

Human Laminin ELISA Kit ab119599

3 References 2 Images

Overview

Product name	Human Laminin ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Cell Lysate
Assay type	Sandwich (quantitative)
Sensitivity	< 10 pg/ml
Range	156 pg/ml - 10000 pg/ml
Assay duration	Multiple steps standard assay
Species reactivity	<b>Reacts with:</b> Human
Product overview	<p>Human Laminin <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Human Laminin in cell culture supernatants, serum and plasma (heparin, EDTA).</p> <p>A Laminin specific rabbit polyclonal antibody has been precoated onto 96-well plates. Standards and test samples are added to the wells and incubated. A biotinylated detection polyclonal antibody from rabbit specific for Laminin is then added followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates were washed away with PBS or TBS buffer. TMB is then used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the Human Laminin amount of sample captured in plate.</p>
Platform	Microplate

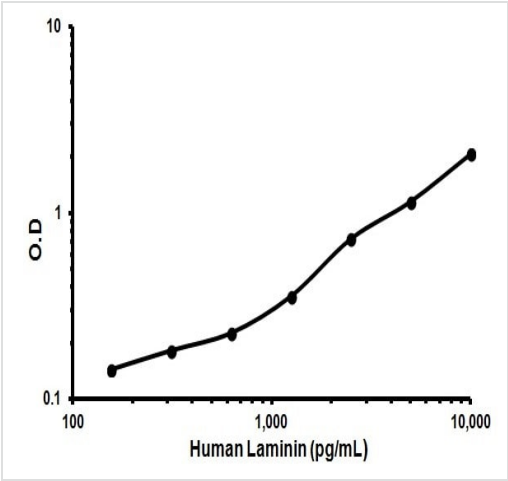
Properties

Storage instructionsStore at -20°C. Please refer to protocols.

Components	Identifier	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml

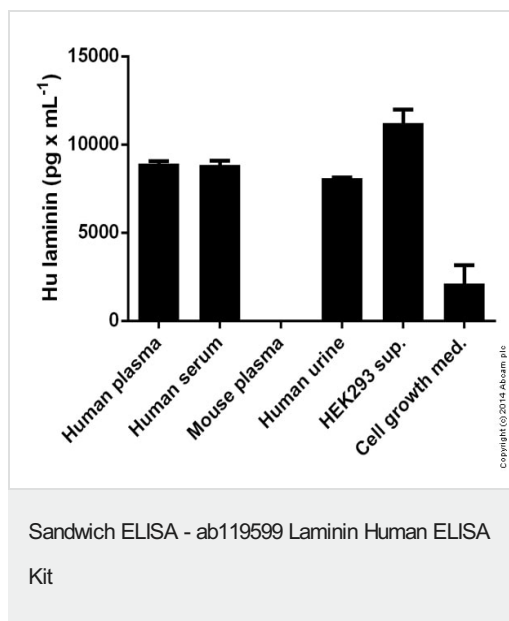
Components	Identifier	1 x 96 tests	1 x 96 tests
Anti-Human Laminin antibody Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti-Human Laminin antibody		1 x 100µl	1 x 100µl
Lyophilized Human Laminin standard		2 x 10ng	2 x 10ng
Plate Seal		1 x 4 units	1 x 4 units
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Amber Bottle	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

Images



Standard curve: mean of duplicates (+/-SD) with background readings subtracted

Sandwich ELISA - ab119599 Laminin Human ELISA Kit



Hu laminin measured in various biological fluids showing quantity (pg) per mL of tested sample

**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