

# PRODUCT INFORMATION

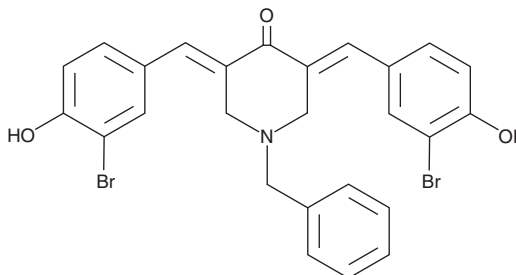


## PRMT4/CARM1 Inhibitor

Item No. 11033

**CAS Registry No.:** 1020399-49-8  
**Formal Name:** 3,5-bis[(3-bromo-4-hydroxyphenyl)methylene]-1-(phenylmethyl)-4-piperidinone  
**Synonym:** Protein Arginine Methyltransferase 4/Coactivator-Associated Arginine Methyltransferase 1 Inhibitor

**MF:** C<sub>26</sub>H<sub>21</sub>Br<sub>2</sub>NO<sub>3</sub>  
**FW:** 555.3  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 250, 367 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

PRMT4/CARM1 inhibitor is supplied as a crystalline solid. A stock solution may be made by dissolving the PRMT4/CARM1 inhibitor in the solvent of choice, which should be purged with an inert gas. PRMT4/CARM1 inhibitor is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PRMT4/CARM1 inhibitor in ethanol is approximately 0.25 mg/ml and approximately 30 mg/ml in DMSO and DMF.

PRMT4/CARM1 inhibitor is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PRMT4/CARM1 inhibitor should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. PRMT4/CARM1 inhibitor has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

PRMT4/CARM1 Inhibitor selectively blocks the activity of PRMT4/CARM1 with an IC<sub>50</sub> value of 7.1 μM.<sup>1</sup> It displays very low affinity towards PRMT1 and SET7 with IC<sub>50</sub> values of 63 and 943 μM, respectively.<sup>1</sup> In human prostate cancer LNCaP cells, this compound reduced the prostate-specific antigen promoter activity at 8-10 μM.<sup>1</sup>

### Reference

1. Cheng, D., Valente, S., Castellano, S., *et al.* Novel 3,5-bis(bromohydroxybenzylidene)piperidin-4-ones as coactivator-associated arginine methyltransferase 1 inhibitors: Enzyme selectivity and cellular activity. *J. Med. Chem.* **54**(13), 4982-4932 (2011).

#### 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, 11/02/2022

#### 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