

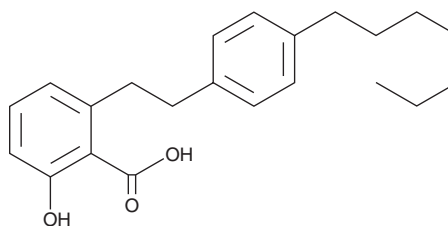
# PRODUCT INFORMATION



## MG149

Item No. 22135

**CAS Registry No.:** 1243583-85-8  
**Formal Name:** 2-[2-(4-heptylphenyl)ethyl]-6-hydroxy-benzoic acid  
**MF:** C<sub>22</sub>H<sub>28</sub>O<sub>3</sub>  
**FW:** 340.5  
**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

MG149 is supplied as a crystalline solid. A stock solution may be made by dissolving the MG149 in the solvent of choice, which should be purged with an inert gas. MG149 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of MG149 in these solvents is approximately 30 mg/ml.

MG149 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MG149 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. MG149 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

MG149 is a histone acetyltransferase (HAT) inhibitor with IC<sub>50</sub> values of 74 and 47 μM for human recombinant Tip60 and males absent on the first (MOF) HATs, respectively.<sup>1,2</sup> MG149 targets the acetyl-CoA binding site and inhibits acetyltransferase activity in the nuclear extract from murine hippocampus, amygdala, and prefrontal cortex.<sup>3</sup> MG149 also inhibits the p53 and NF-κB pathways.

### References

1. Legartová, S., Stixová, L., Strnad, H., *et al.* Basic nuclear processes affected by histone acetyltransferases and histone deacetylase inhibitors. *Epigenomics* **5(4)**, 379-396 (2013).
2. Su, J., Wang, F., Cai, Y., *et al.* The functional analysis of histone acetyltransferase MOF in tumorigenesis. *Int. J. Mol. Sci.* **17(1)**, (2016).
3. Ghizzoni, M., Wu, J., Gao, T., *et al.* 6-alkylsalicylates are selective Tip60 inhibitors and target the acetyl-CoA binding site. *Eur. J. Med. Chem.* **47(1)**, 337-344 (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