

# Mouse BMP2 ELISA Kit ab119582

2 References 画像数 1

医薬用外毒物

## 製品の概要

製品名	Mouse BMP2 ELISA Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Other biological fluids, Tissue Extracts
アッセイタイプ	Sandwich (quantitative)
検出感度	< 2 pg/ml
検出範囲	62.5 pg/ml - 4000 pg/ml
ステップ	Multiple steps standard assay
種交差性	交差種: Mouse
製品の概要	<p>Abcam's mouse BMP2 <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of mouse BMP2 in bone tissue, cell culture supernatants and serum.</p> <p>A BMP2 specific rat monoclonal 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 goat, specific for BMP2 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 mouse BMP2 amount of sample captured in plate.</p>
アプリケーション	適用あり: Sandwich ELISA
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
Anti-mouse BMP2 antibody Microplate (12 x 8 wells)		1 unit

内容	ラベル	1 x 96 tests
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 130µl
Biotinylated anti-mouse BMP2 antibody		1 x 130µl
Lyophilized recombinant mouse BMP2 standard		2 x 10ng
Plate Seal		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

<b>機能</b>	Induces cartilage and bone formation.
<b>組織特異性</b>	Particularly abundant in lung, spleen and colon and in low but significant levels in heart, brain, placenta, liver, skeletal muscle, kidney, pancreas, prostate, ovary and small intestine.
<b>配列類似性</b>	Belongs to the TGF-beta family.
<b>細胞内局在</b>	Secreted.

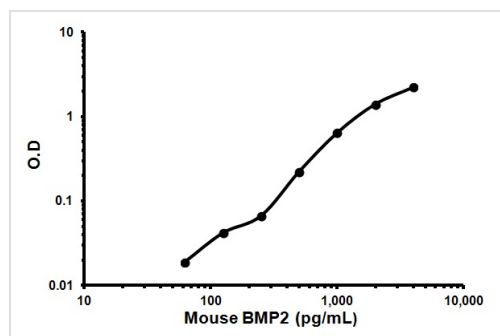
## アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab119582** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

アプリケーション	Abreviews	特記事項
Sandwich ELISA		Use at an assay dependent concentration.

## 画像



Representative Standard Curve using ab119582.

Typical Standard Curve

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

## **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