

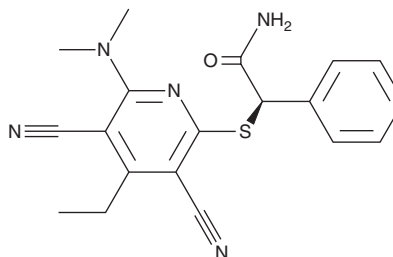
PRODUCT INFORMATION



GSK3484862

Item No. 38897

CAS Registry No.: 2170136-65-7
Formal Name: αR-[[3,5-dicyano-6-(dimethylamino)-4-ethyl-2-pyridinyl]thio]-benzeneacetamide
MF: C₁₉H₁₉N₅OS
FW: 365.5
Purity: ≥95%
UV/Vis.: λ_{max}: 270, 304, 347 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

GSK3484862 is supplied as a solid. A stock solution may be made by dissolving the GSK3484862 in the solvent of choice, which should be purged with an inert gas. GSK3484862 is soluble in DMSO.

Description

GSK3484862 is an inhibitor of DNA methyltransferase 1 (DNMT1; IC₅₀ = 0.23 μM).¹ It is selective for DNMT1 over DNMT3A/3L and DNMT3B/3L (IC₅₀s = >50 μM for both). GSK3484862 (100-1,000 nM) inhibits the growth of MV4-11 leukemia cells. It also decreases promoter DNA methylation in MV4-11 cells.

Reference

1. Pappalardi, M.B., Keenan, K., Cockerill, M., et al. Discovery of a first-in-class reversible DNMT1-selective inhibitor with improved tolerability and efficacy in acute myeloid leukemia. *Nat. Cancer* **2(10)**, 1002-1017 (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, 09/07/2023

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