

PRODUCT INFORMATION



GDC-1971

Item No. 39884

CAS Registry No.: 2377352-49-1

Formal Name: (3R)-1'-[3-(3,4-dihydro-1,5-naphthyridin-1(2H)-yl)-1H-pyrazolo[3,4-b]pyrazin-6-yl]-spiro[benzofuran-2(3H),4'-piperidin]-3-amine

Synonym: Migoprotafib

MF: C₂₅H₂₆N₈O

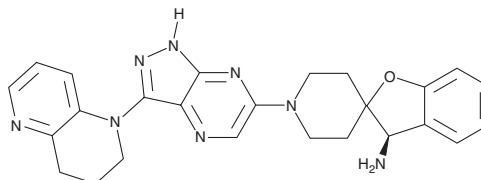
FW: 454.5

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

GDC-1971 is supplied as a solid. A stock solution may be made by dissolving the GDC-1971 in the solvent of choice, which should be purged with an inert gas. GDC-1971 is sparingly soluble (1-10 mg/ml) in methanol and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

GDC-1971 is an inhibitor of Src homology region 2 domain-containing phosphatase 2 (SHP-2; IC₅₀ = <0.001 μM).¹ It is selective for SHP-2 over SHP-2 containing a glutamate-to-lysine substitution at position 76 (SHP-2^{E76K}; IC₅₀ = 0.17 μM). GDC-1971 reduces the levels of phosphorylated ERK in a reporter assay using KYSE-520 esophageal cancer cells (IC₉₀ = 0.012 μM) and inhibits proliferation of the same cells (IC₅₀ = 0.32 μM). *In vivo*, GDC-1971 (30 or 100 mg/kg two times per day) reduces tumor volume in KYSE-520 and NCI H358 non-small cell lung cancer (NSCLC) mouse xenograft models.

Reference

1. Taylor, A.M., Williams, B.R., Giordanetto, F., et al. Identification of GDC-1971 (RLY-1971), a SHP2 inhibitor designed for the treatment of solid tumors. *J. Med. Chem.* **66**(19), 13384-13399 (2023).

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, 05/31/2024

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