

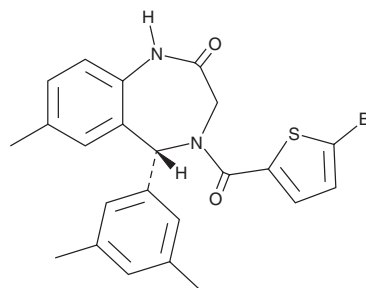
PRODUCT INFORMATION



SGC6870

Item No. 33712

CAS Registry No.: 2561471-27-8
Formal Name: 4-[(5R-bromo-2-thienyl)carbonyl]-5-(3,5-dimethylphenyl)-1,3,4,5-tetrahydro-7-methyl-2H-1,4-benzodiazepin-2-one
MF: C₂₃H₂₁BrN₂O₂S
FW: 469.4
Purity: ≥98%
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

SGC6870 is supplied as a solid. A stock solution may be made by dissolving the SGC6870 in the solvent of choice, which should be purged with an inert gas. SGC6870 is soluble in ethanol and DMSO.

Description

SGC6870 is an inhibitor of protein arginine methyltransferase 6 (PRMT6; IC₅₀ = 77 nM).¹ It is selective for PRMT6 over a panel of 76 methyltransferases, G protein-coupled receptors (GPCRs), ion channels, receptors, and other enzymes at 1 μM. SGC6870 decreases the levels of dimethylated arginine 2 on histone 3 (H3R2me2) and dimethylated arginine 3 on histone 4 (H4R3me2) in HEK293T cells (IC₅₀s = 0.9 and 0.6 μM, respectively). See the Structural Genomics Consortium (SGC) website for more information.

Reference

1. Shen, Y., Li, F., Szewczyk, M.M., et al. A first-in-class, highly selective and cell-active allosteric inhibitor of protein arginine methyltransferase 6. *J. Med. Chem.* **64**(7), 3697-3706 (2021).

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, 06/19/2024

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