

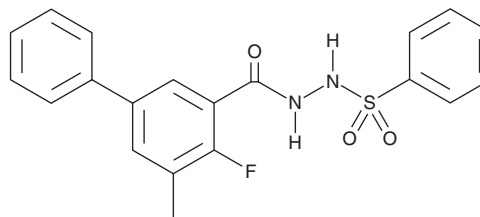
# PRODUCT INFORMATION



## MOZ-IN-3

Item No. 27402

**CAS Registry No.:** 2055397-18-5  
**Formal Name:** 4-fluoro-5-methyl-[1,1'-biphenyl]-3-carboxylic acid, 2-(phenylsulfonyl)hydrazide  
**Synonym:** WM-8014  
**MF:** C<sub>20</sub>H<sub>17</sub>FN<sub>2</sub>O<sub>3</sub>S  
**FW:** 384.4  
**Purity:** ≥98%  
**Supplied as:** A crystalline 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

MOZ-IN-3 is supplied as a crystalline solid. A stock solution may be made by dissolving the MOZ-IN-3 in the solvent of choice, which should be purged with an inert gas. MOZ-IN-3 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of MOZ-IN-3 in these solvents is approximately 2, 33, and 50 mg/ml, respectively.

MOZ-IN-3 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MOZ-IN-3 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. MOZ-IN-3 has a solubility of approximately 0.1 mg/ml in a 1:7 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

MOZ-IN-3 is an inhibitor of lysine acetyltransferase 6A (KAT6A/MOZ) and KAT6B/MORF (IC<sub>50</sub>s = 8 and 28 nM, respectively).<sup>1</sup> It is selective for KAT6A and KAT6B over KAT5 and KAT7 (IC<sub>50</sub>s = 224 and 342 nM, respectively). MOZ-IN-3 induces cell cycle arrest in and inhibits proliferation of mouse embryonic fibroblasts (IC<sub>50</sub> = 2.4 μM) and inhibits proliferation of EMRK1184 lymphoma cells (IC<sub>50</sub> = 2.3 μM). It induces genetic and morphological changes associated with cellular senescence without inducing DNA damage, apoptosis, or necrosis. MOZ-IN-3 (10 μM) decreases global levels of acetylated histone H3 lysine 14 (H3K14Ac) by 49% and reduces H3K9Ac levels at the transcription start sites of *Ezh2*, *Mek*, and *E2f2*.

### Reference

1. Baell, J.B., Leaver, D.J., Hermans, S.J., *et al.* Inhibitors of histone acetyltransferases KAT6A/B induce senescence and arrest tumour growth. *Nature* **560(7717)**, 253-257 (2018).

#### 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, 12/19/2022

#### 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