

PRODUCT INFORMATION



GSK2983559

Item No. 31458

CAS Registry No.: 1579965-12-0

Formal Name: 2-[[4-(5-benzothiazolylamino)-6-[(1,1-dimethylethyl)sulfonyl]-7-quinazolinyl]oxy]-ethanol, 1-(dihydrogen phosphate)

MF: C₂₁H₂₃N₄O₇PS₂

FW: 538.5

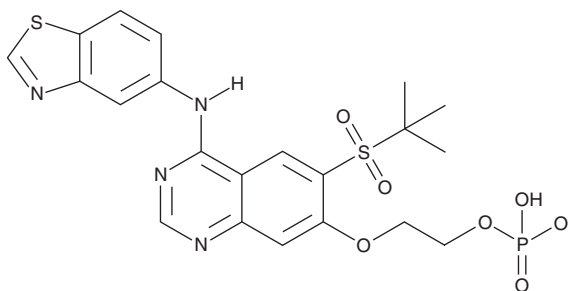
Purity: ≥98%

UV/Vis.: λ_{max}: 218, 245, 344 nm

Supplied as: A crystalline 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

GSK2983559 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK2983559 in the solvent of choice, which should be purged with an inert gas. GSK2983559 is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml. GSK2983559 is also slightly soluble in dimethyl formamide.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of GSK2983559 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of GSK2983559 in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

GSK2983559 is a receptor-interacting protein kinase 2 (RIPK2) inhibitor.¹ It inhibits RIPK2 by 65% in a kinase assay when used at a concentration of 10 μM. GSK2983559 (10 μM) also inhibits VEGFR3 by greater than 90%, as well as 14 additional kinases by 60 to 89% in a panel of 344 kinases. *In vivo*, GSK2983559 (7.5 and 145 mg/kg twice per day) reduces colonic damage in a mouse model of TNBS-induced colitis. GSK2983559 is also a prodrug that is cleaved to an active metabolite in the gastrointestinal tract that inhibits RIPK2 more potently than GSK2983559.

Reference

1. Haile, P.A., Casillas, L.N., Votta, B.J., *et al.* Discovery of a first-in-class receptor interacting protein 2 (RIP2) kinase specific clinical candidate, 2-((4-(benzo[d]thiazol-5-ylamino)-6-(*tert*-butylsulfonyl)quinazolin-7-yl)oxy)ethyl dihydrogen phosphate, for the treatment of inflammatory diseases. *J. Med. Chem.* **62**(14), 6482-6494 (2019).

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, 11/09/2022

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