



EIA for Quantitative Determination of anti-Thrombin-Receptor-like-1-Antibodies

Introduction

The Thrombin-Receptor-like-1 (PAR-2, F2RL1) is a seven transmembrane, G-protein- coupled receptor which is located in thrombocytes and t-lymphocytes. The protease trypsin splits off a N-terminal end of the receptor. It is activated by the arised domain. The occurrence of auto- antibodies against thrombin-receptor-like-1 is associated with the existence of different vascular diseases.

The CellTrend anti-Thrombin-Receptor-like-1-Antibody-EIA is designed for the determination of Antibodies (IgG) against the Thrombin-Receptor in serum and plasma.

Principle of the test

The CellTrend anti-Thrombin-Receptor-like-1-Antibody-EIA is an antibody screening test. The Thrombin-Receptor-like-1 has been pre-coated onto a microtiter plate. During the first incubation the anti-Thrombin-Receptor-like-1-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-like-1-Antibody.

Performance Characteristics

Standard curve:

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

cut off:

20 U/ml

Sample materials:

Serum, Plasma

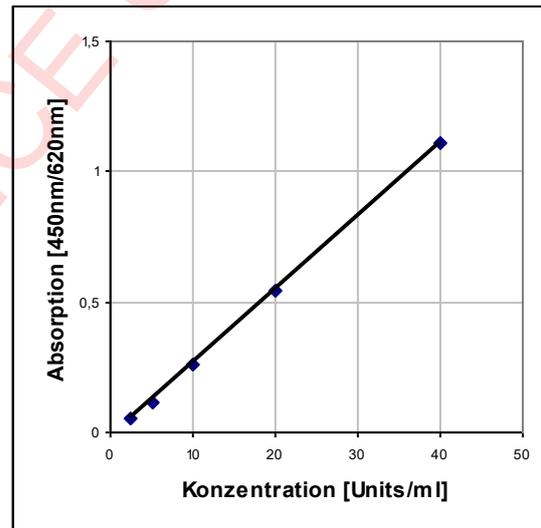
Intraassay-Precision:

7,83%

Interassay-Precision:

9,12%

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-2, 1x96 determ.	12300	1100,-

FOR REFERENCE USE ONLY