

Rat CINC-3 ELISA Kit (CXCL2) ab155463

1 References 2 Images

Overview

[illegible]

Abcam's CINC-3 (CXCL2) Rat ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay designed for the quantitative measurement of rat CINC-3 (CXCL2) in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for CINC-3 coated on a 96-well plate. Standards and samples are pipetted into the wells and CINC-3 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Rat CINC-3 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-3 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Platform Microplate

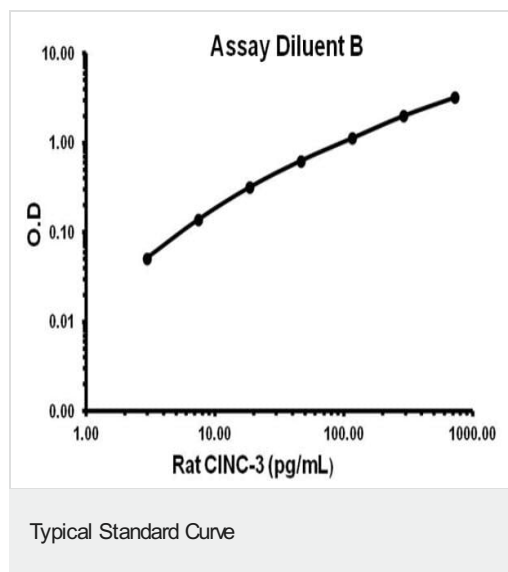
Properties

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

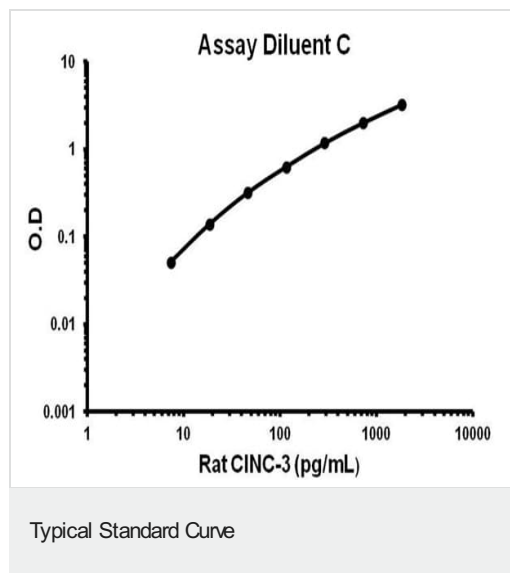
Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
600X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent C	1 x 30ml
Biotinylated anti-rat CINC-3	2 vials
CINC-3 Microplate (12 x 8 wells)	1 unit
Recombinant Rat CINC-3 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Cellular localization Secreted

Images



Representative standard curve using ab155463 - Assay Diluent B.



Representative standard curve using ab155463 - Assay Diluent C.

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