

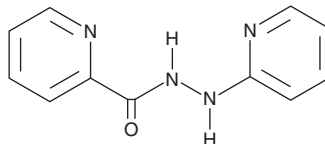
PRODUCT INFORMATION



PFI-90

Item No. 40120

CAS Registry No.: 53995-62-3
Formal Name: 2-(2-pyridinyl)hydrazide,
2-pyridinecarboxylic acid
MF: $C_{11}H_{10}N_4O$
FW: 214.2
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

PFI-90 is supplied as a solid. A stock solution may be made by dissolving the PFI-90 in the solvent of choice, which should be purged with an inert gas. PFI-90 is soluble in methanol.

Description

PFI-90 is an inhibitor of lysine demethylases (KDMs).¹ It inhibits KDM3B, KDM4B, KDM5A, and KDM6B by 46.8 to 87.1% when used at a concentration of 10 μ M. PFI-90 induces cell death in SJSA-1 osteosarcoma and TC-32 Ewing's sarcoma cancer cells (IC_{50} s = 2 and 1 μ M, respectively).

Reference

1. Khan, J., Hawley, R.G., Woldemichael, G.M., *et al.* Inhibitors of histone demethylases (PFI-63 and PFI-90) for the treatment of cancer and for the inhibition of histone demethylase in cells. **WO 2021/101929 A1** (2021).

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, 02/06/2024

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