

# PRODUCT INFORMATION



## BAY-299

Item No. 19777

**CAS Registry No.:** 2080306-23-4  
**Formal Name:** 2-(2,3-dihydro-1,3,6-trimethyl-2-oxo-1H-benzimidazol-5-yl)-6-(3-hydroxypropyl)-1H-benz[de]isoquinoline-1,3(2H)-dione

**MF:** C<sub>25</sub>H<sub>23</sub>N<sub>3</sub>O<sub>4</sub>

**FW:** 429.5

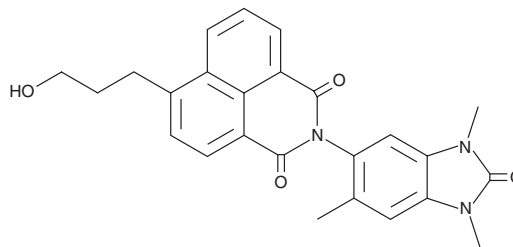
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 215, 236, 294, 342 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

BAY-299 is supplied as a crystalline solid. A stock solution may be made by dissolving the BAY-299 in the solvent of choice, which should be purged with an inert gas. BAY-299 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of BAY-299 in these solvents is approximately 20 and 30 mg/ml, respectively.

### Description

BAY-299 is a potent and selective inhibitor of the bromodomain and PHD finger-containing (BRPF) family protein BRD1 (IC<sub>50</sub> = 6 nM), also known as BRPF2, and the second bromodomain of transcription initiation factor TFIID subunits 1 (TAF1; IC<sub>50</sub> = 13 nM).<sup>1</sup> BAY-299 is >30-fold selective over BRPF1, BRPF3, BRD9, and ATAD2 and is >300-fold selective over BRD4. See the Structural Genomics Consortium (SGC) website for more information.

### Reference

1. Pérez-Salvia, M., and Esteller, M. Bromodomain inhibitors and cancer therapy: From structures to applications. *Epigenetics* (2016).

#### 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, 12/22/2022

#### 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