

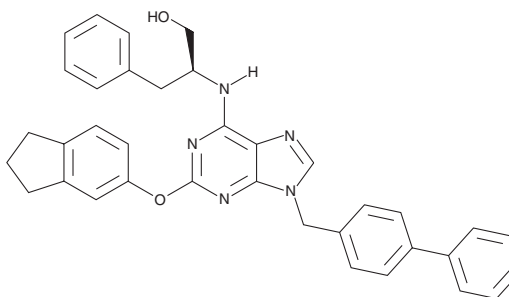
PRODUCT INFORMATION



QS11

Item No. 21247

CAS Registry No.: 944328-88-5
Formal Name: β S-[[9-([1,1'-biphenyl]-4-ylmethyl)-2-[(2,3-dihydro-1H-inden-5-yl)oxy]-9H-purin-6-yl]amino]-benzenepropanol
MF: C₃₆H₃₃N₅O₂
FW: 567.7
Purity: \geq 98%
UV/Vis.: λ_{max} : 272 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

QS11 is supplied as a crystalline solid. A stock solution may be made by dissolving the QS11 in the solvent of choice, which should be purged with an inert gas. QS11 is soluble in organic solvents such as ethanol and DMSO. The solubility of QS11 in these solvents is approximately 100 mM.

Description

QS11 is an inhibitor of GTPase activating protein of ADP-ribosylation factors (ARFGAPs) that increases activity of ARFs in NIH3T3 cells at a concentration of 1 μ M.¹ It activates Wnt/ β -catenin activity in a reporter assay using HEK293 cells (EC_{50} = 0.5 μ M). QS11 exhibits synergy with Wnt-3a ligand-conditioned media, inducing a 200-fold increase in reporter activity compared to 40- and 2-fold increases observed with Wnt-3a ligand-conditioned media alone or QS11 treatment alone, respectively. It inhibits ARFGAP1 with an EC_{50} value of 1.5 μ M in an enzyme assay and inhibits the migration of MDA-MB-231 breast cancer cells that overexpress AMAP1, an ARFGAP essential to cellular invasiveness.^{1,2}

References

1. Zhang, Q., Major, M.B., Takanashi, S., *et al.* Small-molecule synergist of the Wnt/ β -catenin signaling pathway. *Proc. Natl. Acad. Sci. U.S.A.* **104**(18), 7444-7448 (2007).
2. Singh, M.K., Gao, H., Sun, W., *et al.* Structure-activity relationship studies of QS11, a small molecule Wnt synergistic agonist. *Bioorg. Med. Chem. Lett.* **25**(21), 4838-4842 (2015).

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, 12/15/2022

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