

Calretinin ELISA Kit ab233616

Recombinant SimpleStep ELISA

6 Images

Overview

Product name	Calretinin ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Brain	6			4.5%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Brain	3			10.5%
Sample type	Milk, Tissue Extracts, Mouse brain extract				
Assay type	Sandwich (quantitative)				
Sensitivity	4.2 pg/ml				
Range	234 pg/ml - 15000 pg/ml				
Recovery	Sample specific recovery				
	Sample type		Average %	Range	
	Milk		111	103% - 115%	
	Tissue Extracts		107	98% - 114%	
	Mouse brain extract		106	97% - 115%	
Assay time	1h 30m				
Assay duration	One step assay				
Species reactivity	Reacts with: Mouse, Rat, Human				
Product overview	Calretinin ELISA Kit (ab233616) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Calretinin protein in milk, mouse brain extract, and tissue extracts. It				

uses our proprietary SimpleStep ELISA® technology. Quantitate Calretinin with 4.2 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

**Notes**

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

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

## Properties

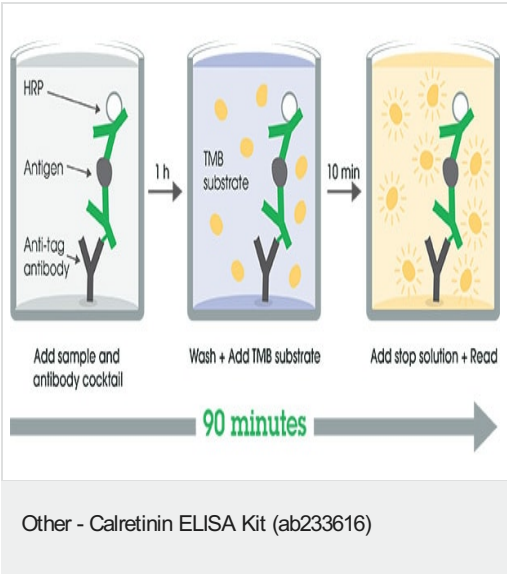
**Storage instructions** Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Calretinin Capture Antibody	1 x 600µl
10X Calretinin Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
50X Cell Extraction Enhancer Solution (ab193971)	1 x 1ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml
Antibody Diluent 5BR	1 x 6ml
Calretinin Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 12ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit

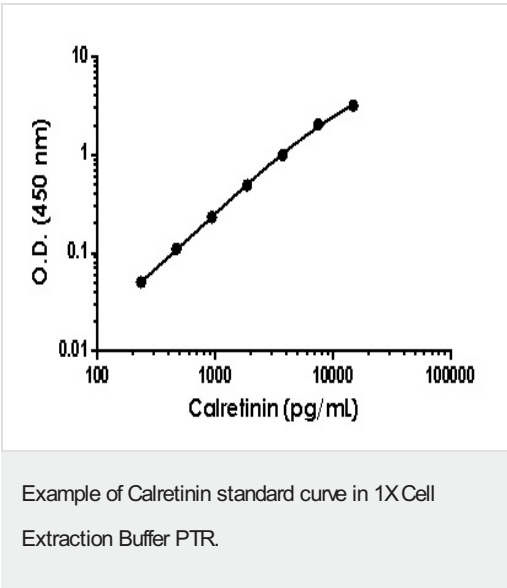
Components	1 x 96 tests
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

Function	Calretinin is a calcium-binding protein which is abundant in auditory neurons.
Tissue specificity	Brain.
Sequence similarities	Belongs to the calbindin family. Contains 6 EF-hand domains.

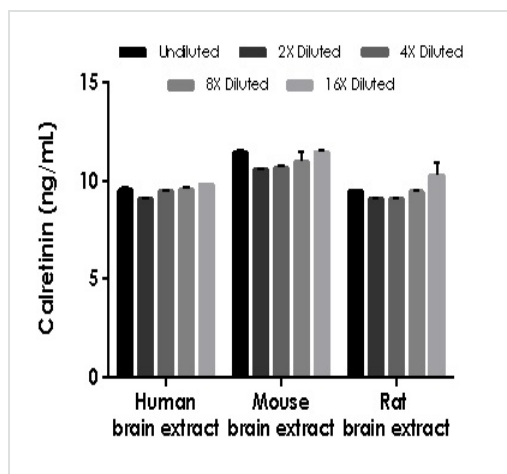
Images



SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.

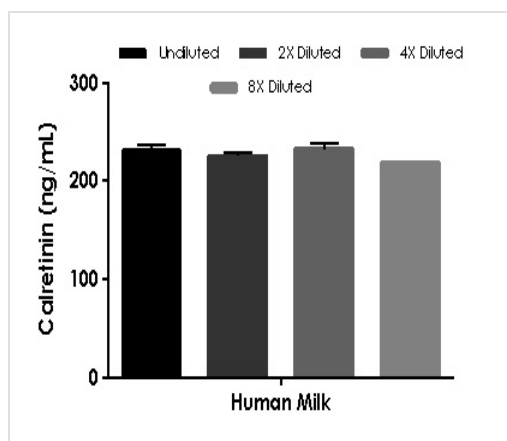


The Calretinin standard curve was prepared as described in Section 10.



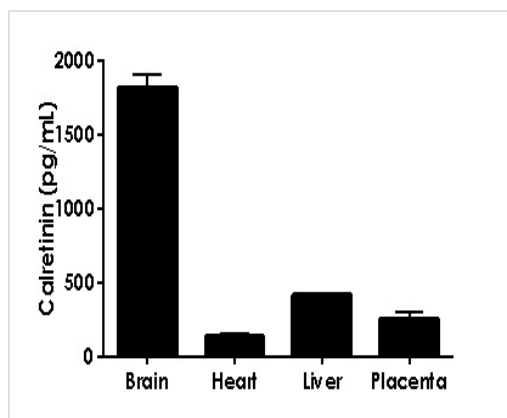
Interpolated concentrations of native Calretinin in human, mouse, and rat brain extract samples.

The concentrations of Calretinin were measured in duplicates, interpolated from the Calretinin standard curves and corrected for sample dilution. Undiluted samples are as follows: brain extract 12.5 µg/mL. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2).



Interpolated concentrations of native Calretinin in human breast milk samples.

The concentrations of Calretinin were measured in duplicates, interpolated from the Calretinin standard curves and corrected for sample dilution. Undiluted samples are as follows: milk 1.25%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Calretinin concentration was determined to be 227 ng/mL in neat human breast milk.



Relative expression of calretinin in normal tissue extracts

Calretinin expression in non-neural and non-malignant tissues are expected to be significantly lower than neural tissues. Interpolated concentrations of native Calretinin in human brain, heart, liver, and placenta extract samples. The concentrations of Calretinin were measured in duplicates, interpolated from the Calretinin standard curves and corrected for sample dilution. Undiluted samples are as follows: brain extract 3 µg/mL, heart, liver, and placenta extract 100 µg/mL. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2).

Powered by  
recombinant antibodies



**Research with confidence**  
Consistent and reproducible results



**Long-term and scalable supply**  
Recombinant technology



**Success from the first experiment**  
Confirmed specificity



**Ethical standards compliant**  
Animal-free production

Sandwich ELISA - Calretinin ELISA Kit (ab233616)

To learn more about the advantages of recombinant antibodies see [here](#).

**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