

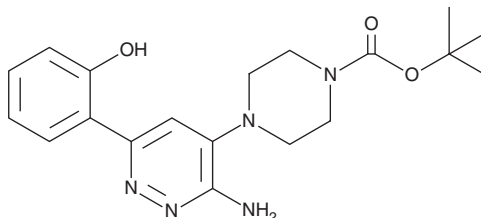
# PRODUCT INFORMATION



## SGC-SMARCA-BRDVIII

Item No. 33567

**CAS Registry No.:** 1997319-84-2  
**Formal Name:** 4-[3-amino-6-(2-hydroxyphenyl)-4-pyridazinyl]-1-piperazinecarboxylic acid, 1,1-dimethylethyl ester  
**MF:** C<sub>19</sub>H<sub>25</sub>N<sub>5</sub>O<sub>3</sub>  
**FW:** 371.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 219, 265, 313 nm  
**Supplied as:** A 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

SGC-SMARCA-BRDVIII is supplied as a solid. A stock solution may be made by dissolving the SGC-SMARCA-BRDVIII in the solvent of choice, which should be purged with an inert gas. SGC-SMARCA-BRDVIII is soluble in DMSO.

### Description

SGC-SMARCA-BRDVIII is a pan-inhibitor of family VIII bromodomain-containing proteins (BRDs), which are components of switch/sucrose nonfermentable (SWI/SNF) chromatin remodeling complexes.<sup>1</sup> It selectively binds to polybromo-1D (PBRM1) bromodomain 5, BRM, also known as SMARCA2, and BRG1, also known as SMARCA4 (K<sub>d</sub>s = 13, 35, and 36 nM, respectively), over PBRM1 bromodomain 2, PBRM1 bromodomain 3, and PBRM1 bromodomain 4 (K<sub>d</sub>s = 3,655, 1,963, and >10,000 nM, respectively). SGC-SMARCA-BRDVIII (1 μM) decreases lipid accumulation induced by the PPARγ agonist rosiglitazone (Item Nos. 71740 | 11884 | 71742) in 3T3-L1 mouse fibroblasts differentiated into adipocytes by a combination of the phosphodiesterase (PDE) inhibitor IBMX (Item No. 13347), the glucocorticoid dexamethasone (Item No. 11015), and insulin. It decreases expression of the genes encoding PPARγ, C/EBPα, and fatty acid binding protein 4 (FABP4), markers of adipocytes, in the same cells when used at a concentration of 0.5 μM. See the Structural Genomics Consortium (SGC) website for more information.

### Reference

1. Wanior, M., Preuss, F., Ni, X., *et al.* Pan-SMARCA/PB1 bromodomain inhibitors and their role in regulating adipogenesis. *J. Med. Chem.* **63**(23), 14680-14699 (2020).

**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, 04/20/2023

### 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