

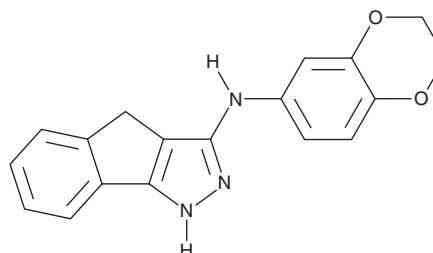
PRODUCT INFORMATION



GN44028

Item No. 37289

CAS Registry No.: 1421448-26-1
Formal Name: N-(2,3-dihydro-1,4-benzodioxin-6-yl)-1,4-dihydro-indeno[1,2-c]pyrazol-3-amine
MF: C₁₈H₁₅N₃O₂
FW: 305.3
Purity: ≥98%
UV/Vis.: λ_{max}: 264 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

GN44028 is supplied as a solid. A stock solution may be made by dissolving the GN44028 in the solvent of choice, which should be purged with an inert gas. GN44028 is soluble in methanol.

Description

GN44028 is an inhibitor of HIF-1α transcriptional activity.¹ It reduces HIF-1α reporter gene expression in HeLa cervical cancer cells (IC₅₀ = 14 nM). GN44028 induces cytotoxicity in HCT116 colorectal, HepG2 hepatic, PC3 prostate, and HeLa cancer cells (IC₅₀s = 2.1, 3.7, 25.4, and 2.1 μM, respectively). It reduces the levels of VEGF, but not HIF-1α or HIF-1α-HIF-1β dimer levels, in HeLa cells when used at concentrations of 0.1 and 1 μM. *In vivo*, GN44028 (5 mg/kg twice per week), alone or in combination with 5-fluorouracil (Item. No. 14416), decreases tumor volume and the number of lung metastases and increases survival percentage in a CT26 murine colon cancer model.²

References

1. Minegishi, H., Fukushima, S., Ban, H.S., *et al.* Discovery of indenopyrazoles as a new class of hypoxia inducible factor (HIF)-1 inhibitors. *ACS Med. Chem. Lett.* **4**(2), 297-301 (2013).
2. Liu, C., Zhang, W., Wang, J., *et al.* Tumor-associated macrophage-derived transforming growth factor-β promotes colorectal cancer progression through HIF1-TRIB3 signaling. *Cancer Sci.* **112**(10), 4198-4207 (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, 10/09/2023

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