **117**, 173-177 (2016).

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, 12/08/2022

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