

Mouse TCA-3 ELISA Kit (CCL1) ab155460

2 Images

Overview

Product name	Mouse TCA-3 ELISA Kit (CCL1)				
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, Serum, Plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	< 30 pg/ml				
Range	0.033 ng/ml - 8 ng/ml				
Recovery	Sample specific recovery				
	Sample type		Average %	Range	
	Cell culture supernatant		91.6	77% - 103%	
	Serum		127.7	119% - 135%	
	Plasma		126.3	109% - 139%	
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Mouse				
Product overview	Abcam's TCA-3 (CCL1) Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i>				

the immobilized antibody. The wells are washed and biotinylated anti-mouse TCA-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 colour develops in proportion to the amount of TCA-3 bound. The Stop Solution changes the colour from blue to yellow, and the intensity of the colour is measured at 450 nm.

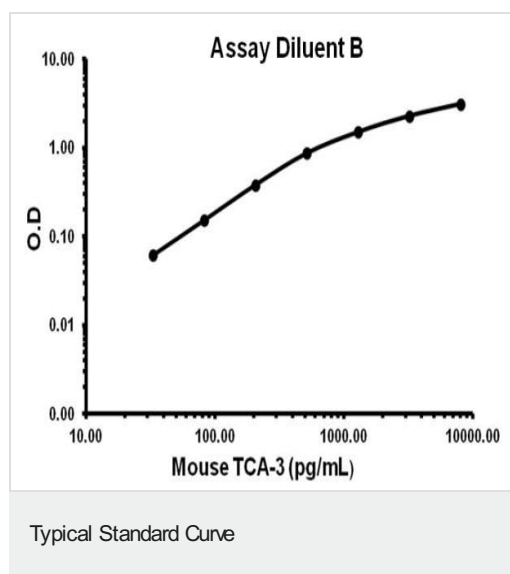
Platform Microplate

Properties

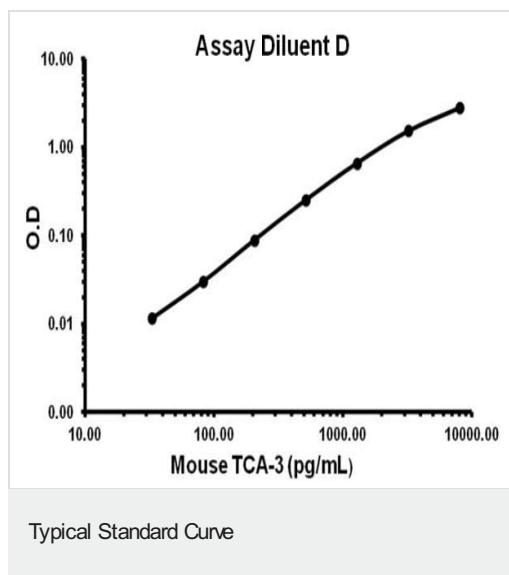
Storage instructions Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
5X Assay Diluent D	1 x 15ml
700X HRP-Streptavidin Concentrate	1 x 200µl
Biotinylated anti-mouse TCA-3	2 vials
Recombinant Mouse TCA-3 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TCA-3 Microplate (12 x 8 wells)	1 unit
TMB One-Step Substrate Reagent	1 x 12ml

Images



Representative standard curve using ab155460 - Assay Diluent B.



Representative standard curve using ab155460 - Assay Diluent D.

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