



EIA for Quantitative Determination of anti-Muscarinic Cholinergic Receptor 1 (M1)-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. Autoantibodies against M1 AChR are significantly elevated in patients suffering from Lambert-Eaton myasthenic syndrome and from myasthenia gravis.

The CellTrend anti-muscarinic cholinergic receptor 1 (M1)-antibody EIA is designed for the determination of antibodies against the muscarinic cholinergic receptor 1 (M1) in serum.

Principle of the assay

The CellTrend muscarinic cholinergic receptor 1 (M1)-EIA is an antibody screening test. M1 receptor has been pre-coated onto a microtiter plate. During the first incubation the anti-muscarinic cholinergic receptor 1-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 1-antibody.

Performance Characteristics

Standard curve:

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

cut off: > 9.0 Units/ml positive

< 9.0 Units/ml negative

Sample materials:

Serum, Plasma

Intraassay-Precision:

(n=10)

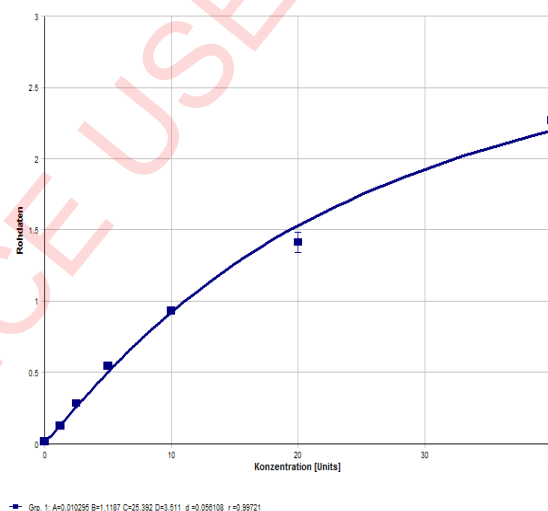
Sample 1 (17.2 U/ml): 4.1%

Interassay-Precision:

(n=10)

Sample 1 (16.5 U/ml): 8.6%

Typical Standard Curve



Assay Procedure

Incubation of samples/ standards/ controls	100 μ l	2 hrs, 4°C
Wash		
Incubation of detection antibody	100 μ l	1 hr, room temperature
Wash		
Substrate incubation	100 μ l/well	20 min, room temperature
Add Stopp solution	100 μ l/well	
Read at 450nm		

Order informations

Product	Catalog number	Price (€)
EIA for Quantitative Determination of anti-M1-AB, 1x96 determin.	15100	1100.-