



# EIA for Quantitative Determination of anti-ACE-II-Antibodies

distributed in the US/Canada by:  
**EAGLE BIOSCIENCES, INC.**  
 20A NW Blvd, Suite 112, Nashua, NH 03063  
 Phone: 617-419-2019 FAX: 617-419-1110  
[www.EagleBio.com](http://www.EagleBio.com) • [info@eaglebio.com](mailto:info@eaglebio.com)



## Introduction

Angiotensin-Converting-Enzyme 2 (ACE-II) is a zinc containing metalloenzyme expressed on endothelial and other cells. ACE-II is an enzyme expressed by cells in the lungs, arteries, heart, kidney, and intestines. ACE-II lowers blood pressure by catalysing the hydrolysis of angiotensin II (Ang-II) into angiotensin (1-7). ACE-II counters the activity of the related angiotensin-converting enzyme by reducing Ang-II and increasing Ang(1-7). The CellTrend anti-ACE-II-Antibody EIA is designed for the determination of IgG antibodies against the ACE-II in serum and plasma.

## Principle of the test

The CellTrend Angiotensin-converting enzyme 2 (ACE-II)-EIA is an antibody screening test. ACE-II enzyme has been pre-coated onto a microtiter plate. During the first incubation the anti-ACE-II-antibodies of the samples are immobilized on the plate. The autoantibodies are detected with a HRP 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-ACE-II antibodies.

## Performance Characteristics

### Standard curve:

6 standards between 1.25 U/ml and 40 U/ml

### Healthy controls:

Mean: 12.9 Units  
SEM: 1.9 Units/ml

### Sample materials:

Serum, Plasma

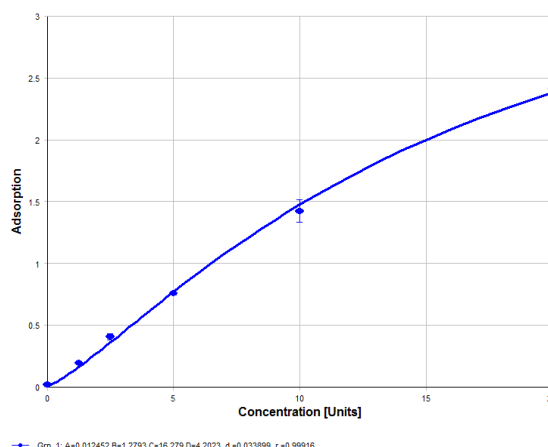
### Intraassay-Precision (n=10):

Sample 1 (11.1 U/ml): 6.8%

### Interassay-Precision (n=10):

Sample 1 (12.0 U/ml): 11.2%

## Typical Standard Curve



## Assay Procedure

Incubation of samples/ standards/ controls	100 µl	120 min, 4°C
Wash		
Incubation of detection antibody	100 µl	60 min, room temperature
Wash		
Substrate incubation	100 µl/well	20 min, room temperature
Add Stopp solution	100 µl/well	
Read at 450nm		

## Order information

Product	Bestell-Nummer	Preis (€)
EIA Quantitative Determination of anti-ACEII-IgG-Antibodies, 1x96 Best.	16000	950,-