

Product datasheet

Mouse CXCL1 Matched Antibody Pair Kit ab213742

1 Image

Overview

---

<b>Product name</b>	Mouse CXCL1 Matched Antibody Pair Kit
<b>Detection method</b>	Colorimetric
<b>Assay type</b>	ELISA set
<b>Range</b>	7.8 pg/ml - 500 pg/ml
<b>Species reactivity</b>	<b>Reacts with:</b> Mouse
<b>Product overview</b>	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 CXCL1. Both capture and detector antibodies are rabbit monoclonal antibodies.

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.

Additional protocol information and tip on the use of the Matched Antibody Pair kits for sandwich ELISA can be found on our [website](#).

Buffer information: The capture antibody is glycerol free. The detector antibody contains glycerol.

**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.**

<b>Platform</b>	Microplate
-----------------	------------

Properties

---

<b>Storage instructions</b>	Store at -20°C. Please refer to protocols.
-----------------------------	--

Components	10 x 96 tests	2 x 96 tests
Mouse CXCL1 Capture Antibody	1 x 100µg	1 x 20µg
Mouse CXCL1 Detector Antibody	1 x 25µg	1 x 5µg
Mouse CXCL1 Lyophilized Protein	1 vial	1 vial

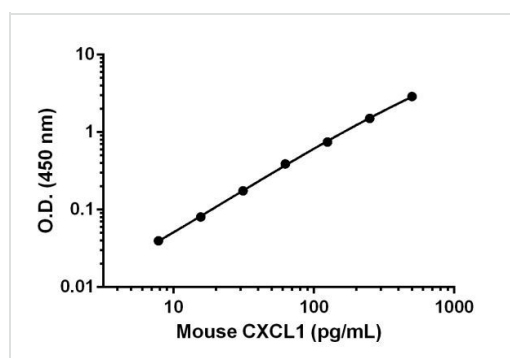
**Function** Has chemotactic activity for neutrophils. May play a role in inflammation and exerts its effects on endothelial cells in an autocrine fashion. In vitro, the processed forms GRO-alpha(4-73), GRO-alpha(5-73) and GRO-alpha(6-73) show a 30-fold higher chemotactic activity.

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

**Post-translational modifications** N-terminal processed forms GRO-alpha(4-73), GRO-alpha(5-73) and GRO-alpha(6-73) are produced by proteolytic cleavage after secretion from peripheral blood monocytes.

**Cellular localization** Secreted.

## Images



Standard calibration curve. Background subtracted values are graphed. Data provided for demonstration purposes only.

Mouse CXCL1 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