

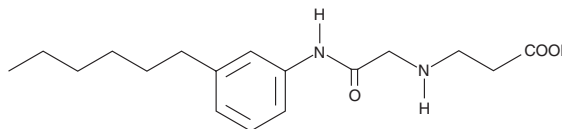
PRODUCT INFORMATION



W123

Item No. 10010992

CAS Registry No.: 1345982-24-2
Formal Name: 3-(2-(3-hexylphenylamino)-2-oxoethylamino)propanoic acid
MF: C₁₇H₂₆N₂O₃
FW: 306.4
Purity: ≥98%
UV/Vis.: λ_{max}: 244 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

W123 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of W123 in these solvents is approximately 5 mg/ml.

W123 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of W123 should be diluted with the aqueous buffer of choice. W123 has a solubility of approximately 0.3 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method.

Description

Sphingosine-1-phosphate (S1P) modulates skin, vascular endothelial, and immune cells via five distinct G protein-coupled receptors. The S1P type 1 receptor, S1P₁, also known as EDG-1, mediates endothelial, lymphoid, and neuronal cell responses to S1P, as well as the synthetic agonist FTY720. W123 is an analog of FTY720 that is a competitive antagonist of S1P₁, as measured by GTPγS activation, mitogen-activated protein kinase recruitment, cell migration, and ligand-induced receptor internalization.¹ It has a K_i value of 0.69 μM when assessed by SEW2871-induced activation of S1P₁.¹ Its specificity for other S1P receptor subtypes remains to be determined.

Reference

1. Wei, S.H., Rosen, H., Matheu, M.P., *et al.* Sphingosine-1-phosphate type 1 receptor agonism inhibits transendothelial migration of medullary T cells to lymphatic sinuses. *Nature Immunology* **6**(12), 1228-1235 (2005).

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, 02/28/2019

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