

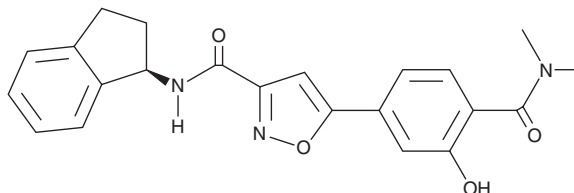
PRODUCT INFORMATION



PFI-6

Item No. 41162

CAS Registry No.: 2675452-91-0
Formal Name: N-[(1R)-2,3-dihydro-1H-inden-1-yl]-5-[4-[(dimethylamino)carbonyl]-3-hydroxyphenyl]-3-isoxazolecarboxamide
MF: C₂₂H₂₁N₃O₄
FW: 391.4
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

PFI-6 is supplied as a solid. A stock solution may be made by dissolving the PFI-6 in the solvent of choice, which should be purged with an inert gas. PFI-6 is slightly soluble (0.1-1 mg/ml) in acetonitrile and methanol, and sparingly soluble (1-10 mg/ml) in DMSO.

Description

PFI-6 is a chemical probe for the YEATS domain-containing proteins MLLT1 and MLLT3 (IC₅₀s = 140 and 160 nM, respectively).¹ It is selective for MLLT1 and MLLT3 over YEATS domain-containing 2 (YEATS2) and YEATS4 (IC₅₀s = >30,000 nM for both). See the Structural Genomics Consortium (SGC) website for more information.

Reference

1. Raux, B., Buchan, K.A., Bennett, J., *et al.* Discovery of PFI-6, a small-molecule chemical probe for the YEATS domain of MLLT1 and MLLT3. *Bioorg. Med. Chem. Lett.* **98**, 129546 (2024).

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, 08/07/2024

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