

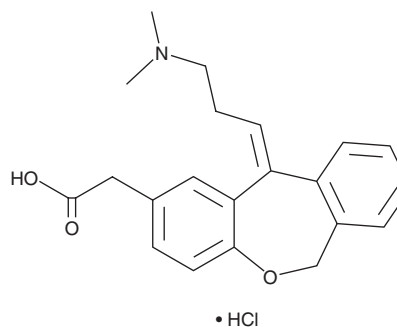
# PRODUCT INFORMATION



## Olopatadine (hydrochloride)

Item No. 11999

**CAS Registry No.:** 140462-76-6  
**Formal Name:** 11Z-[3-(dimethylamino)propylidene]-6,11-dihydro-dibenz[b,e]oxepin-2-acetic acid, monohydrochloride  
**Synonyms:** ALO 4943A, KW 4679  
**MF:**  $C_{21}H_{23}NO_3 \cdot HCl$   
**FW:** 373.9  
**Purity:**  $\geq 98\%$   
**UV/Vis.:**  $\lambda_{max}$ : 206, 301 nm  
**Supplied as:** A crystalline solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

Olopatadine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the olopatadine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Olopatadine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of olopatadine (hydrochloride) in these solvents is approximately 0.25, 3, and 5 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of olopatadine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of olopatadine (hydrochloride) in PBS (pH 7.2) is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Olopatadine is a histamine  $H_1$  receptor antagonist ( $K_i = 41$  nM).<sup>1</sup> It is 1,059- and 4,177-fold selective for histamine  $H_1$  over  $H_2$  and  $H_3$  receptors, respectively. Olopatadine inhibits histamine-induced phosphoinositide turnover in isolated human conjunctival epithelial cells, isolated human corneal fibroblasts, and human trabecular meshwork (TM3) cells ( $IC_{50}$ s = 9.5, 19, and 39.9 nM, respectively). *In vivo*, olopatadine inhibits passive cutaneous anaphylaxis in rats ( $ED_{50}$  = 49  $\mu$ g/kg) and IgG1-mediated bronchoconstriction in ovalbumin-sensitized guinea pigs ( $ED_{50}$  = 30  $\mu$ g/kg).<sup>2</sup> Formulations containing olopatadine have been used in the treatment of allergic rhinitis and conjunctivitis, as well as in the treatment of itch in patients with well-controlled urticaria.

### References

1. Sharif, N.A., Xu, S.X., and Yanni, J.M. Olopatadine (AL-4943A): Ligand binding and functional studies on a novel, long acting  $H_1$ -selective histamine antagonist and anti-allergic agent for use in allergic conjunctivitis. *J. Ocul. Pharmacol. Ther.* **12**(4), 401-407 (1996).
2. Ohshima, E., Otaki, S., Sato, H., et al. Synthesis and antiallergic activity of 11-(aminoalkylidene)-6,11-dihydrodibenz[b,e]oxepin derivatives. *J. Med. Chem.* **35**(11), 2074-2084 (1992).

#### 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, 11/08/2022

#### 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