

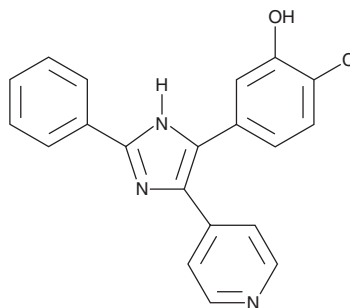
PRODUCT INFORMATION



L-779,450

Item No. 17807

CAS Registry No.: 303727-31-3
Formal Name: 2-chloro-5-[2-phenyl-4-(4-pyridinyl)-1H-imidazol-5-yl]-phenol
Synonym: Raf Kinase Inhibitor IV
MF: C₂₀H₁₄ClN₃O
FW: 347.8
Purity: ≥98%
UV/Vis.: λ_{max}: 251, 295 nm
Supplied as: A crystalline 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

L-779,450 is supplied as a crystalline solid. A stock solution may be made by dissolving the L-779,450 in the solvent of choice, which should be purged with an inert gas. L-779,450 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of L-779,450 in ethanol is approximately 2 mg/ml and approximately 30 mg/ml in DMSO and DMF.

L-779,450 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, L-779,450 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. L-779,450 has a solubility of approximately 0.12 mg/ml in a 1:7 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

L-779,450 is an ATP-competitive B-Raf inhibitor (IC₅₀ = 10 nM; K_d = 2.4 nM) that displays >7, >30, and >70-fold selectivity over p38α, GSK3β, and Lck, respectively.¹ It has been shown to inhibit cell proliferation both in B-Raf mutated and wild-type melanoma cell lines, as well as to enhance tumor necrosis factor-related apoptosis-inducing ligand (TRAIL)-mediated apoptosis in these cells.²

References

1. Takle, A.K., Brown, M.J.B., Savies, S., *et al.* The identification of potent, selective and CNS penetrant furan-based inhibitors of B-Raf kinase. *Bioorg. Med. Chem. Lett.* **18(15)**, 4373-4376 (2008).
2. Berger, A., Quast, S.-A., Plötz, M., *et al.* RAF inhibition overcomes resistance to TRAIL-induced apoptosis in melanoma cells. *J. Invest. Dermatol.* **134(2)**, 430-440 (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