

Mouse CXCL15 ELISA Kit ab193720

1 Image

Overview

Product name	Mouse CXCL15 ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 10%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 12%
Sample type	Cell culture supernatant, Serum, Plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	0.8 ng/ml				
Range	0.8 ng/ml - 200 ng/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		120.1		102% - 128%
	Plasma		134		118% - 150%
	Cell culture media		135.5		124% - 144%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Mouse				
Product overview	Abcam's CXCL15 mouse ELISA Kit (ab193720) is an <i>in vitro</i> enzyme-linked immunosorbent				

CXCL15 present in the samples. The wells are washed and biotinylated anti-mouse CXCL15 antibody is added. After washing away any unbound biotinylated antibody, an HRP-conjugated streptavidin is pipetted to the wells. After incubation, the wells are again washed, followed by the addition of a TMB substrate solution to the wells. Color will develop in proportion to the amount of CXCL15 bound in each well. Addition of the Stop Solution will change the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**Platform** Pre-coated microplate (12 x 8 well strips)

## Properties

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

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent D	1 x 15ml
800X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Biotinylated Mouse CXCL15 detection antibody	2 vials
mouse CXCL15 Standards (lyophilized)	2 vials
Pre-coated mouse CXCL15 Microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

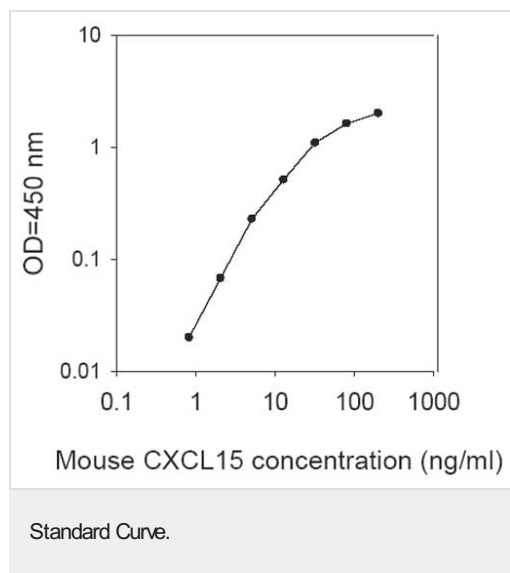
**Function** Chemotactic for neutrophils. Involved in lung-specific neutrophil trafficking during normal and inflammatory conditions.

**Tissue specificity** Expression restricted to the lung, produced by bronchoepithelial cells and is released into the airways. Expressed at low levels in fetal lung.

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

**Cellular localization** Secreted.

## Images



Typical Standard Curve using ab193720 CXCL15 Mouse ELISA Kit.

**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