

Lipid Peroxidation (4-HNE) Assay Kit ab238538

[33 References](#) [1 Image](#)

Overview

Product name	Lipid Peroxidation (4-HNE) Assay Kit
Detection method	Colorimetric
Sample type	Serum, Plasma, Cell Lysate, Purified protein, Tissue Lysate
Assay type	Competitive
Species reactivity	Reacts with: Vertebrata
Product overview	Lipid Peroxidation (4-HNE) Assay Kit (ab238538) is designed for the rapid detection and quantitation of 4-hydroxynonenal (4-HNE)-protein adducts.

4-HNE is a well known by-product of lipid peroxidation and is widely accepted as a stable marker for oxidative stress. This Competitive ELISA kit will allow the quantitation of 4-HNE adduct in protein samples determined by comparing its absorbance with that of a known 4-HNE-BSA standard curve.

Platform Microplate reader

Properties

Storage instructions Please refer to protocols.

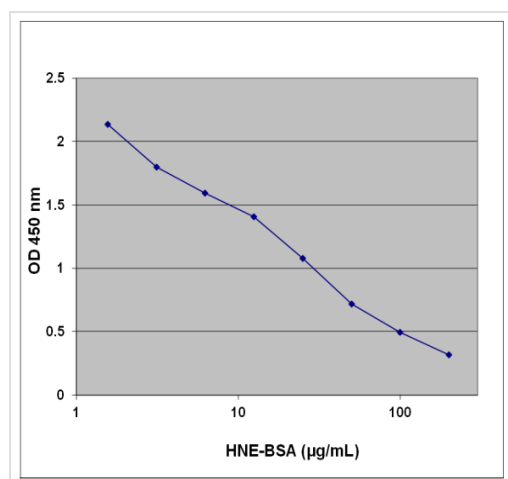
Components	96 tests	32 tests
100X Conjugate Diluent	1 x 300µl	1 x 100µl
10X Wash Buffer	1 x 100ml	1 x 30ml
Protein Binding Strip Well Plate	1 unit	1 unit
Anti-HNE Antibody (1000X)	1 x 10µl	1 x 5µl
Assay Diluent	1 x 50ml	1 x 20ml
HNE Conjugate	1 x 50µl	1 x 20µl
HNE-BSA Standard	1 x 250µl	1 x 100µl

Components	96 tests	32 tests
Secondary Antibody, HRP Conjugate (1000X)	1 x 20µl	1 x 20µl
Stop Solution	1 x 12ml	1 x 4ml
Substrate Solution	1 x 12ml	1 x 4ml

Relevance Aldehydic products of lipid peroxidation, such as 4 hydroxynonenal (4 HNE), have been implicated in the etiology of pathological changes under oxidative stress as a key mediator of oxidative stress induced cell death. It is a stable product of lipid peroxidation, is proarrhythmic and may contribute to the cytotoxic effects of oxidative stress.

Cellular localization Cytoplasmic

Images



HNE-BSA Competitive ELISA Standard Curve.

The following figures demonstrate typical HNE Adduct Competitive ELISA results. One should use the data below for reference only. This data should not be used to interpret actual results.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors