

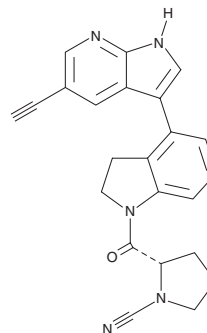
PRODUCT INFORMATION



IMP-1710

Item No. 31391

CAS Registry No.: 2383117-96-0
Formal Name: (2S)-2-[[4-(5-ethynyl-1H-pyrrolo[2,3-b]pyridin-3-yl)-2,3-dihydro-1H-indol-1-yl]carbonyl]-1-pyrrolidinecarbonitrile
Synonym: Click Tag™ IMP-1710
MF: C₂₃H₁₉N₅O
FW: 381.4
Purity: ≥98%
UV/Vis.: λ_{max}: 249 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

IMP-1710 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as chloroform purged with an inert gas can be used. The solubility of IMP-1710 in chloroform is approximately 10 mg/ml.

Description

IMP-1710 is an inhibitor of ubiquitin C-terminal hydrolase L1 (UCH-L1; IC₅₀ = 38 nM in a fluorescence polarization assay) that contains an alkyne moiety for use in click chemistry reactions.¹ It is selective for UCH-L1 over a panel of 20 deubiquitinating enzymes at 1 μM. IMP-1710 inhibits TGF-β1-induced fibroblast-to-myofibroblast transition (FMT) in primary lung fibroblasts isolated from patients with idiopathic pulmonary fibrosis (IC₅₀ = 740 nM). It has been used for labeling UCH-L1 in cell-based assays, followed by click reactions with azide-modified TAMRA and biotin capture reagents.

Reference

1. Panyain, N., Godinat, A., Lanyon-Hogg, T., *et al.* Discovery of a potent and selective covalent inhibitor and activity-based probe for the deubiquitylating enzyme UCHL1, with antifibrotic activity. *J. Am. Chem. Soc.* **142(28)**, 12020-12026 (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, 08/24/2023

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