

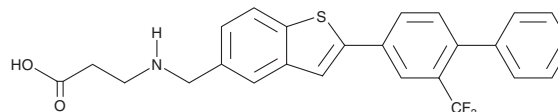
PRODUCT INFORMATION



AUY954

Item No. 9000548

CAS Registry No.: 820240-77-5
Formal Name: N-[[2-[2-(trifluoromethyl)[1,1'-biphenyl]-4-yl]benzo[b]thien-5-yl]methyl]-β-alanine
MF: C₂₅H₂₀F₃NO₂S
FW: 455.5
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 263, 311 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

AUY954 is supplied as a crystalline solid. A stock solution may be made by dissolving the AUY954 in the solvent of choice, which should be purged with an inert gas. AUY954 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AUY954 in these solvents is approximately 30 mg/ml.

AUY954 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AUY954 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AUY954 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

AUY954 is an orally bioavailable and selective agonist of the sphingosine-1-phosphate receptor 1 (S1P₁; EC₅₀ = 1.2 nM for stimulating GTPγS binding to S1P₁ in CHO cells).¹ It is highly selective for the S1P₁ receptor with EC₅₀ values of >10,000, 1,210, >1,000, and 340 nM for S1P₂, S1P₃, S1P₄, and S1P₅ receptors, respectively. AUY954 reduces circulating lymphocytes and prevents rejection of rat heart allografts. In a rat model of experimental autoimmune neuritis, chronic treatment with AUY954 (10 mg/kg), starting on the day of immunization, almost completely prevents peripheral paralysis.² In the same model, it also decreases T cell, B cell, and macrophage infiltration and perivascular demyelination in sciatic nerves.

References

1. Pan, S., Mi, Y., Pally, C., *et al.* A monoselective sphingosine-1-phosphate receptor-1 agonist prevents allograft rejection in a stringent rat heart transplantation model. *Chem. Biol.* **13(11)**, 1227-1234 (2006).
2. Zhang, Z.Y., Zhang, Z., Zug, C., *et al.* AUY954, a selective S1P₁ modulator, prevents experimental autoimmune neuritis. *J. Neuroimmunol.* **216(1-2)**, 59-65 (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, 12/12/2022

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