

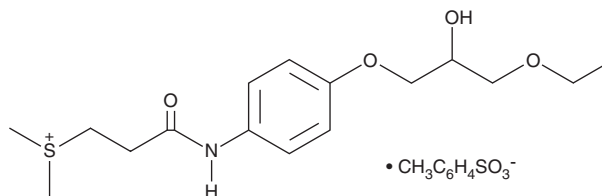
PRODUCT INFORMATION



Suplatast (tosylate)

Item No. 30316

CAS Registry No.: 94055-76-2
Formal Name: [3-[[4-(3-ethoxy-2-hydroxypropoxy)phenyl]amino]-3-oxopropyl]dimethyl-sulfonium, 4-methylbenzenesulfonate
MF: C₁₆H₂₆NO₄S • C₇H₇O₃S
FW: 499.6
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 253 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

Suplatast (tosylate) is supplied as a crystalline solid. A stock solution may be made by dissolving the suplatast (tosylate) in the solvent of choice, which should be purged with an inert gas. Suplatast (tosylate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of suplatast (tosylate) in these solvents is approximately 30 mg/ml.

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 suplatast (tosylate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of suplatast (tosylate) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Suplatast is an antiallergic agent.^{1,2} It inhibits IL-4 and IL-5 production in conalbumin-stimulated D10.G4.1 murine T helper 2 (Th2) cells in a concentration-dependent manner.¹ Suplatast (100 mg/kg) inhibits ovalbumin-induced increases in eosinophil and total cell numbers, as well as IL-4, IL-5, and IL-13, but not IFN-γ, levels in bronchoalveolar lavage fluid (BALF) in an ovalbumin-sensitized mouse model of asthma.² It also inhibits ovalbumin-induced increases in ovalbumin-specific IgE in serum and bronchial hyperresponsiveness to methacholine in the ovalbumin-sensitized mouse model of asthma.

References

1. Yamaya, H., Basaki, Y., Togawa, M., *et al.* Down-regulation of Th2 cell-mediated murine peritoneal eosinophilia by antiallergic agents. *Life Sci.* **56(19)**, 1647-1654 (1995).
2. Zhao, G.D., Yokoyama, A., Kohno, N., *et al.* Effect of suplatast tosilate (IPD-1151T) on a mouse model of asthma: Inhibition of eosinophilic inflammation and bronchial hyperresponsiveness. *Int. Arch. Allergy. Immunol.* **121(2)**, 116-122 (2000).

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, 05/20/2020

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