

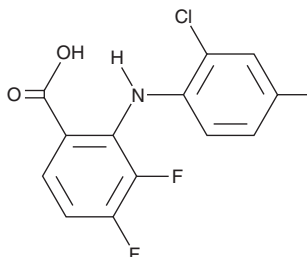
PRODUCT INFORMATION



Zapnometinib

Item No. 38517

CAS Registry No.: 303175-44-2
Formal Name: 2-[(2-chloro-4-iodophenyl)amino]-3,4-difluoro-benzoic acid
Synonyms: ATR-002, PD 0184264
MF: C₁₃H₇ClF₂INO₂
FW: 409.6
Purity: ≥95%
UV/Vis.: λ_{max}: 280, 325 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

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

Description

Zapnometinib is a MEK inhibitor (IC₅₀ = 5.73 nM) and an active metabolite of CI-1040 (Item No. 11580).^{1,2} It reduces viral titers in A549 and MDCK cells infected with influenza A (EC₅₀s = 6.356 and 4.498 μM, respectively).² Zapnometinib (50 mg/kg per day) increases survival in H1N1pdm09 influenza A-infected mice.

References

1. Teclé, H., Shao, J., Li, Y., *et al.* Beyond the MEK-pocket: Can current MEK kinase inhibitors be utilized to synthesize novel type III NCKIs? Does the MEK-pocket exist in kinases other than MEK? *Bioorg. Med. Chem. Lett.* **19(1)**, 226-229 (2009).
2. Laure, M., Hamza, H., Koch-Heier, J., *et al.* Antiviral efficacy against influenza virus and pharmacokinetic analysis of a novel MEK-inhibitor, ATR-002, in cell culture and in the mouse model. *Antiviral Res.* **178**, 104806 (2020).

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, 06/08/2023

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