

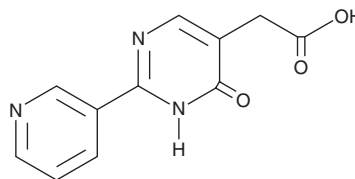
PRODUCT INFORMATION



CAY10781

Item No. 33778

CAS Registry No.: 1210360-87-4
Formal Name: 3,4-dihydro-4-oxo-2-(3-pyridinyl)-5-pyrimidineacetic acid
MF: C₁₁H₉N₃O₃
FW: 231.2
Purity: ≥90%
UV/Vis.: λ_{max}: 288 nm
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

CAY10781 is supplied as a solid. A stock solution may be made by dissolving the CAY10781 in the solvent of choice, which should be purged with an inert gas. CAY10781 is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Description

CAY10781 is an inhibitor of the protein-protein interaction between neuropilin-1 (NRP-1) and VEGF-A.¹ It inhibits the interaction by 43% when used at a concentration of 12.5 μM. CAY10781 also inhibits VEGF-A-induced phosphorylation of VEGFR2 in catecholamine A-differentiated (CAD) cells when used at the same concentration.

Reference

1. Perez-Miller, S., Patek, M., Moutal, A., *et al.* In silico identification and validation of inhibitors of the interaction between neuropilin receptor 1 and SARS-CoV-2 Spike protein. *bioRxiv Preprint*, 308783 (2020).

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, 11/15/2022

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