

# PRODUCT INFORMATION



## S-Geranylgeranyl-L-glutathione

Item No. 9004116

CAS Registry No.: 2364639-42-7

Synonyms: GGG

MF:  $C_{30}H_{49}N_3O_6S$

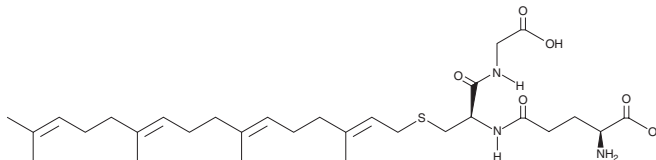
FW: 579.8

Purity:  $\geq 98\%$

Supplied as: A solid

Storage:  $-20^{\circ}C$

Stability:  $\geq 4$  years



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

### Laboratory Procedures

S-Geranylgeranyl-L-glutathione is supplied as a solid. A stock solution may be made by dissolving the S-geranylgeranyl-L-glutathione in the solvent of choice, which should be purged with an inert gas. S-Geranylgeranyl-L-glutathione is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml (warmed and vortexed).

### Description

S-Geranylgeranyl-L-glutathione is a ligand of the orphan G protein-coupled receptor (GPCR) P2RY8.<sup>1</sup> It selectively induces internalization of P2RY8 over sphingosine-1-phosphate receptor 2 (S1P<sub>2</sub>), GPR55, cysteinyl leukotriene receptor 1 (CysLT<sub>1</sub> receptor), and CysLT<sub>2</sub> receptor at 100 nM. S-Geranylgeranyl-L-glutathione (10 nM) inhibits the migration of P2RY8-expressing WEHI-231 B cell lymphoma cells and isolated human tonsil germinal center B cells induced by chemokine (C-X-C motif) ligand 12 (WEHI-231).

### Reference

1. Lu, E., Wolfreys, F.D., Muppidi, J.R., *et al.* S-Geranylgeranyl-L-glutathione is a ligand for human B cell-confinement receptor P2RY8. *Nature* **567(7724)**, 244-248 (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, 10/17/2023

#### 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