

Mouse mTOR ELISA Kit ab206311

リコンビナント SimpleStep ELISA

★★★★★ 1 Abreviews 画像数 6

製品の概要

製品名	Mouse mTOR 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%
	cell lysate	8			= 3.1%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	cell lysate	3			= 6.9%
サンプルの種類	Cell culture supernatant, Cell culture extracts, Tissue Extracts, Tissue Homogenate				
アッセイタイプ	Sandwich (quantitative)				
検出感度	16.7 pg/ml				
検出範囲	39.06 pg/ml - 2500 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	
	Cell culture media	97		94% - 101%	
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 mTOR ELISA Kit (ab206311) 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

Mammalian target of rapamycin (mTOR) is a serine/threonine protein kinase part of two distinct signaling complexes, mTORC1 and mTORC2. These two complexes share four proteins (mTOR, mLST8, DEPTOR, Tti1/tel2), with only mTORC1 containing Raptor and PRAS40 and mTORC2 containing Rictor, mSin1 and Protor1/2. The complex mTORC1 (rapamycin sensitive complex) coordinates inputs from growth factors, stress, energy status, oxygen and amino acids levels to control processes such as protein and lipid synthesis and autophagy. The complex mTORC2 is insensitive to nutrients and rapamycin, but it responds to insulin signaling. It also controls ion transport and cell shape by targeting serum/glucocorticoid protein kinase (SGK1) and protein kinase (PKC- $\alpha$ ) respectively.

The canonical regulation of mTORC1 occurs through the TSC/Rheb pathway which receives signals from AKT, AMPK and IKK $\beta$  to activate the complex. Phosphorylation of mTOR at Ser2448 is carried out directly by AKT kinase as well as p70S6 kinase acting as a feedback signal. Phosphorylation at this site is a biomarker for the activation state of the PI-3 kinase pathway as well as the activation status of mTOR. Activation of mTOR leads to phosphorylation of PRAS40, raptor and DEPTOR and the consequential activation of mTORC1. Deregulated signaling of mTOR has been implicated in diseases such as cancer, metabolic syndrome, neurodegeneration and aging. Constitutive activation of PI3K-mTORC1 signaling in cancer cells inhibits autophagy, deregulates protein synthesis via 4E-BP1/eIF4E and increases de novo lipid synthesis via SREBP1. Similarly mTOR signaling is a key factor in the regulation of tissue metabolism in the normal and nutrient overload state affecting the hypothalamus, adipose tissue, the liver, skeletal muscle and pancreas. Notably, rat and human mTOR are 99.5% and 98.9% identical to mouse mTOR, respectively.

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.

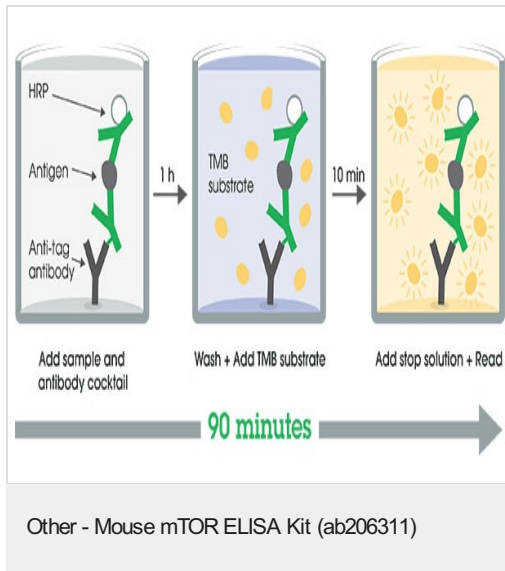
内容	1 x 96 tests
10X Wash Buffer PT (ab206977)	1 x 20ml

内容	1 x 96 tests
50X Cell Extraction Enhancer Solution (ab193971)	1 x 1ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml
Antibody Diluent 4BI	1 x 6ml
Mouse mTOR Capture Antibody (lyophilized)	1 vial
Mouse mTOR Detector Antibody (lyophilized)	1 vial
Mouse mTOR Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

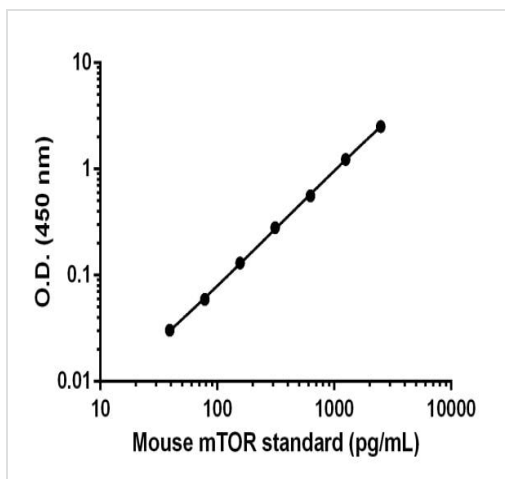
<b>機能</b>	<p>Kinase subunit of both mTORC1 and mTORC2, which regulates cell growth and survival in response to nutrient and hormonal signals. mTORC1 is activated in response to growth factors or amino-acids. Growth factor-stimulated mTORC1 activation involves AKT1-mediated phosphorylation of TSC1-TSC2, which leads to the activation of the RHEB GTPase that potentially activates the protein kinase activity of mTORC1. Amino-acid-signaling to mTORC1 requires its relocalization to the lysosomes mediated by the Ragulator complex and the Rag GTPases. Activated mTORC1 up-regulates protein synthesis by phosphorylating key regulators of mRNA translation and ribosome synthesis. mTORC1 phosphorylates EIF4EBP1 and releases it from inhibiting the elongation initiation factor 4E (eIF4E). mTORC1 phosphorylates and activates S6K1 at 'Thr-421', which then promotes protein synthesis by phosphorylating PDCD4 and targeting it for degradation. Phosphorylates MAF1 leading to attenuation of its RNA polymerase III-repressive function. mTORC2 is also activated by growth factors, but seems to be nutrient-insensitive. mTORC2 seems to function upstream of Rho GTPases to regulate the actin cytoskeleton, probably by activating one or more Rho-type guanine nucleotide exchange factors. mTORC2 promotes the serum-induced formation of stress-fibers or F-actin. mTORC2 plays a critical role in AKT1 'Ser-473' phosphorylation, which may facilitate the phosphorylation of the activation loop of AKT1 on 'Thr-308' by PDK1 which is a prerequisite for full activation. mTORC2 regulates the phosphorylation of SGK1 at 'Ser-422'. mTORC2 also modulates the phosphorylation of PRKCA on 'Ser-657'.</p>
<b>組織特異性</b>	Expressed in numerous tissues, with highest levels in testis.
<b>配列類似性</b>	<p>Belongs to the PI3/PI4-kinase family.</p> <p>Contains 1 FAT domain.</p> <p>Contains 1 FATC domain.</p> <p>Contains 7 HEAT repeats.</p> <p>Contains 1 PI3K/PI4K domain.</p>
<b>翻訳後修飾</b>	Autophosphorylated; when part of mTORC1 or mTORC2.
<b>細胞内局在</b>	Endoplasmic reticulum membrane. Golgi apparatus membrane. Mitochondrion outer membrane.

