

# Histamine . dihydrochloride

Histamine triggers the inflammatory response.

Citations: 7

[View Online »](#)

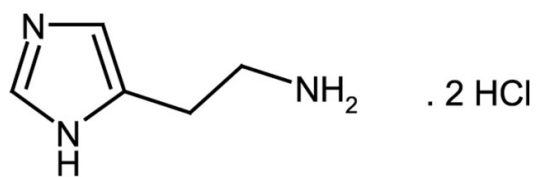
## Ordering Information

[Order Online »](#)

|                  |    |
|------------------|----|
| ALX-550-132-G005 | 5g |
|------------------|----|

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Long Term Storage** +4°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 1H-Imidazole-4-ethanamine . 2HCl

**Appearance** White to off-white solid.

**CAS** 56-92-8

**Couple Target** Histamine receptor

**Couple Type** Activator, Ligand

**Formula**  $C_5H_9N_3 \cdot 2HCl$

**Identity** Determined by IR.

**MI** **14:** 4719

**MW** 111.1 . 73.0

**Purity** ≥98% (HPLC)

**RTECS** MS1575000

**Solubility** Soluble in water (50mg/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)