

PRODUCT INFORMATION



M-110

Item No. 39165

CAS Registry No.: 1395048-49-3

Formal Name: 2-[[3-(4-morpholinyl)propyl] amino]-4-pyridinecarboxylic acid, 2-[1-(4-chloro-2-hydroxyphenyl) propylidene]hydrazide

MF: C₂₂H₂₈ClN₅O₃

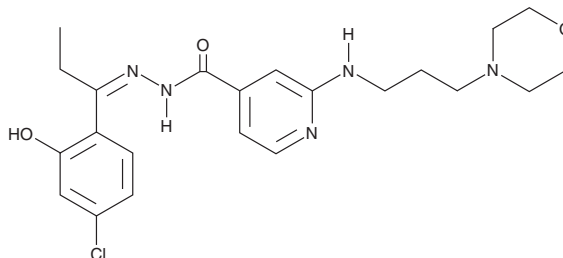
FW: 445.9

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

M-110 is supplied as a solid. A stock solution may be made by dissolving the M-110 in the solvent of choice, which should be purged with an inert gas. M-110 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO

Description

M-110 is a Pim-3 kinase inhibitor (IC₅₀ = 0.05 μM).¹ It is selective for Pim-3 over Pim-1 and Pim-2 (IC₅₀s = 2.5 μM for both). It inhibits the growth of DU145 and MIA PaCa-2 prostate cancer cells (IC₅₀s = 0.9 and 3.9 μM, respectively). M-110 (10 μM) inhibits IL-6-induced STAT3 phosphorylation in DU145 cells.

Reference

1. Chang, M., Kanwar, N., Feng, E., *et al.* PIM kinase inhibitors downregulate STAT3^{Tyr705} phosphorylation. *Mol. Cancer Ther.* **9**(9), 2478-2487 (2010).

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, 10/15/2024

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