Lysosome. Cytoplasm. Nucleus > PML body. Shuttles between cytoplasm and nucleus. Accumulates in the nucleus in response to hypoxia (By similarity). Targeting to lysosomes depends on amino acid availability and RRAGA and RRAGB.

## 画像

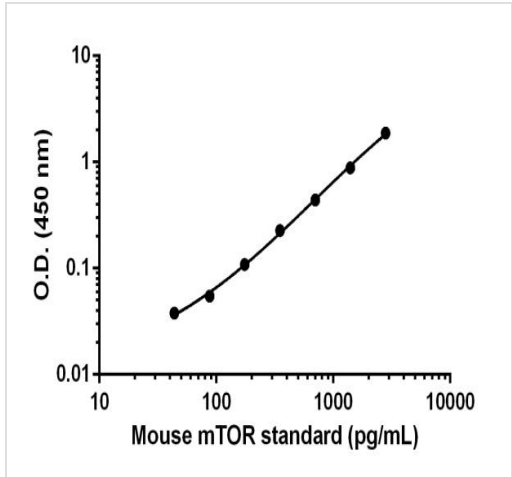


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



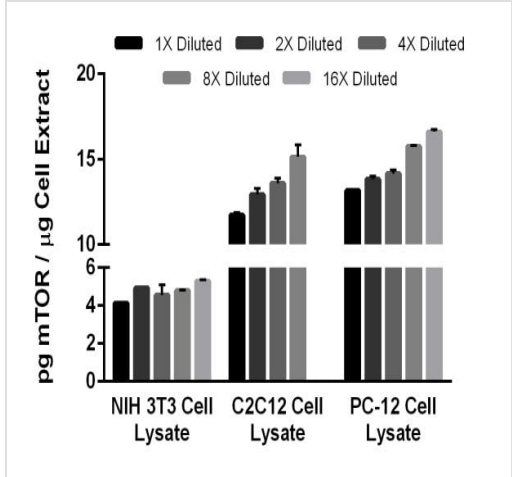
Background-subtracted data values (mean  $\pm$  SD) are graphed.

Example of the mouse mTOR standard curve in Sample Diluent NS.



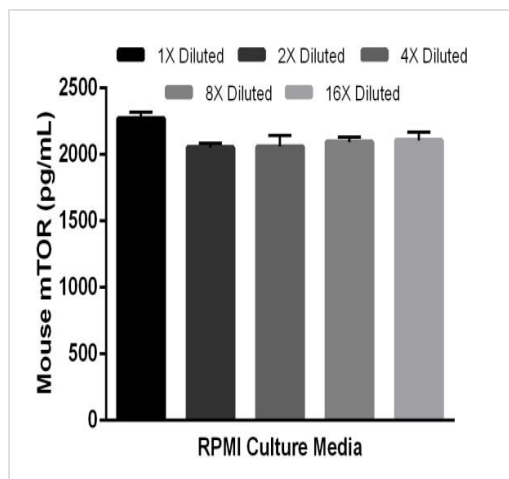
Background-subtracted data values (mean +/- SD) are graphed.

Example of the mouse mTOR standard curve in 1X Cell Extraction Buffer PTR.



Native mouse mTOR was measured in 600 µg/mL NIH 3T3 cell extract and 200 µg/mL C2C12 cell extract diluted in a 2-fold dilution series in 1X Cell Extraction Buffer PTR. Native rat mTOR was measured in 200 µg/mL PC-12 cell extract diluted in a 2-fold dilution series in 1X Cell Extraction Buffer PTR. The concentrations of mouse and rat mTOR were measured in duplicate and interpolated from the mouse mTOR standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are graphed (mean +/- SD).

Native linearity of dilution mTOR in cell extracts.



Linearity of dilution of mouse mTOR in RPMI culture media.

Recombinant mouse mTOR was spiked into 10% RPMI culture media and diluted in a 2-fold dilution series in Sample Diluent NS. The concentrations of mTOR were measured in duplicate and interpolated from the mouse mTOR standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are graphed (mean +/- SD).

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 mTOR ELISA Kit  
(ab206311)

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