

Mouse CT-1 ELISA Kit ab155453

1 Image

Overview

|                    |  |   |           |    |            |
|--------------------|--|---|-----------|----|------------|
| Product name       | Mouse CT-1 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        | < 0.2 ng/ml  |   |           |    |            |
| Range              | 0.25 ng/ml - 60 ng/ml  |   |           |    |            |
| Recovery           | Sample specific recovery   |   |           |    |            |
|                    | Sample type  |   | Average % |    | Range      |
|                    | Serum  |   | 78.13     |    | 70% - 88%  |
|                    | Plasma   |   | 83.28     |    | 73% - 96%  |
|                    | Cell culture media   |   | 92.96     |    | 84% - 100% |
| Assay duration     | Multiple steps standard assay  |   |           |    |            |
| Species reactivity | Reacts with: Mouse   |   |           |    |            |
| Product overview   | <p>Abcam's CT-1 Mouse ELISA Kit (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay designed for the quantitative measurement of Mouse CT-1 in serum, plasma and cell culture supernatants.</p> <p>This assay employs an antibody specific for CT-1 coated on a 96-well plate. Standards and samples are pipetted into the wells and CT-1 present in a sample is bound to the wells by the</p> |   |           |    |            |

immobilized antibody. The wells are washed and biotinylated anti-Mouse CT-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 CT-1 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 |
|---|--------------|
| 200X HRP-Streptavidin Concentrate             | 1 x 200µl    |
| 20X Wash Buffer                               | 1 x 25ml     |
| 5X Assay Diluent B                            | 1 x 15ml     |
| Assay Diluent C                               | 1 x 30ml     |
| Biotinylated anti-mouse CT-1                  | 2 vials      |
| CT-1 Microplate (12 x 8 wells)                | 1 unit       |
| Recombinant Mouse CT-1 Standard (lyophilized) | 2 vials      |
| Stop Solution                                 | 1 x 8ml      |
| TMB One-Step Substrate Reagent                | 1 x 12ml     |

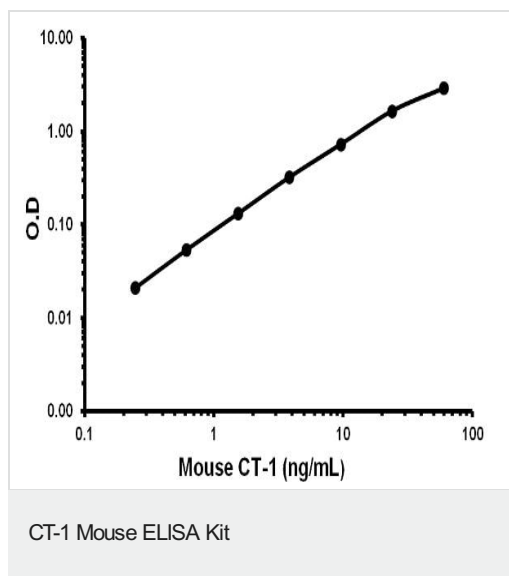
**Function** Induces cardiac myocyte hypertrophy in vitro. Binds to and activates the ILST/gp130 receptor.

**Tissue specificity** Highly expressed in heart, skeletal muscle, prostate and ovary. Lower levels in lung, kidney, pancreas, thymus, testis and small intestine. Little or no expression in brain, placenta, liver, spleen, colon or peripheral blood leukocytes.

**Sequence similarities** Belongs to the IL-6 superfamily.

**Cellular localization** Secreted.

## Images



Representative standard curve using ab155453

**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