

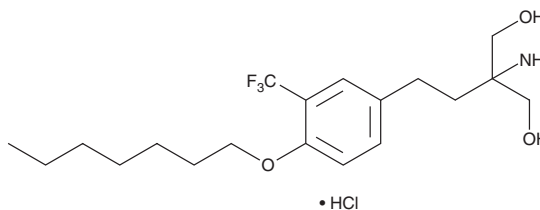
PRODUCT INFORMATION



Amiselimod (hydrochloride)

Item No. 20970

CAS Registry No.: 942398-84-7
Formal Name: 2-amino-2-[2-[4-(heptyloxy)-3-(trifluoromethyl)phenyl]ethyl]-1,3-propanediol, monohydrochloride
Synonym: MT-1303
MF: C₁₉H₃₀F₃NO₃ • HCl
FW: 413.9
Purity: ≥95%
UV/Vis.: λ_{max}: 225, 283 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

Amiselimod (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the amiselimod (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Amiselimod (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of amiselimod (hydrochloride) in these solvents is approximately 15, 20, and 25 mg/ml, respectively.

Amiselimod (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, amiselimod (hydrochloride) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Amiselimod (hydrochloride) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Amiselimod is a prodrug form of the sphingosine-1-phosphate receptor 1 (S1P₁) antagonist amiselimod phosphate (amiselimod-P).¹ *In vivo*, amiselimod (0.3 mg/kg per day) inhibits the development of chronic colitis induced by adoptive transfer of CD4⁺CD45RB^{high} T cells in mice.²

References

1. Sugahara, K., Maeda, Y., Shimano, K., *et al.* Amiselimod, a novel sphingosine 1-phosphate receptor-1 modulator, has potent therapeutic efficacy for autoimmune diseases, with low bradycardia risk. *Br. J. Pharmacol.* **174(1)**, 15-27 (2017).
2. Shimano, K., Maeda, Y., Kataoka, H., *et al.* Amiselimod (MT-1303), a novel sphingosine 1-phosphate receptor-1 functional antagonist, inhibits progress of chronic colitis induced by transfer of CD4⁺CD45RB^{high} T cells. *PLoS One* **14(12)**, e0226154 (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, 12/14/2022

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