

Mouse BMP2 ELISA Kit ab119582

8 References 1 图像

概述

产品名称	小鼠BMP2 ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Other biological fluids, Tissue Extracts
检测类型	Sandwich (quantitative)
灵敏度	< 2 pg/ml
范围	62.5 pg/ml - 4000 pg/ml
实验步骤	Multiple steps standard assay
种属反应性	与反应: Mouse
产品概述	<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>
平台	Microplate

性能

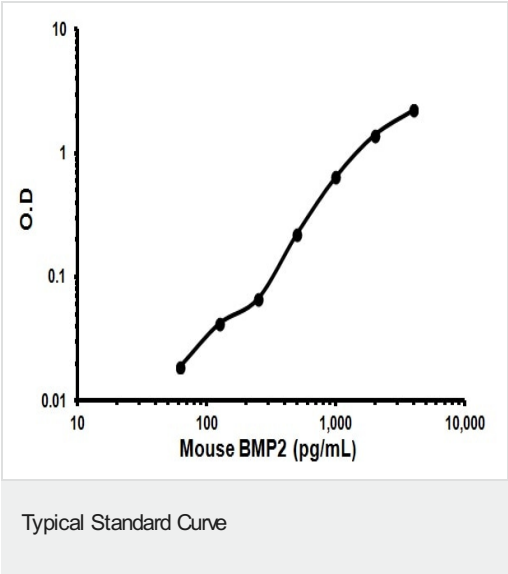
存放说明Store at -20°C. Please refer to protocols.

组件	标识符	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
Anti-mouse BMP2 antibody Microplate (12 x 8 wells)		1 unit	1 unit

组件	标识符	1 x 96 tests	1 x 96 tests
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti-mouse BMP2 antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant mouse BMP2 standard		2 x 10ng	2 x 10ng
Plate Seal		4 units	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

功能	Induces cartilage and bone formation.
组织特异性	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.
序列相似性	Belongs to the TGF-beta family.
细胞定位	Secreted.

图片



Representative Standard Curve using ab119582.

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.cn/abpromise> or contact our technical team.

Terms and conditions

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