

Rat Laminin ELISA Kit ab119573

画像数 1

製品の概要

製品名	Rat Laminin ELISA Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma, Other biological fluids, Tissue Extracts
アッセイタイプ	Sandwich (quantitative)
検出感度	< 10 pg/ml
検出範囲	156 pg/ml - 10000 pg/ml
ステップ	Multiple steps standard assay
種交差性	交差種: Rat
製品の概要	Abcam's rat Laminin <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of rat Laminin in cell culture supernatants, serum, cell lysates and plasma (heparin, EDTA).

A Laminin specific mouse monoclonal antibody has been precoated onto 96-well plates. Standards (form rat sarcoma basement membrane) and test samples are added to the wells and incubated. A biotinylated detection polyclonal antibody from goat, specific for Laminin is then added followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are 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 rat Laminin amount of sample captured in plate.

試験プラットフォーム	Microplate
------------	------------

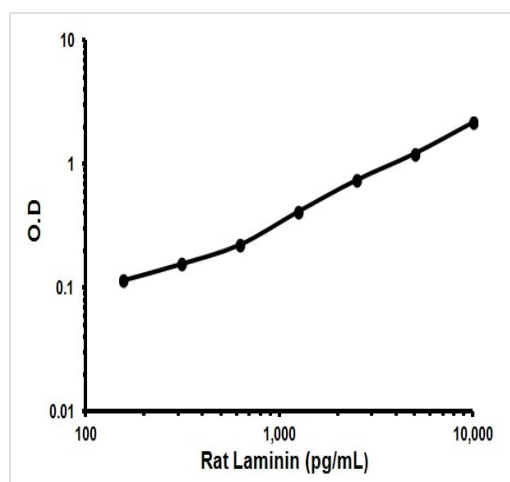
製品の特性

保存方法	Store at -20°C. Please refer to protocols.
------	--

内容	ラベル	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml

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

画像



Representative Standard Curve using ab119573

Typical 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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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