

PRODUCT INFORMATION



Dihydro Montelukast

Item No. 22374

CAS Registry No.: 142147-67-9
Formal Name: 1-[[[(1R)-1-[3-[2-(7-chloro-2-quinolinyl)ethyl]phenyl]-3-[2-(1-hydroxy-1-methylethyl)phenyl]propyl]thio]methyl]-cyclopropaneacetic acid

MF: C₃₅H₃₈ClNO₃S

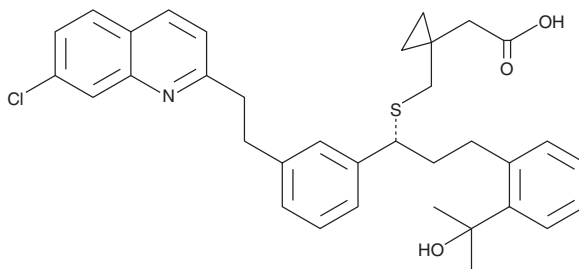
FW: 588.2

Purity: ≥98%

Supplied as: A 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

Dihydro montelukast is supplied as a solid. A stock solution may be made by dissolving the dihydro montelukast in the solvent of choice, which should be purged with an inert gas. Dihydro montelukast is soluble in chloroform.

Description

Dihydro montelukast is a potential impurity found in commercial montelukast preparations. Montelukast (sodium salt) (Item No. 10008318) is a potent and selective cysteinyl leukotriene 1 (CysLT₁) receptor antagonist. Formulations containing montelukast are used for the treatment of asthma as well as for the symptoms associated with allergic rhinitis.¹⁻⁴

References

1. Leff, J. A., Busse, W. W., Pearlman, D., *et al.* Montelukast, a leukotriene-receptor antagonist, for the treatment of mild asthma and exercise-induced bronchoconstriction. *N. Engl. J. Med.* **339(3)**, 147-152 (1998).
2. Reiss, T. F., Chervinsky, P., Dockhorn, R. J., *et al.* Montelukast, a once-daily leukotriene receptor antagonist, in the treatment of chronic asthma. A multicenter, randomized, double-blind trial. *Arch. Intern. Med.* **158(11)**, 1213-1220 (1998).
3. Reiss, T. F., Altman, L. C., Chervinsky, P., *et al.* Effects of montelukast (MK-0476), a new potent cysteinyl leukotriene (LTD₄) receptor antagonist, in patients with chronic asthma. *J. Allergy Clin. Immunol.* **98(3)**, 528-534 (1996).
4. Peters-Golden, M. and Henderson, W.R. The role of leukotrienes in allergic rhinitis. *Ann. Allergy Asthma Immunol.* **94(6)**, 609-618 (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, 11/14/2022

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