

PRODUCT INFORMATION



RLY-2608

Item No. 44472

CAS Registry No.: 2733573-94-7
Formal Name: N-[(3R)-3-(2-chloro-5-fluorophenyl)-6-(5-cyano[1,2,4] triazolo[1,5-a]pyridin-6-yl)-2,3-dihydro-1-oxo-1H-isoindol-4-yl]-3-fluoro-5-(trifluoromethyl)-benzamide

MF: C₂₉H₁₄ClF₅N₆O₂

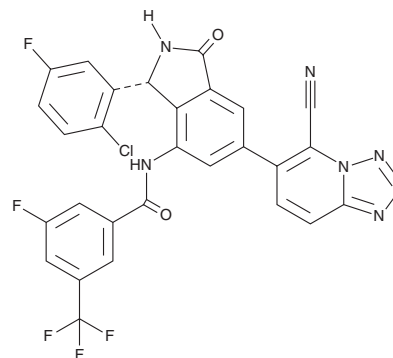
FW: 608.9

Purity: ≥95%

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

RLY-2608 is supplied as a solid. A stock solution may be made by dissolving the RLY-2608 in the solvent of choice, which should be purged with an inert gas. RLY-2608 is soluble (≥10 mg/ml) in DMSO.

Description

RLY-2608 is an inhibitor of PI3Kα containing a histidine-to-arginine mutation at position 1,047 (PI3Kα^{H1047R}; IC₅₀ = 4 nM).¹ It is selective for PI3Kα^{H1047R} over wild-type PI3Kα (IC₅₀ = 48 nM). RLY-2608 (25 and 100 mg/kg) reduces tumor volume in T47D and ST1056 PI3Kα^{H1047R} mutant-containing breast cancer mouse xenograft models.

Reference

1. Varkaris, A., Pazolli, E., Gunaydin, H., *et al.* Discovery and clinical proof-of-concept of RLY-2608, a first-in-class mutant-selective allosteric PI3Kα inhibitor that decouples antitumor activity from hyperinsulinemia. *Cancer Discov.* **14**(2), 240-257 (2024).

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, 10/22/2025

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