

PRODUCT INFORMATION



BH400

Item No. 41892

Formal Name: 5-(((4'-fluoro-[1,1'-biphenyl]-4-yl)methyl)amino)-2-hydroxybenzoic acid

MF: C₂₀H₁₆FNO₃

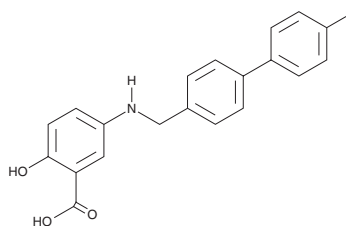
FW: 337.4

Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

BH400 is supplied as a solid. A stock solution may be made by dissolving the BH400 in the solvent of choice, which should be purged with an inert gas. BH400 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

BH400 is an agonist of peroxisome proliferator-activated receptor α (PPAR α) and an inhibitor of stimulator of interferon genes (STING; IC₅₀ = 8.1 μ M).¹ It induces transactivation of PPAR α in a reporter assay using THP-1 cells (EC₅₀ = 1.2 μ M). BH400 (10 μ M) inhibits decreases in viability induced by 4-hydroxy nonenal (4-HNE; Item No. 32100) or hydrogen peroxide (H₂O₂) in 661W retinal ganglion cells. It prevents LPS-induced increases in phosphorylated TANK-binding kinase 1 (TBK1) and IL-6 levels and LPS-induced decreases in peroxisome proliferator-activated receptor- γ coactivator-1 α (PGC-1 α) levels in HMC3 microglial cells when used at a concentration of 10 μ M.

Reference

1. Hu, B., Cui, Y., Lee, J.J., *et al.* Design and assessment of first-generation heterobifunctional PPAR α /STING modulators. *ACS Med. Chem. Lett.* **15**(8), 1279-1286 (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, 10/24/2024

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