

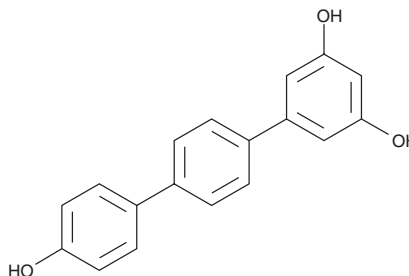
PRODUCT INFORMATION



CAY10503

Item No. 10008872

CAS Registry No.: 890854-82-7
Formal Name: 4-(4-hydroxyphenyl)phenyl-3,5-benzenediol
MF: C₁₈H₁₄O₃
FW: 278.3
Purity: ≥98%
UV/Vis.: λ_{max}: 207, 291 nm
Supplied as: A crystalline 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

CAY10503 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10503 in the solvent of choice, which should be purged with an inert gas. CAY10503 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10503 in ethanol and DMF is approximately 30 mg/ml and approximately 20 mg/ml in DMSO.

CAY10503 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10503 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. CAY10503 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

CAY10503 is a proapoptotic, antiproliferative compound that is able to arrest cell cycle progression in the G₀-G₁ phase.¹ CAY10503 inhibits the growth of HL60, multidrug resistant HL60R, and K562 cells with IC₅₀ values of 7.0, 23, and 20 μM, respectively. Accumulation of HL60 cells in G₀-G₁ occurred within eight hours following treatment with 50 μM CAY10503, whereas 10 μM CAY10503 required 96 hours for cell cycle arrest to appear. CAY10503 also induces differentiation of HL60 cells into the following positive cells: CD14 (monocytic marker) as well as CD11b and CD11c (granulocytic markers). Approximately 60% of HL60 cells exposed to 10 μM CAY10503 expressed these markers within 72 hours.¹

Reference

1. Roberti, M., Pizzirani, D., Recanatini, M., *et al.* Identification of a terphenyl derivative that blocks the cell cycle in the G₀-G₁ phase and induces differentiation in leukemia cells. *J. Med. Chem.* **49(10)**, 3012-3018 (2006).

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, 06/02/2023

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