

Product datasheet

Mouse CXCL11 Matched Antibody Pair Kit ab212066

1 Image

Overview

Product name	Mouse CXCL11 Matched Antibody Pair Kit
Detection method	Colorimetric
Assay type	ELISA set
Range	62.5 pg/ml - 4000 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 CXCL11. 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® (ab204519) 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 CXCL11 Capture Antibody	1 x 100µg	1 x 20µg
Mouse CXCL11 Detector Antibody	1 x 25µg	1 x 5µg
Mouse CXCL11 Lyophilized Protein	1 vial	1 vial

Function Chemotactic for interleukin-activated T-cells but not unstimulated T-cells, neutrophils or monocytes. Induces calcium release in activated T-cells. Binds to CXCR3. May play an important role in CNS diseases which involve T-cell recruitment. May play a role in skin immune responses.

Tissue specificity High levels in peripheral blood leukocytes, pancreas and liver astrocytes. Moderate levels in thymus, spleen and lung. Low levels in placenta, prostate and small intestine. Also found in epidermal basal layer keratinocytes in skin disorders.

Sequence similarities Belongs to the intercrine alpha (chemokine CxC) family.

Cellular localization Secreted.

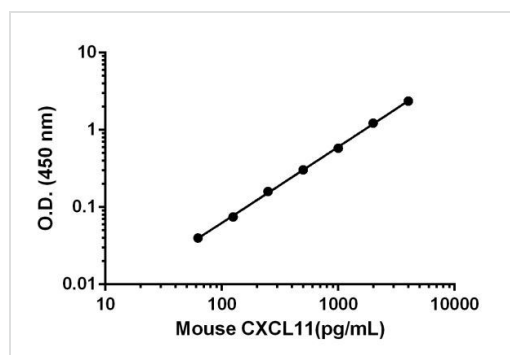
Applications

Our [Abpromise guarantee](#) covers the use of **ab212066** 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.

Mouse CXCL11 standard curve

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