

PRODUCT INFORMATION

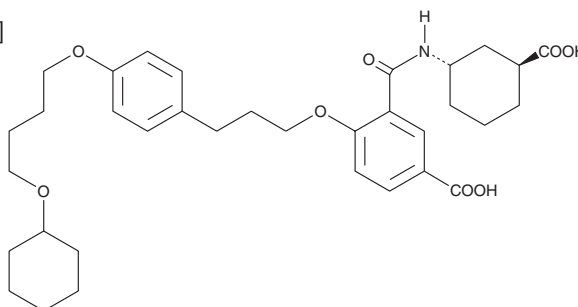


HAMI3379

Item No. 10580

CAS Registry No.: 1245653-57-9
Formal Name: 3-[[[(1S,3S)-3-carboxycyclohexyl]amino]carbonyl]-4-[3-[4-(4-(cyclohexyloxy)butoxy)phenyl]propoxy]-benzoic acid

MF: C₃₄H₄₅NO₈
FW: 595.7
Purity: ≥90%
UV/Vis.: λ_{max}: 220 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

HAMI3379 is supplied as a crystalline solid. A stock solution may be made by dissolving the HAMI3379 in the solvent of choice, which should be purged with an inert gas. HAMI3379 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of HAMI3379 in ethanol is approximately 5 mg/ml and approximately 20 mg/ml in DMSO and DMF.

HAMI3379 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, HAMI3379 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. HAMI3379 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

HAMI3379 is a cysteinyl leukotriene 2 (CysLT₂) receptor antagonist (IC₅₀ = 37.9 nM in a radioligand binding assay).¹ It is selective for CysLT₂ over CysLT₁ (IC₅₀ = >30 μM in a radioligand binding assay). HAMI3379 inhibits calcium mobilization induced by leukotriene D₄ (LTD₄; Item No. 20310) or leukotriene C₄ (LTC₄; Item No. 20210) in CHO cells expressing human CysLT₂ (IC₅₀s = 3.8 and 4.4 nM, respectively). It reverses the LTC₄-induced increases in perfusion pressure and decreases in contractility in isolated Langendorff-perfused guinea pig hearts in a concentration-dependent manner. HAMI3379 (0.1 mg/kg) prevents decreases in body weight and neurological deficit scores in a rat model of *S. pneumoniae*-induced meningitis, as well as reduces infarct volume in a rat model of cerebral ischemia brain injury induced by middle cerebral artery occlusion (MCAO).^{2,3} It also prevents increases in airway hyperresponsiveness in a mouse model of ovalbumin-induced asthma when administered at a dose of 10 mg/kg.⁴

References

1. Wunder, F., Tinel, H., Kast, R., *et al.* Pharmacological characterization of the first potent and selective antagonist at the cysteinyl leukotriene 2 (CysLT₂) receptor. *Br. J. Pharmacol.* **160**(2), 399-409 (2010).
2. Yu, S., Chen, X., Li, X., *et al.* Neuroprotective effects of CysLTR antagonist on *Streptococcus pneumoniae*-induced meningitis in rats. *Exp. Ther. Med.* **24**(1), 443 (2022).
3. Shi, Q.J., Wang, H., Liu, Z.X., *et al.* HAMI 3379, a CysLT₂R antagonist, dose- and time-dependently attenuates brain injury and inhibits microglial inflammation after focal cerebral ischemia in rats. *Neuroscience* **291**, 53-69 (2015).
4. Trinh, H.K.T., Suh, D.-H., Nguyen, T.V.T., *et al.* Characterization of cysteinyl leukotriene-related receptors and their interactions in a mouse model of asthma. *Prostaglandins Leukot. Essent. Fatty Acids* **141**, 17-23 (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, 02/06/2023

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