

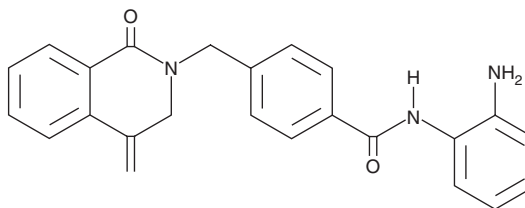
PRODUCT INFORMATION



MI-192

Item No. 18288

CAS Registry No.: 1415340-63-4
Formal Name: N-(2-aminophenyl)-4-[(3,4-dihydro-4-methylene-1-oxo-2(1H)-isoquinolinyl)methyl]-benzamide
MF: C₂₄H₂₁N₃O₂
FW: 383.4
Purity: ≥98%
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 203, 234 nm
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

MI-192 is supplied as a crystalline solid. A stock solution may be made by dissolving the MI-192 in the solvent of choice. MI-192 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of MI-192 in these solvents is approximately 5 and 1 mg/ml, respectively.

MI-192 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MI-192 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. MI-192 has a solubility of approximately 0.1 mg/ml in a 1:7 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

MI-192 is an inhibitor of histone deacetylases (HDACs) that preferentially inhibits HDAC2 (IC₅₀ = 30 nM) and HDAC3 (IC₅₀ = 16 nM) over HDAC1, 4, 6, 7, and 8 (IC₅₀s = 4.8, 5, >10, 4.1, and >10 μM, respectively).¹ At 0.1-0.4 μM, it induces differentiation and promotes apoptosis of acute myeloid leukemic cell lines U937, HL60, and Kasumi-1.¹

Reference

1. Boissinot, M., Inman, M., Hempshall, A., *et al.* Induction of differentiation and apoptosis in leukaemic cell lines by the novel benzamide family histone deacetylase 2 and 3 inhibitor MI-192. *Leuk. Res.* **36**, 1304-1310 (2012).

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, 12/02/2022

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