

PRODUCT INFORMATION



GSK1904529A

Item No. 29155

CAS Registry No.: 1089283-49-7
Formal Name: N-(2,6-difluorophenyl)-5-[3-[2-[[5-ethyl-2-methoxy-4-[4-[4-(methylsulfonyl)-1-piperazinyl]-1-piperidinyl]phenyl]amino]-4-pyrimidinyl]imidazo[1,2-a]pyridin-2-yl]-2-methoxybenzamide

MF: C₄₄H₄₇F₂N₉O₅S

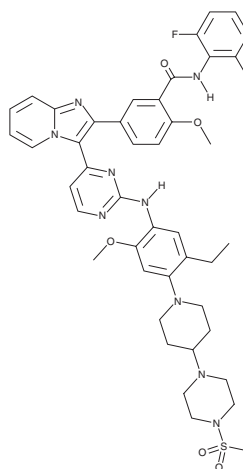
FW: 852.0

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.

Description

GSK1904529A is a dual inhibitor of the insulin receptor (InsR) and insulin-like growth factor 1 receptor (IGF-1R) kinases (IC₅₀s = 25 and 27 nM, respectively, in a cell-free assay).¹ It selectively inhibits InsR and IGF-1R kinases over a panel of 45 serine/threonine and tyrosine kinases (IC₅₀s = >1 μM for all). GSK1904529A decreases proliferation of 13 cancer cell lines, including multiple myeloma, sarcoma, colon, and breast cancer cells, with IC₅₀ values ranging from 35 to 189 nM in a panel of 35 cell lines. It reduces tumor growth in NIH3T3-LISN, COLO 205, HT-29, and BxPC3 mouse xenograft models when administered at a dose of 30 mg/kg for 21 days.

Reference

1. Sabbatini, P., Rowand, J.L., Groy, A., *et al.* Antitumor activity of GSK1904529A, a small-molecule inhibitor of the insulin-like growth factor-I receptor tyrosine kinase. *Clin. Cancer Res.* **15(9)**, 3058-3067 (2009).

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, 10/04/2022

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