

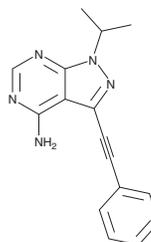
PRODUCT INFORMATION



SPP 86

Item No. 37390

CAS Registry No.: 1357349-91-7
Formal Name: 1-(1-methylethyl)-3-(2-phenylethynyl)-1H-pyrazolo[3,4-d]pyrimidin-4-amine
MF: C₁₆H₁₅N₅
FW: 277.3
Purity: ≥98%
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

SPP 86 is supplied as a solid. A stock solution may be made by dissolving the SPP 86 in the solvent of choice, which should be purged with an inert gas. SPP 86 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

SPP 86 is a RET inhibitor ($IC_{50} = 0.008 \mu\text{M}$).¹ It is selective for RET over a panel of 44 serine/threonine/tyrosine kinases at 10 μM but does inhibit nine by greater than 75%. SPP 86 is also selective for RET over a panel of 63 receptor and non-receptor tyrosine kinases at 0.4 μM but does inhibit six, including RET, by greater than 75%. SPP 86 reduces the proliferation of HCT116 colon cancer and TPC-1 thyroid carcinoma cells (EC_{50} s = 2.4 and 3.3 μM , respectively). It inhibits viral replication in BHK-21 cells infected with foot-and-mouth disease virus (FMDV) when used at a concentration of 5 μM .²

References

1. Dinér, P., Alao, J.P., Söderlund, J., *et al.* Preparation of 3-substituted-1-isopropyl-1H-pyrazolo[3,4-d]pyrimidin-4-amines as RET kinase inhibitors. *J. Med. Chem.* **55(10)**, 4872-4876 (2012).
2. Wang, X., Liao, Y., Abdullah, S.W., *et al.* FGFR1-mediated enhancement of foot-and-mouth disease virus entry. *Vet. Microbiol.* **298:110237**, (2024).

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, 06/13/2025

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