

## Product datasheet

# Human TACE ELISA Kit (ADAM17) ab113321

[1 References](#) [1 Image](#)

### Overview

<b>Product name</b>	Human TACE ELISA Kit (ADAM17)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell culture supernatant, Serum, Plasma
<b>Assay type</b>	Sandwich (quantitative)
<b>Sensitivity</b>	< 70 pg/ml
<b>Range</b>	78.15 pg/ml - 5000 pg/ml
<b>Recovery</b>	100 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	98.65	89% - 124%
Serum	112.3	85% - 132%
Plasma	101.1	72% - 128%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Product overview** Abcam's TACE Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human TACE in serum, plasma and cell culture supernatants.

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

**Notes** Optimization may be required with urine samples.

The standard is affinity purified.

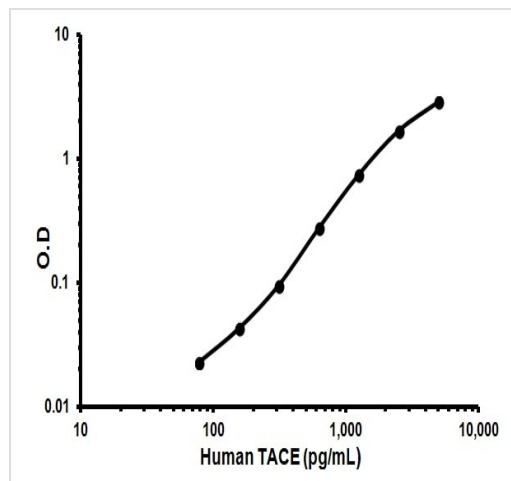
**Platform**

Microplate

**Properties****Storage instructions**

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

Components	Identifier	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-Human TACE	Item F	2 vials
Recombinant Human TACE Standard (lyophilized)		2 vials
Stop Solution		1 x 8ml
TACE Microplate (12 x 8 wells)	Item A	1 unit
TMB One-Step Substrate Reagent		1 x 12ml

**Images**

Representative standard curve using ab113321

Typical Standard Curve

**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