

PRODUCT INFORMATION



SGC3027

Item No. 25325

CAS Registry No.: 2624313-13-7
Formal Name: N-(4-(((2S,3S,4R,5R)-5-(6-amino-9H-purin-9-yl)-3,4-dihydroxytetrahydrofuran-2-yl)methyl)thio)butyl)-N-((4'-chloro-[1,1'-biphenyl]-3-yl)methyl)-3-methyl-3-(2,4,5-trimethyl-3,6-dioxocyclohexa-1,4-dien-1-yl)butanamide

MF: C₄₁H₄₇ClN₆O₆S

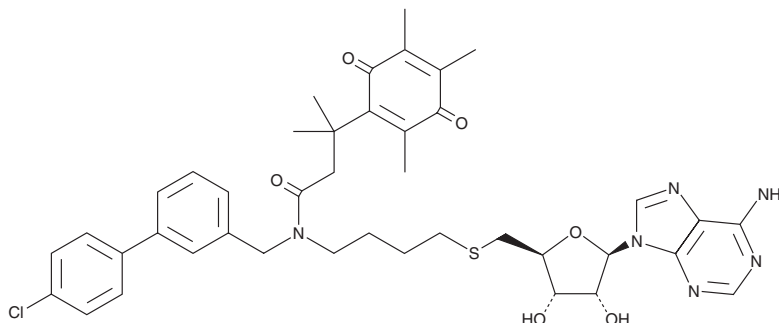
FW: 787.4

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

SGC3027 is supplied as a solid. A stock solution may be made by dissolving the SGC3027 in the solvent of choice, which should be purged with an inert gas. SGC3027 is soluble in the organic solvent DMSO.

Description

SGC3027 is a prodrug form of the protein arginine methyltransferase 7 (PRMT7) inhibitor SGC8158. It is converted by intracellular reductases to SGC8158, which inhibits PRMT7 with an IC₅₀ value of less than 2.5 nM for inhibition of histone H2B (23-37) methylation. SGC3027 inhibits methylation of heat shock protein 70 (Hsp70) *in vitro* (IC₅₀ = 2.4 μM). See the Structural Genomics Consortium (SGC) website for more information.

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, 01/04/2023

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