

# Ketotifen . fumarate

Histamine receptor ligand

Selective histamine H<sub>1</sub> antagonist. Inhibits histamine release from monocytes and lymphocytes at 0.1-10 µM.

Citations: 3

[View Online »](#)

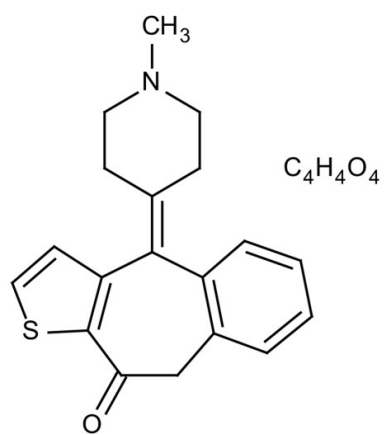
Ordering Information

[Order Online »](#)

|                |    |
|----------------|----|
| BML-NH106-0001 | 1g |
|----------------|----|

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** Ambient

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Appearance** White solid.

**CAS** 34580-14-8

**Couple Target** Histamine receptor

**Couple Type** Ligand

**Formula**  $C_{19}H_{19}NOS \cdot C_4H_4O_4$

**MW** 425.5

**Purity**  $\geq 98\%$  (HPLC)

**Solubility** Soluble in water (10 mg/ml).

**Last modified:** May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)