

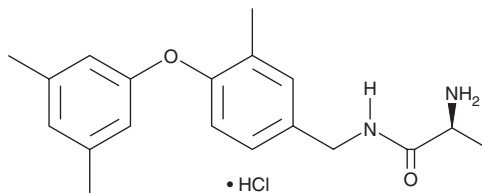
# PRODUCT INFORMATION



## SGC2085 (hydrochloride)

Item No. 21879

**CAS Registry No.:** 1821908-49-9  
**Formal Name:** 2S-amino-N-[[4-(3,5-dimethylphenoxy)-3-methylphenyl]methyl]-propanamide, monohydrochloride  
**MF:** C<sub>19</sub>H<sub>24</sub>N<sub>2</sub>O<sub>2</sub> • HCl  
**FW:** 348.9  
**Purity:** ≥98%  
**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

SGC2085 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the SGC2085 (hydrochloride) in the solvent of choice. SGC2085 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of SGC2085 (hydrochloride) in these solvents is approximately 30 mg/ml.

SGC2085 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SGC2085 (hydrochloride) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. SGC2085 (hydrochloride) has a solubility of approximately 0.1 mg/ml in a 1:10 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

SGC2085 is an inhibitor of protein arginine methyltransferase 4 (PRMT4/CARM1; IC<sub>50</sub> = 50 nM).<sup>1</sup> It is selective for PRMT4/CARM1 over PRMT8 at concentrations up to 50 μM and PRMT1, PRMT3, PRMT5, and PRMT7 up to 100 μM. SGC2085 also weakly inhibits PRMT6 (IC<sub>50</sub> = 5.2 μM). It has no activity in HEK293 cells at concentrations up to 10 μM.

### Reference

1. Ferreira de Freitas, R., Eram, M.S., Smil, D., *et al.* Discovery of a potent and selective coactivator associated arginine methyltransferase 1 (CARM1) inhibitor by virtual screening. *J. Med. Chem.* **59**(14), 6838-6847 (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