

Histamine . dihydrochloride

Histamine triggers the inflammatory response.

Citations: 7

[View Online »](#)

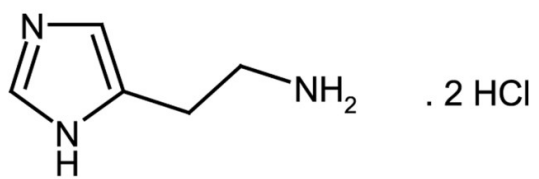
Ordering Information

[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)