



EIA for Quantitative Determination of anti-Muscarinic Cholinergic Receptor 2 (M2)-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 complex regional pain syndrome (Morbus Sudeck). The CellTrend anti-muscarinic cholinergic receptor 2 (M2)-antibody EIA is designed for the determination of antibodies against the muscarinic cholinergic receptor 2 (M2) in serum.

Principle of the assay

The CellTrend muscarinic cholinergic receptor 2 (M2)-EIA is an antibody screening test. M2 receptor has been pre-coated onto a microtiter plate. During the first incubation the anti-muscarinic cholinergic receptor 2-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 2-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)

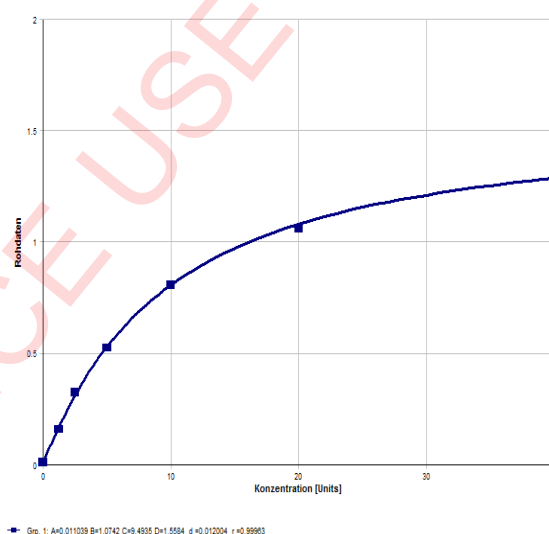
Sample 1 (19.5 U/ml): 9.9%

Interassay-Precision:

(n=10)

Sample 1 (17.8 U/ml): 15.2%

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-M2-AB, 1x96 determin.	15200	1100.-