



EIA for Quantitative Determination of anti-Muscarinic Cholinergic Receptor 5 (M5)-Antibodies

Introduction

Muscarinic cholinergic receptors, or mAChRs, are acetylcholine receptors that form G protein-receptor complexes in the cell membranes of certain neurons and other cells.

The CellTrend anti-muscarinic cholinergic receptor 5 (M5)-antibody EIA is designed for the determination of antibodies against the muscarinic cholinergic receptor 5 (M5) in serum.

Principle of the assay

The CellTrend muscarinic cholinergic receptor 5 (M5)-EIA is an antibody screening test. M5 receptor has been pre-coated onto a microtiter plate. During the first incubation the anti-muscarinic cholinergic receptor 5-antibodies of the samples are immobilised on the plate. The autoantibodies 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-muscarinic cholinergic receptor 5-antibody.

Performance Characteristics

Standard curve:

5 standards between 2.5 U/ml and 40 U/ml

cut off: > 14.2 Units/ml positive

< 14.2 Units/ml negative

Sample materials:

Serum, Plasma

Intraassay-Precision:

(n=10)

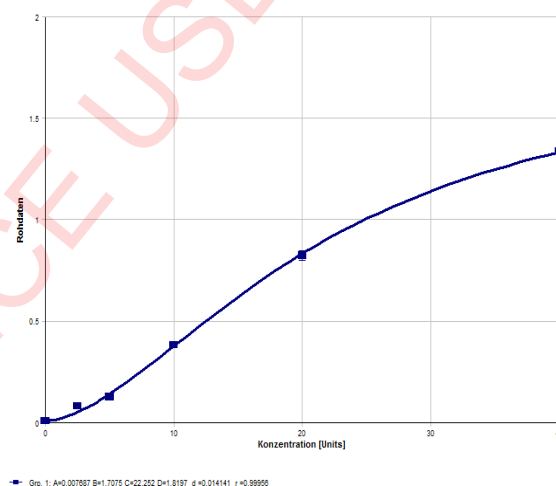
Sample 1 (14.5 U/ml): 3.9%

Interassay-Precision:

(n=10)

Sample 1 (13.8 U/ml): 3.6%

Typical Standard Curve



Assay Procedure

Incubation of samples/ standards/ controls	100 µl	2 hrs, 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-M5-AB, 1x96 determin.	15500	1100.-