

Product datasheet

Hyaluronic Acid ELISA Kit ab287799

1 Image

Overview

Product name Hyaluronic Acid ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 8%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Sample type Serum, Plasma, Other biological fluids, Tissue Homogenate

Assay type Competitive

Sensitivity < 0.938 ng/ml

Range 1.563 ng/ml - 100 ng/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	95	85% - 103%
Hep Plasma	91	85% - 99%
EDTA Plasma	94	87% - 104%

Assay duration Multiple steps standard assay

Product overview

Hyaluronic acid (HA) is an anionic, nonsulfated glycosaminoglycan distributed widely throughout connective, epithelial, and neural tissues. It is unique among glycosaminoglycans in that it is nonsulfated, forms in the plasma membrane instead of the Golgi apparatus, and can be very large. One of the chief components of the extracellular matrix, hyaluronan contributes significantly to cell proliferation and migration, and may also be involved in the progression of some malignant tumors. The Hyaluronic Acid ELISA kit ab287799 is a competitive ELISA assay for the

quantitative measurement of Hyaluronic Acid in serum, plasma and cell culture supernatants. The density of color is proportional to the amount of Hyaluronic Acid captured from samples.

Platform Microplate (12 x 8 well strips)

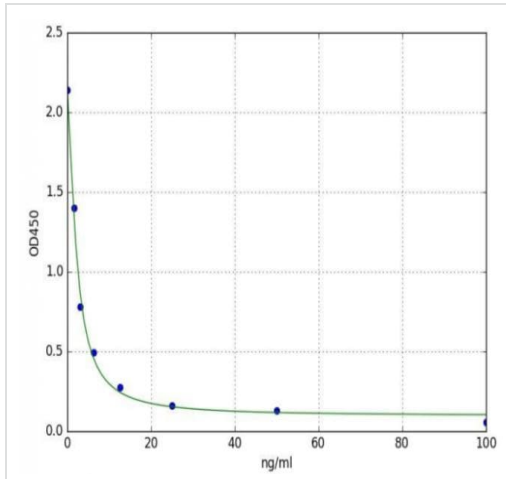
Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	96 tests
25X Wash buffer	1 x 30ml
Antibody dilution buffer	1 x 10ml
Biotin-Detection antibody (Concentrated)	1 x 60µl
HRP-Streptavidin Conjugate (SABC)	1 x 120µl
Lyophilized Standard	2 vials
Micro ELISA Plate	1 unit
Plate Sealer	5 units
SABC dilution buffer	1 x 10ml
Sample / Standard Dilution Buffer	1 x 20ml
Stop Solution	1 x 10ml
TMB Substrate	1 x 10ml

Relevance Hyaluronic acid is a mucopolysaccharide which is an integral part of the gel-like substance of animal connective tissue that serves as a lubricating and protective agent.

Images



Typical Standard Curve: These standard curves are for demonstration only. A standard curve must be run with each assay.

Hyaluronic Acid (HA) ELISA Kit Standard Curve

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