

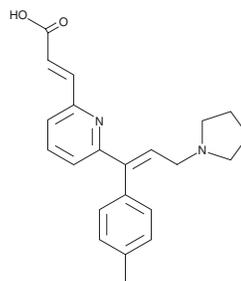
# PRODUCT INFORMATION



## Acrivastine

Item No. 24021

**CAS Registry No.:** 87848-99-5  
**Formal Name:** (2E)-3-[6-[(1E)-1-(4-methylphenyl)-3-(1-pyrrolidinyl)-1-propen-1-yl]-2-pyridinyl]-2-propenoic acid  
**Synonyms:** BW 0270C, BW A825C  
**MF:** C<sub>22</sub>H<sub>24</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 348.4  
**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

Acrivastine is supplied as a solid. A stock solution may be made by dissolving the acrivastine in the solvent of choice, which should be purged with an inert gas. Acrivastine is slightly soluble in DMSO and methanol.

### Description

Acrivastine is a histamine H<sub>1</sub> receptor antagonist with a K<sub>i</sub> value of 10 nM in COS-7 cells expressing the human receptor.<sup>1</sup> *In vivo*, acrivastine (1 mg/kg) completely inhibits response to histamine in guinea pigs.<sup>2</sup> Formulations containing acrivastine have been used for the treatment of seasonal allergies and hay fever.

### References

1. Wieland, K., Laak, A.M., Smit, M.J., *et al.* Mutational analysis of the antagonist-binding site of the histamine H<sub>1</sub> receptor. *J. Biol. Chem.* **274**(42), 29994-30000 (1999).
2. Hoshiko, K., Chapman, I.D., and Morley, J. Histamine(H1) antagonists and airway hyperreactivity in the guinea-pig. *Agents Actions Suppl.* **34**, 323-333 (1991).

#### 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, 11/21/2022

#### 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