

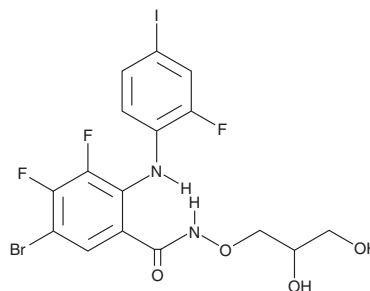
PRODUCT INFORMATION



PD 318088

Item No. 30265

CAS Registry No.: 391210-00-7
Formal Name: 5-bromo-N-(2,3-dihydroxypropoxy)-3,4-difluoro-2-[(2-fluoro-4-iodophenyl)amino]-benzamide
MF: C₁₆H₁₃BrF₃IN₂O₄
FW: 561.1
Purity: ≥98%
UV/Vis.: λ_{max}: 239, 284 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

PD 318088 is supplied as a crystalline solid. A stock solution may be made by dissolving the PD 318088 in the solvent of choice, which should be purged with an inert gas. PD 318088 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PD 318088 in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

PD 318088 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PD 318088 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. PD 318088 has a solubility of approximately 0.20 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

PD 318088 is an allosteric inhibitor of MEK1 and an analog of the MEK inhibitor PD 184352 (CI-1040; Item No. 11580).^{1,2} PD318088 binds to MEK1 concurrently with Mg-ATP and stabilizes an inactive conformation of MEK1, as well as prevents formation of tetramers and higher-order oligomers.

References

- Ohren, J.F., Chen, H., Pavlovsky, A., *et al.* Structures of human MAP kinase kinase 1 (MEK1) and MEK2 describe novel noncompetitive kinase inhibition. *Nat. Struct. Mol. Biol.* **11(12)**, 1192-1197 (2004).
- Shang, J., Lu, S., Jiang, Y., *et al.* Allosteric modulators of MEK1: Drug design and discovery. *Chem. Biol. Drug Des.* **88(4)**, 485-497 (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, 05/14/2020

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