

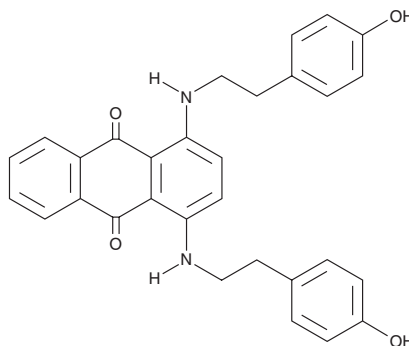
PRODUCT INFORMATION



JFD00244

Item No. 14648

CAS Registry No.: 96969-83-4
Formal Name: 1,4-bis[[2-(4-hydroxyphenyl)ethyl]amino]-9,10-anthracenedione
Synonym: BML-266
MF: C₃₀H₂₆N₂O₄
FW: 478.5
Purity: ≥95%
UV/Vis.: λ_{max}: 226, 258, 596, 642 nm
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

JFD00244 is supplied as a crystalline solid. A stock solution may be made by dissolving the JFD00244 in the solvent of choice, which should be purged with an inert gas. JFD00244 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of JFD00244 in ethanol is approximately 0.3 mg/ml and approximately 30 mg/ml in DMSO and DMF.

JFD00244 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JFD00244 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. JFD00244 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Human sirtuin 2 (SIRT2) is a NAD⁺-dependent protein deacetylase with diverse targets, including α-tubulin, phosphoenolpyruvate carboxykinase, and forkhead box protein O1. JFD00244 is an inhibitor of SIRT2 (IC₅₀ = 56.7 μM).¹ At 5 μM, JFD00244 induces granulocytic differentiation in the acute promyelocytic leukemia cell line NB4.²

References

1. Kiviranta, P.H., Leppänen, J., Kyrylenko, S., *et al.* N,N'-Bisbenzylidenebenzene-1,4-diamines and N,N'-bisbenzylidenenaphthalene-1,4-diamines as sirtuin type 2 (SIRT2) inhibitors. *J. Med. Chem.* **49**(26), 7907-7911 (2006).
2. Sunami, Y., Araki, M., Hironaka, Y., *et al.* Inhibition of the NAD-dependent protein deacetylase SIRT2 induces granulocytic differentiation in human leukemia cells. *PLoS One* **8**(2), 57633 (2013).

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, 11/11/2022

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