

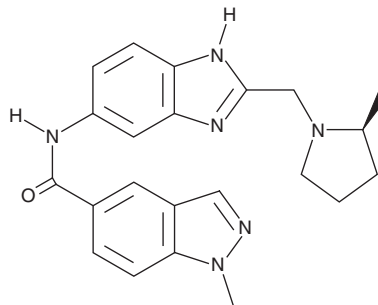
PRODUCT INFORMATION



SGC-iMLLT

Item No. 28113

CAS Registry No.: 2255338-25-9
Formal Name: 1-methyl-N-[2-[[[(2S)-2-methyl-1-pyrrolidinyl]methyl]-1H-benzimidazol-6-yl]-1H-indazole-5-carboxamide
MF: C₂₂H₂₄N₆O
FW: 388.5
Purity: ≥95%
UV/Vis.: λ_{max}: 234, 295 nm
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

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

Description

SGC-iMLLT is an inhibitor of the YEATS domain-containing proteins MLLT1 and MLLT3.¹ It binds to the MLLT1 YEATS1 and MLLT3 YEATS3 domains (K_ds = 129 and 77 nM, respectively).¹ It is selective for YEATS1 and -3 over YEATS2 and -4 (IC₅₀s = >10 μM), as well as a panel of 48 bromodomains at 50 μM. SGC-iMLLT inhibits the interaction of histone H3.3 with MLLT3 (IC₅₀ = 400 nM in a reporter assay) and decreases expression of the tumorigenic genes MYC and DDN in MV-4-11 leukemia cells when used at a concentration of 1 μM.

Reference

1. Moustakim, M., Christott, T., Monteiro, O.P., *et al.* Discovery of an MLLT1/3 YEATS domain chemical probe. *Angew. Chem. Int. Ed. Engl.* **57**(50), 16302-16307 (2018).

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, 02/27/2026

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