

Mouse LH Beta ELISA Kit ab235648

リコンビナント SimpleStep ELISA

★★★★★ 1 Abreviews 1 References 画像数 4

製品の概要

製品名	Mouse LH Beta ELISA Kit				
検出方法	Colorimetric				
再現性	Intra-Assay (同時再現性)				
	サンプル	N	平均値	SD	CV%
	Pituitary	8			4.9%
サンプルの種類	Tissue Extracts				
アッセイタイプ	Sandwich (quantitative)				
検出感度	4.7 pg/ml				
検出範囲	117 pg/ml - 7500 pg/ml				
添加回収試験	特定サンプルでの回収試験				
	サンプルの種類		平均 %	測定範囲	
	Tissue Extracts		97	% - %	
全工程の試験時間	1h 30m				
ステップ	One step assay				
種交差性	交差種: Mouse				
製品の概要	Mouse LH Beta ELISA Kit (ab235648) is a single-wash 90 min sandwich ELISA designed for the				

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate ([ab203359](#)) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

## 特記事項

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

## 試験プラットフォーム

Pre-coated microplate (12 x 8 well strips)

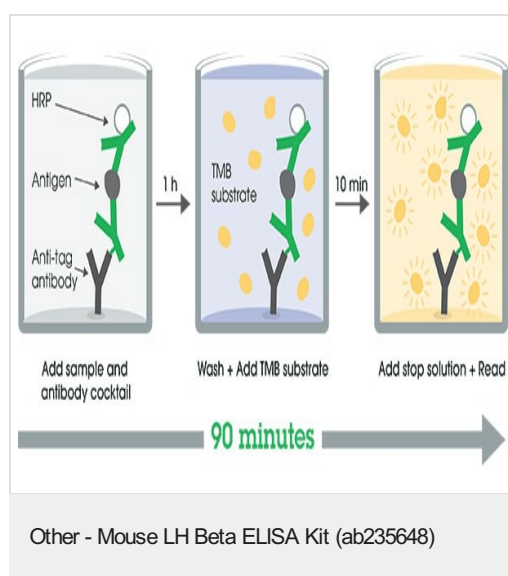
## 製品の特性

### 保存方法

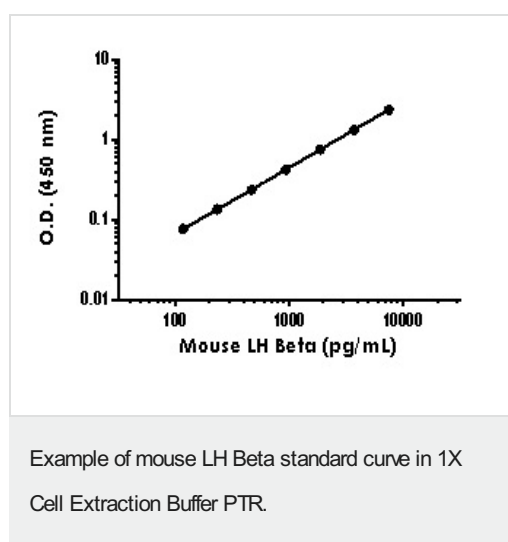
Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
10X Mouse LH Beta Capture Antibody	1 x 600µl
10X Mouse LH Beta Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
50X Cell Extraction Enhancer Solution (ab193971)	1 x 1ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml
Antibody Diluent 5BI	1 x 6ml
Mouse LH Beta Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 12ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

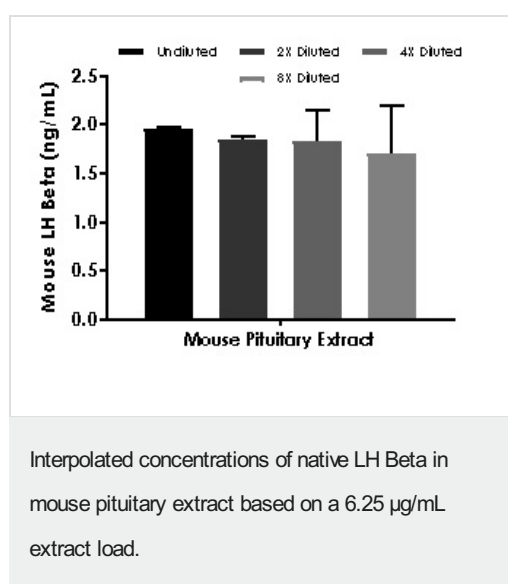
## 画像



SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



Background-subtracted data values (mean  $\pm$  SD) are graphed.



The concentrations of LH Beta were measured in duplicate and interpolated from the LH Beta standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean  $\pm$  SD,  $n=2$ ). The mean LH Beta concentration was determined to be 1.84 ng/mL in mouse pituitary extract.

Powered by  
recombinant antibodies



Research with  
confidence  
Consistent and  
reproducible results



Long-term and  
scalable supply  
Recombinant  
technology



Success from the  
first experiment  
Confirmed  
specificity



Ethical standards  
compliant  
Animal-free  
production

Sandwich ELISA - Mouse LH Beta ELISA Kit  
(ab235648)

To learn more about the advantages of recombinant antibodies see [here](#).

**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