

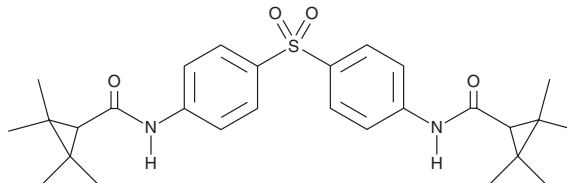
PRODUCT INFORMATION



BRD7116

Item No. 20570

CAS Registry No.: 329059-55-4
Formal Name: N,N'-(sulfonyldi-4,1-phenylene)
bis[2,2,3,3-tetramethyl-
cyclopropanecarboxamide]
MF: C₂₈H₃₆N₂O₄S
FW: 496.7
Purity: ≥98%
UV/Vis.: λ_{max}: 263, 292 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

BRD7116 is supplied as a crystalline solid. A stock solution may be made by dissolving the BRD7116 in the solvent of choice, which should be purged with an inert gas. BRD7116 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of BRD7116 in these solvents is approximately 10 mg/ml.

BRD7116 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BRD7116 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BRD7116 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BRD7116 is a *bis*-aryl sulfone that acts as a selective inhibitor of leukemia stem cell activity. It exhibits an EC₅₀ value of 200 nM for leukemia stem cells isolated from the bone marrow of leukemic animals in co-culture but does not affect normal hematopoietic stem cells or AML cell lines (EC₅₀s = 20 μM).¹ Its exact mechanism is not well understood but is thought to involve impairment of the stroma's ability to support leukemia stem cells by inducing transcriptional changes consistent with myeloid differentiation.¹

Reference

1. Hartwell, K.A., Miller, P.G., Mukherjee, S., *et al.* Niche-based screening identifies small-molecule inhibitors of leukemia stem cells. *Nat. Chem. Biol.* **9**(12), 840-848 (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, 12/16/2022

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