

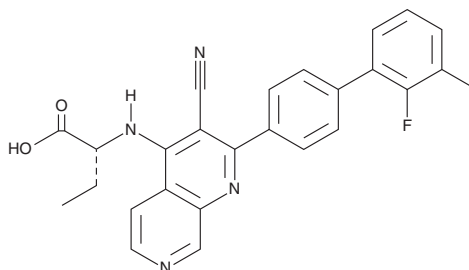
PRODUCT INFORMATION



BAY-091

Item No. 43634

CAS Registry No.: 2922280-34-8
Formal Name: (2R)-2-[[3-cyano-2-(2'-fluoro-3'-methyl[1,1'-biphenyl]-4-yl)-1,7-naphthyridin-4-yl]amino]-butanoic acid
MF: C₂₆H₂₁FN₄O₂
FW: 440.5
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

BAY-091 is supplied as a solid. A stock solution may be made by dissolving the BAY-091 in the solvent of choice, which should be purged with an inert gas. BAY-091 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

BAY-091 is an inhibitor of phosphatidylinositol 5-phosphate 4-kinase type-2 α (PIP4K2A; IC₅₀ = 1 nM).¹

Reference

1. Wortmann, L., Bräuer, N., Holton, S.J., *et al.* Discovery and characterization of the potent and highly selective 1,7-naphthyridine-based inhibitors BAY-091 and BAY-297 of the kinase PIP4K2A. *J. Med. Chem.* **64**(21), 15883-15911 (2021).

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, 08/21/2025

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