

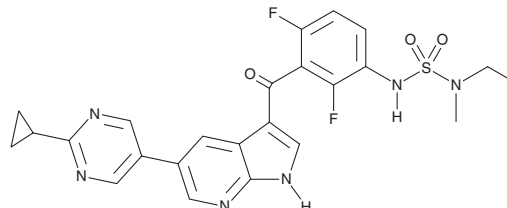
PRODUCT INFORMATION



PLX7904

Item No. 20710

CAS Registry No.: 1393465-84-3
Formal Name: N'-[3-[[5-(2-cyclopropyl-5-pyrimidinyl)-1H-pyrrolo[2,3-b]pyridin-3-yl]carbonyl]-2,4-difluorophenyl]-N-ethyl-N-methyl-sulfamide
MF: C₂₄H₂₂F₂N₆O₃S
FW: 512.5
Purity: ≥98%
UV/Vis.: λ_{max}: 258, 305 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

PLX7904 is supplied as a crystalline solid. A stock solution may be made by dissolving the PLX7904 in the solvent of choice, which should be purged with an inert gas. PLX7904 is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

Description

PLX7904 is a RAF inhibitor (IC₅₀s = 2.4, 140, and 91 nM for mutant B-RAF^{V600E}, wild-type B-RAF, and C-RAF, respectively).¹ It inhibits phosphorylation of ERK in A375 and COLO 829 cells (IC₅₀s = 16 and 18 nM, respectively). Unlike PLX4032 (Item No. 10618), BAY 43-9006 (Item No. 10009644), and dabrafenib (Item No. 16989), PLX7904 does not induce paradoxical pERK activation and proliferation of cancer cell lines (EC₅₀s = >200 μM). PLX7904 (25 mg/kg twice per day) inhibits tumor growth in a mouse COLO 205 colon cancer xenograft model.

Reference

1. Zhang, C., Spevak, W., Zhang, Y., et al. RAF inhibitors that evade paradoxical MAPK pathway activation. *Nature* **526**(7574), 583-586 (2015).

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, 12/01/2022

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