

PRODUCT INFORMATION



MU1210

Item No. 41308

CAS Registry No.: 2275601-87-9

Formal Name: 5-(1-methyl-1H-pyrazol-4-yl)-3-[3-(4-pyridinyl)phenyl]-furo[3,2-b]pyridine

MF: C₂₂H₁₆N₄O

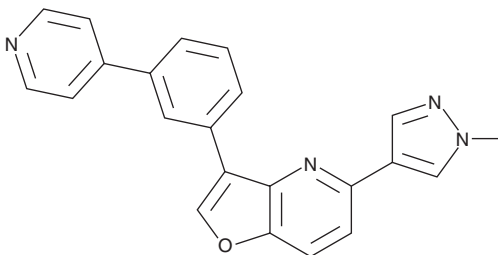
FW: 352.4

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

MU1210 is supplied as a solid. A stock solution may be made by dissolving the MU1210 in the solvent of choice, which should be purged with an inert gas. MU1210 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile and methanol.

Description

MU1210 is an inhibitor of CDC-like kinase 1 (Clk1), Clk2, and Clk4 (IC₅₀s = 8, 20, and 12 nM, respectively).¹ It is selective for these kinases over Clk3 (IC₅₀ = >3,000 nM), as well as dual-specificity tyrosine phosphorylation-regulated kinase 1A (DYRK1A), DYRK1B, DYRK2, homeodomain-interacting protein kinase 1 (HIPK1), and HIPK3 (IC₅₀s = 213, 956, 1,309, 187, and 159 nM, respectively) but does not inhibit HIPK2 (IC₅₀ = 23 nM).

Reference

1. Moyano, P.M., Némec, V., and Paruch, K. Cdc-Like kinases (CLKs): Biology, chemical probes, and therapeutic potential. *Int. J. Mol. Sci.* **21**(20), 7549 (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, 08/20/2024

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