

# Human Thrombomodulin ELISA Set (without plates) ab47358

[1 Image](#)

### Overview

<b>Product name</b>	Human Thrombomodulin ELISA Set (without plates)
<b>Sample type</b>	Cell culture supernatant, Serum, Plasma
<b>Assay type</b>	Sandwich (quantitative)
<b>Sensitivity</b>	< 0.31 ng/ml
<b>Range</b>	0.62 ng/ml - 20 ng/ml
<b>Assay time</b>	1h 45m
<b>Assay duration</b>	Multiple steps standard assay
<b>Species reactivity</b>	<b>Reacts with:</b> Human
<b>Product overview</b>	<p>Abcam's Thrombomodulin Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of Thrombomodulin in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>This ELISA Set is intended to measure human Thrombomodulin in cell -culture supernatants, as well as in serum and plasma samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.</p> <p>These reagents are intended for research use only, not for diagnostic purposes.</p> <p>The Thrombomodulin ELISA Set includes both the capture and the detection antibodies as well as Horse-Radish peroxidase conjugated Streptavidin (HRP-Strep) and TMB which are the reagents used in the colour development step.</p> <p>Recognizes both natural and recombinant human Thrombomodulin.</p>
<b>Tested applications</b>	<b>Suitable for:</b> Sandwich ELISA
<b>Platform</b>	Reagents

### Properties

**Storage instructions**

Store at +4°C. Please refer to protocols.

Components	10 x 96 tests	5 x 96 tests	15 x 96 tests
Chromogen TMB Substrate Solution	4 vials	2 vials	6 vials
Detection biotinylated anti-Thrombomodulin antibody	2 x 0.55ml	1 x 0.55ml	3 x 0.55ml
Streptavidin-HRP	2 vials	1 vial	3 vials
Thrombomodulin Capture Antibody	2 x 0.1ml	1 x 0.1ml	3 x 0.1ml
Thrombomodulin standard	10 vials	5 vials	15 vials

**Relevance**

Thrombomodulin is a specific endothelial cell receptor that forms a 1:1 stoichiometric complex with thrombin. Endothelial cells are unique in synthesizing thrombomodulin. The thrombomodulin/thrombin complex is responsible for the conversion of protein C to the activated protein C (protein Ca), which degrades clotting factors Va and VIIIa and reduces the amount of thrombin generated. Defects in Thrombomodulin may be a cause of thromboembolic disease.

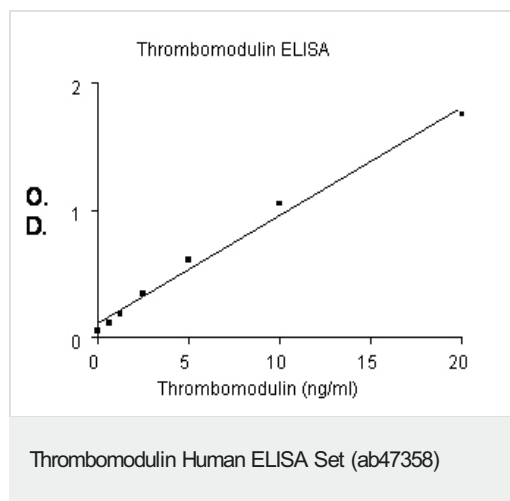
**Cellular localization**

single-pass type I membrane protein

**Applications****The Abpromise guarantee**Our **Abpromise guarantee** covers the use of ab47358 in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
<b>Sandwich ELISA</b>		Use at an assay dependent concentration.

**Images**

Thrombomodulin Standard curve ranging from 0.625 to 20 ng/ml

**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