

PRODUCT INFORMATION



GSK682753A

Item No. 45381

CAS Registry No.: 1334294-76-6
Formal Name: 8-[(2E)-3-(4-chlorophenyl)-1-oxo-2-propen-1-yl]-3-[(3,4-dichlorophenyl)methyl]-1-oxa-3,8-diazaspiro[4.5]decan-2-one

MF: C₂₃H₂₁Cl₃N₂O₃

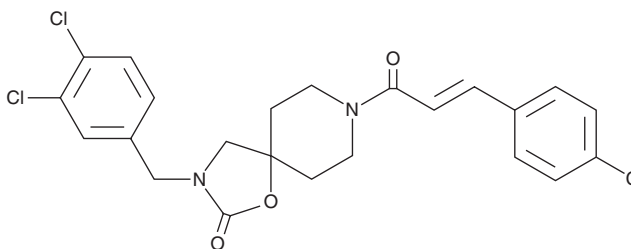
FW: 479.8

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

GSK682753A is supplied as a solid. A stock solution may be made by dissolving the GSK682753A in the solvent of choice, which should be purged with an inert gas. GSK682753A is slightly soluble (0.1-1 mg/ml) in methanol.

Description

GSK682753A is an Epstein-Barr virus-induced G protein-coupled receptor 2 (EBI2) inverse agonist.¹ It inhibits CREB activity in a reporter assay using HEK293 cells transiently transfected with EBI2 and Gqi4myr, a chimeric G_α subunit that recognizes G_{α_i} receptors but activates G_{α_q}-mediated pathways (IC₅₀ = 53.6 nM). GSK682753A (0.001-1 μM) inhibits antibody-induced proliferation of B cells isolated from transgenic mice overexpressing human EBI2.

Reference

1. Benned-Jensen, T., Smethurst, C., Holst, P. J., *et al.* Ligand modulation of the Epstein-Barr virus-induced seven-transmembrane receptor EBI2: Identification of a potent and efficacious inverse agonist. *The Journal of Biological Chemistry* **286**(33), 29292-29302 (2011).

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, 03/18/2026

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