

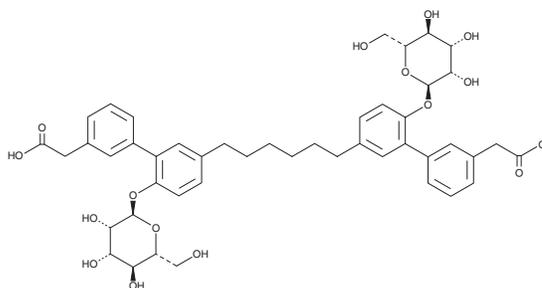
PRODUCT INFORMATION



Bimosiamose

Item No. 30878

CAS Registry No.: 187269-40-5
Formal Name: 3',3'''-(1,6-hexanediyl)bis[6'-(α -D-mannopyranosyloxy)-[1,1'-biphenyl]-3-acetic acid
Synonym: TBC 1269
MF: C₄₆H₅₄O₁₆
FW: 862.9
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

Bimosiamose is supplied as a crystalline solid. A stock solution may be made by dissolving the bimosiamose in the solvent of choice, which should be purged with an inert gas. Bimosiamose is soluble in DMSO.

Description

Bimosiamose is a pan-selectin inhibitor (IC_{50} s = 88, 20, and 86 μ M for E-, P-, and L-selectin, respectively).¹ It reduces adhesion of HL-60 cells to E- and P-selectin when used at a concentration of 129 μ M.² *In vivo*, bimosiamose (200 mg/kg) reduces skin scaliness, induration, and erythema in a patient-derived xenograft mouse model of psoriasis. It decreases hepatic alanine aminotransferase (ALT), aspartate aminotransferase (AST), lactate dehydrogenase (LDH), and myeloperoxidase (MPO) levels and reduces neutrophil infiltration in a rat model of severe hepatic ischemia-reperfusion injury.¹ Bimosiamose also prevents nematode-induced late airway reactivity in sheep and reduces airway hyperresponsiveness in a mouse model of asthma.³

References

1. Palma-Vargas, J.M., Toledo-Pereyra, L., Dean, R.E., *et al.* Small-molecule selectin inhibitor protects against liver inflammatory response after ischemia and reperfusion *J. Am. Coll. Surg.* **185**(4), 365-372 (1997).
2. Friedrich, M., Bock, D., Philipp, S., *et al.* Pan-selectin antagonism improves psoriasis manifestation in mice and man. *Arch. Dermatol. Res.* **297**(8), 345-351 (2008).
3. Aydt, E. and Wolff, G. Development of synthetic pan-selectin antagonists: A new treatment strategy for chronic inflammation in asthma. *Pathobiology* **70**(5), 297-301 (2002).

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, 06/17/2020

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