

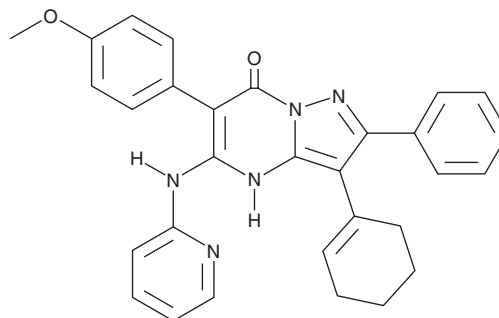
PRODUCT INFORMATION



AG-270

Item No. 36246

CAS Registry No.: 2201056-66-6
Formal Name: 3-(1-cyclohexen-1-yl)-6-(4-methoxyphenyl)-2-phenyl-5-(2-pyridinylamino)-pyrazolo[1,5-a]pyrimidin-7(4H)-one
MF: C₃₀H₂₇N₅O₂
FW: 489.6
Purity: ≥95%
UV/Vis.: λ_{max}: 273, 317 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

AG-270 is supplied as a solid. A stock solution may be made by dissolving the AG-270 in the solvent of choice, which should be purged with an inert gas. AG-270 is slightly soluble in DMSO.

Description

AG-270 is an inhibitor of methionine adenosyltransferase 2A (MAT2A; IC₅₀ = 0.014 μM).¹ It reduces levels of S-(5'-adenosyl)-L-methionine (SAM; Item Nos. 16376 | 17595) in, and inhibits the proliferation of, HCT116 cells with a homozygous deletion of *MTAP*, the gene encoding methylthioadenosine phosphorylase (IC₅₀s = 0.2 and 0.257 μM, respectively). AG-270 (200 mg/kg) reduces intratumor SAM levels and tumor growth in KP4 and HCT116 *MTAP*-null, but not HCT116 *MTAP*-wild-type, mouse xenograft models.

Reference

1. Konteatis, Z., Travins, J., Gross, S., Marjon, K., *et al.* Discovery of AG-270, a first-in-class oral MAT2A inhibitor for the treatment of tumors with homozygous *MTAP* deletion. *J. Med. Chem.* **64**(8), 4430-4449 (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, 08/16/2024

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