

Product datasheet

Rat CINC-1 ELISA Kit (CXCL1) ab155462

[2 Images](#)

Overview

Product name Rat CINC-1 ELISA Kit (CXCL1)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
--------	---	------	----	-----

Intra-assay < 10%

Inter-assay

Sample	n	Mean	SD	CV%
--------	---	------	----	-----

Inter-assay < 12%

Sample type Cell culture supernatant, Plasma

Assay type Sandwich (quantitative)

Sensitivity 0.05 ng/ml

Range 0.047 ng/ml - 15 ng/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	112.5	99% - 122%
Plasma	107.9	97% - 125%

Assay time 5h 00m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Rat

Product overview

Abcam's CINC-1 (CXCL-1) Rat ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay designed for the quantitative measurement of rat CINC-1 in serum, plasma and cell culture supernatants. (Collect plasma using EDTA and heparin as an anticoagulant). Citrate is not recommended as anticoagulants for use in this assay. Rat CINC-1 concentration in normal plasma is very low and it may not be detected in this assay.

This assay employs an antibody specific for rat CINC-1 coated on a 96-well plate. Standards and samples are pipetted into the wells and CINC-1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-rat CINC-1 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of CINC-1 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

We have not been able to detect the endogenous Rat CINC-1 in normal serum with ab155462, only in serum spiked with Rat CINC-1.

Notes Sensitivity of the assay is dependant on the Assay Diluent used. Please refer to Protocol Booklet for more details.

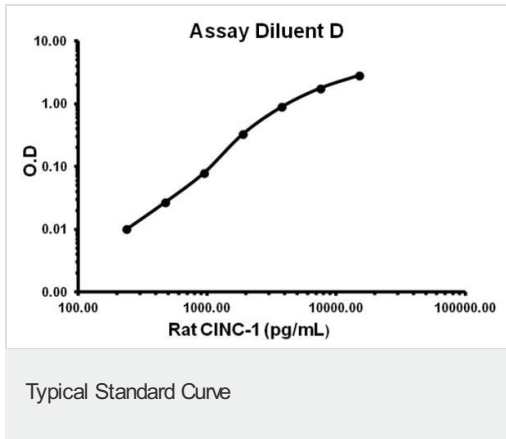
Platform Microplate

Properties

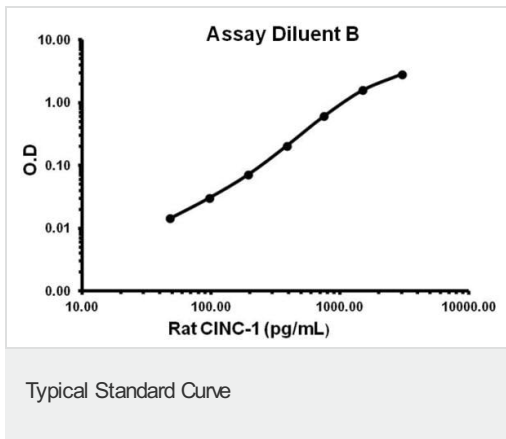
Storage instructions Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer Concentrate	1 x 25ml
5X Assay Diluent B	1 x 15ml
5X Assay Diluent D	1 x 15ml
Biotinylated anti-rat CINC-1	2 vials
CINC-1 Microplate (12 x 8 wells)	1 unit
Recombinant Rat CINC-1 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Images



Representative standard curve using ab155462 - Assay Diluent D.



Representative standard curve using ab155462 - Assay Diluent B.

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