

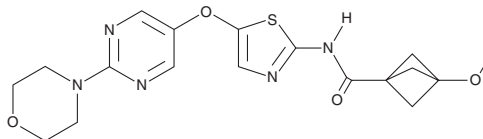
PRODUCT INFORMATION



RGN6024

Item No. 45752

CAS Registry No.: 3055038-20-2
Formal Name: 3-methoxy-N-[5-[[2-(4-morpholinyl)-5-pyrimidinyl]oxy]-2-thiazolyl]-bicyclo[1.1.1]pentane-1-carboxamide
MF: C₁₈H₂₁N₅O₄S
FW: 403.5
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

RGN6024 is supplied as a solid. A stock solution may be made by dissolving the RGN6024 in the solvent of choice, which should be purged with an inert gas. RGN6024 is soluble (≥20 mg/ml) in DMSO.

Description

RGN6024 is a tubulin destabilizer.¹ It reduces tumor growth in a temozolomide-resistant LN-18 glioblastoma mouse xenograft model when administered orally at doses of 7.5 or 15 mg/kg.

Reference

1. Patrón, L.A., Yeoman, H., Ramos, J., *et al.* RGN6024 is a brain-penetrant, small-molecule tubulin destabilizer for the treatment of glioblastoma. *Mol. Cancer Ther.* **24**(8), 1129-1144 (2025).

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, 06/11/2026

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