

Product datasheet

Mouse Osteoprotegerin Matched Antibody Pair Kit
ab215078

1 Image

Overview

Product name	Mouse Osteoprotegerin Matched Antibody Pair Kit
Detection method	Colorimetric
Assay type	ELISA set
Range	15.62 pg/ml - 1000 pg/ml
Species reactivity	Reacts with: Mouse
Product overview	<p>Matched Antibody Pair kit is a titrated unlabeled capture antibody, a titrated biotin-labeled detector and a calibrated protein standard. The Matched Antibody Pair Kit can be used to quantify native and recombinant mouse Osteoprotegerin.</p> <p>Both capture and detector antibodies are rabbit monoclonal antibodies.</p> <p>Optimization of the kit reagents to sample type, immunoassay format or instrumentation may be required. Guidelines for use of this kit in a standard 96-well microplate sandwich ELISA using HRP/TMB system of colorimetric detection is described in this assay procedure for the purposes of quantification.</p> <p>Additional protocol information and tips on the use of the Matched Antibody Pair kits for sandwich ELISA can be found on our website.</p> <p>For additional information on the performance of the antibody pair used in this kit, please see our equivalent SimpleStep ELISA® (ab203365) which uses the same antibody pair.</p> <p>To receive an electronic copy of the Certificate of Analysis, please send an email with "CoA for matched antibody pair kit" in the subject line and the desired product number and lot number in the body of the email.</p> <p>Buffer information: The capture antibody is glycerol free. The detector antibody contains glycerol.</p>
Tested applications	Suitable for: ELISA
Platform	Reagents

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	10 x 96 tests	2 x 96 tests
Mouse Osteoprotegerin Capture Antibody	1 x 100µg	1 x 20µg
Mouse Osteoprotegerin Detector Antibody	1 x 25µg	1 x 5µg
Mouse Osteoprotegerin Lyophilized Protein	1 vial	1 vial

Function Acts as decoy receptor for RANKL and thereby neutralizes its function in osteoclastogenesis. Inhibits the activation of osteoclasts and promotes osteoclast apoptosis in vitro. Bone homeostasis seems to depend on the local RANKL/OPG ratio. May also play a role in preventing arterial calcification. May act as decoy receptor for TRAIL and protect against apoptosis. TRAIL binding blocks the inhibition of osteoclastogenesis.

Tissue specificity Highly expressed in adult lung, heart, kidney, liver, spleen, thymus, prostate, ovary, small intestine, thyroid, lymph node, trachea, adrenal gland, testis, and bone marrow. Detected at very low levels in brain, placenta and skeletal muscle. Highly expressed in fetal kidney, liver and lung.

Involvement in disease Defects in TNFRSF11B are the cause of juvenile Paget disease (JPD) [MIM:239000]; also known as hyperostosis corticalis deformans juvenilis or hereditary hyperphosphatasia or chronic congenital idiopathic hyperphosphatasia. JPD is a rare autosomal recessive osteopathy that presents in infancy or early childhood. The disorder is characterized by rapidly remodeling woven bone, osteopenia, debilitating fractures, and deformities due to a markedly accelerated rate of bone remodeling throughout the skeleton. Approximately 40 cases of JPD have been reported worldwide. Unless it is treated with drugs that block osteoclast-mediated skeletal resorption, the disease can be fatal.

Sequence similarities Contains 2 death domains.
Contains 4 TNFR-Cys repeats.

Post-translational modifications N-glycosylated. Contains sialic acid residues.
The N-terminus is blocked.

Cellular localization Secreted.

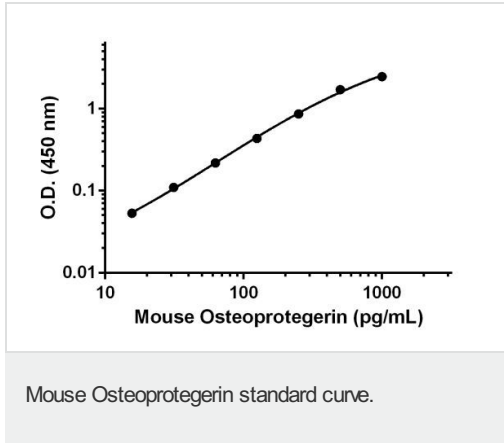
Applications

Our [Abpromise guarantee](#) covers the use of **ab215078** in the following tested applications.

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

Application	Abreviews	Notes
ELISA		Use at an assay dependent concentration.

Images



Standard calibration curve. Background subtracted values are graphed.

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.com/abpromise> or contact our technical team.

Terms and conditions

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