



# Ultra-Sensitive ELISA Kit Detects Histamine in Both Native and Converted States

## Histamine ELISA Kit

**Ultra-sensitive Histamine ELISA kit enabling detection of histamine in serum, plasma, urine and tissue culture media**

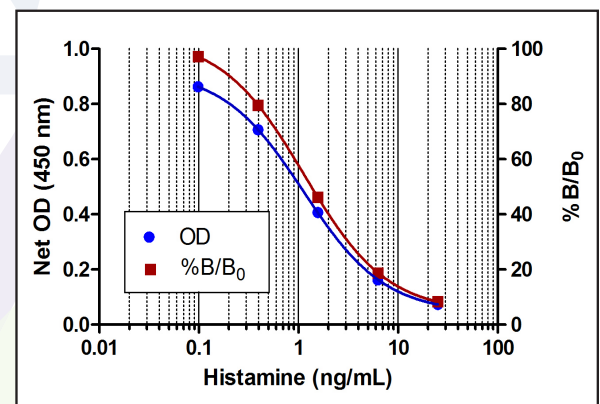
### Histamine ELISA Kit (ENZ-KIT140)

Histamine is an important mediator of immediate-type-allergic reactions. Because of its potent role in the immune response, histamine is present in a very transitory state in the body and can only be measured within minutes of release. The half-life of histamine in a biological system is four minutes before its conversion to n-methyl histamine. N-methyl histamine is a major product of histamine metabolism, is present in urine and has a longer half-life than that of histamine. This assay reacts equally to n-methyl histamine and histamine, making it an excellent choice for measuring histamine levels in urine.

#### Key Features

- Ultra-sensitive measurement of histamine, detecting as little as 0.03 ng/ml
- Does not require an acetylation step
- N-methyl histamine cross-reacts 100% in the assay. Histamine is detected in serum and plasma and n-methyl histamine is detected in urine
- Broad dynamic range suitable for a large variety of samples
- High throughput format with results in 2 hours for up to 40 samples in duplicate

Ultra-sensitive Measurement of Histamine



Standard curve experiment using Histamine ELISA Kit (ENZ-KIT140)

PRODUCT NAME	PRODUCT #	PRODUCT NAME	PRODUCT #
Histamine ELISA Kit	ENZ-KIT140	IL-2 ELISA Kit	ADI-900-118A
Histamine pAb	ENZ-ABS292	IL-4 ELISA Kit	ADI-900-145A
ALDetect™ Lipid Peroxidation Assay Kit	BML-AK170	IL-5 ELISA Kit	ENZ-KIT139
Angiotensin I ELISA Kit	ADI-900-203	IL-6 ELISA Kit	ADI-900-033
Angiotensin II ELISA Kit	ADI-900-204	IL-10 ELISA Kit	ADI-900-036
Bradykinin ELISA Kit	ADI-900-206	LTB <sub>4</sub> ELISA Kit	ADI-900-068
COX-2 ELISA Kit	ADI-900-094	PGE <sub>2</sub> ELISA Kit	ADI-900-001
CRP ELISA Kit	ENZ-KIT102	Serotonin ELISA Kit	ADI-900-175
Endothelin-I ELISA Kit	ADI-900-020A	SOD Activity Kit	ADI-900-157
IL-1β ELISA Kit	ADI-900-130	TNFα ELISA Kit	ADI-900-099

Advance Your Research:  
[www.enzolifesciences.com](http://www.enzolifesciences.com)

#### GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.  
 10 Executive Boulevard  
 Farmingdale, NY 11735  
 Phone: 800.942.0430  
 Fax: 631.694.7501  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

#### EUROPE

Enzo Life Sciences (ELS) AG  
 Industriestrasse 17  
 CH-4415 Lausen, Switzerland  
 Phone: +41 61 926 8989  
 Fax: +41 61 926 8979  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)