

Mouse RANK ELISA Kit (CD265) ab119606

1 图像

概述

产品名称	小鼠RANK ELISA试剂盒(CD265)
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum
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 RANK (CD265) <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of mouse RANK (CD265) in cell culture supernatants and serum.</p> <p>A RANK (CD265) 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 RANK 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 RANK amount of sample captured in plate.</p>
经测试应用	适用于: Sandwich ELISA
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
ABC Diluent Buffer	Blue Cap	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml

组件	标识符	1 x 96 tests
Anti-Mouse RANK antibody Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 130µl
Biotinylated anti-mouse RANK antibody		1 x 130µl
Lyophilized recombinant mouse RANK standard		2 x 10ng
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

功能	Receptor for TNFSF11/RANKL/TRANCE/OPGL; essential for RANKL-mediated osteoclastogenesis. Involved in the regulation of interactions between T-cells and dendritic cells.
组织特异性	Ubiquitous expression with high levels in skeletal muscle, thymus, liver, colon, small intestine and adrenal gland.
疾病相关	<p>Defects in TNFRSF11A are the cause of familial expansile osteolysis (FEO) [MIM:174810]. FEO is a rare autosomal dominant bone disorder characterized by focal areas of increased bone remodeling. The osteolytic lesions develop usually in the long bones during early adulthood. FEO is often associated with early onset deafness and loss of dentition.</p> <p>Defects in TNFRSF11A are a cause of Paget disease of bone type 2 (PDB2) [MIM:602080]; also known as familial Paget disease of bone. PDB2 is a bone-remodeling disorder with clinical similarities to FEO. Unlike FEO, however, affected individuals have involvement of the axial skeleton with lesions in the spine, pelvis and skull.</p> <p>Defects in TNFRSF11A are the cause of osteopetrosis autosomal recessive type 7 (OPTB7) [MIM:612301]; also called osteoclast-poor osteopetrosis with hypogammaglobulinemia. Osteopetrosis is a rare genetic disease characterized by abnormally dense bone, due to defective resorption of immature bone. The disorder occurs in two forms: a severe autosomal recessive form occurring in utero, infancy, or childhood, and a benign autosomal dominant form occurring in adolescence or adulthood. OPTB7 is characterized by paucity of osteoclasts, suggesting a molecular defect in osteoclast development. OPTB7 is associated with hypogammaglobulinemia.</p>
序列相似性	Contains 4 TNFR-Cys repeats.
细胞定位	Membrane.

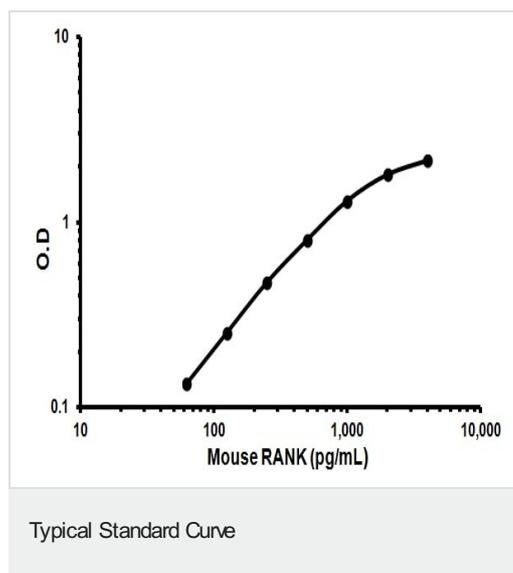
应用

The Abpromise guarantee **Abpromise™** 承诺保证使用 ab119606 于以下的经测试应用

“应用说明”部分 下显示的仅为推荐的起始稀释度;实际最佳的稀释度/浓度应由使用者检定。

应用	Ab评论	说明
Sandwich ELISA		Use at an assay dependent concentration.

图片



Representative standard curve using ab119606

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