



EIA for Quantitative Determination of anti-Thrombin-Receptor-Antibodies

Introduction

The Thrombin-Receptor (PAR-1, F2R) is a seven transmembrane, G-protein- coupled receptor which is located in thrombocytes and endothelial cells. These enzymes cleave the N-terminus of the receptor, which in turn acts as a tethered ligand. In the cleaved state, part of the receptor itself acts as the agonist, causing a physiological response. The occurrence of auto- antibodies against thrombin receptor is associated with the existence of different vascular diseases.

The CellTrend anti-Thrombin-Receptor -Antibody-EIA is designed for the determination of Antibodies (IgG) against the Thrombin-Receptor in serum and plasma

Performance Characteristics

Standard curve:

5 standards between 2,5 U/ml and 40 U/ml

cut off:

> 6.0 U/ml no PE

5.0 – 6.0 U/ml at risk

< 5.0 PE

(Preeclampsia, Dechend et al, ISSHP 2015)

Sample materials:

Serum, Plasma

Intraassay-Precision:

(n=10)

Sample 1 (18.2 U/ml): 1.9%

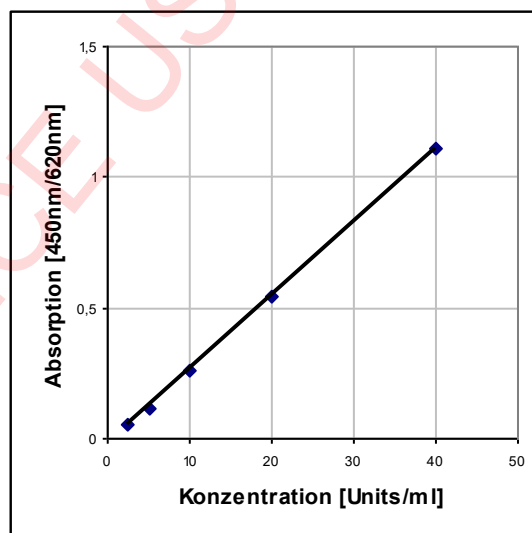
Interassay-Precision:

6.62%

Principle of the test

The CellTrend anti-Thrombin-Receptor-Antibody-EIA is an antibody screening test. The Thrombin-Receptor has been pre-coated onto a microtiter plate. During the first incubation the anti-Thrombin-Receptor-Antibodies of the samples are immobilised on the plate. The auto-antibodies are detected with a POD labeled anti-human IgG antibody. In the following enzymatic substrate reaction the intensity of the colour correlates with the concentration and/ or avidity of anti-Thrombin-Receptor-Antibody.

Typical Standard Curve



Assay Procedure

Incubation of samples/ standards/ controls	100 µl	120 min, 4°C
Wash		
Incubation of detection antibody	100 µl	1 hr, room temperature
Wash		
Substrate incubation	100 µl/well	20 min, room temperature
Add Stopp solution	100 µl/well	
Read at 450nm		

Order informations

Product	Catalog number	Price (€)
EIA for Quantitative Determination of anti-PAR-1-AB, 1x96 determ.	12200	1100,-

CE *In vitro*-Diagnostikum