

Human Myostatin ELISA kit ab267656

[1 References](#) [1 Image](#)

Overview

Product name Human Myostatin ELISA kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 12%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity 0.83 ng/ml

Range 0.819 ng/ml - 200 ng/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	101.9	86% - 123%
Plasma	89.4	67% - 122%
Cell culture media	102	99% - 107%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Human Myostatin ELISA Kit is designed for the quantitative determination of Myostatin in cell culture supernatants, plasma and serum samples.

This assay employs an antibody specific for human Myostatin coated on a 96-well plate. Standards and samples are pipetted into the wells and Myostatin present in a sample is bound to

the wells by the immobilized antibody. The wells are washed and biotinylated anti-human Myostatin antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of Myostatin bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

This ELISA kit shows no cross-reactivity with the following cytokines tested: Human ACE, Activin RIB (ALK-4), ADAM23, Armin, Cardiotrophin-1 (CT-1), Cathepsin V, FABP1 (LFABP), FGF-20, HAI-1, IL-27 R alpha (WSX-1), Insulin R (CD220), Kallikrein 7, LIF R alpha, Lipocalin-1, LTbR (Lymphotoxin beta R), Mesothelin, MFRP, Neuropilin-2, Neurturin, Nidogen-2, Olfactomedin-2 (Noelin-2), p53, PDECGF (Thymidine Phosphorylase), PDGF-CC, Progranulin, Ret, ROBO4, Semaphorin 6B, Serpin F1 (PEDF), SREC-I (SCARF1), SREC-II (SCARF2), TLR1, TLR3, TPP1 (Tripeptidyl-peptidase I), TREM-2, TrkC, TROY (TNFRSF19), Uromodulin, XIAP.

Platform Microplate

Properties

Storage instructions Store at -20°C. Please refer to protocols.

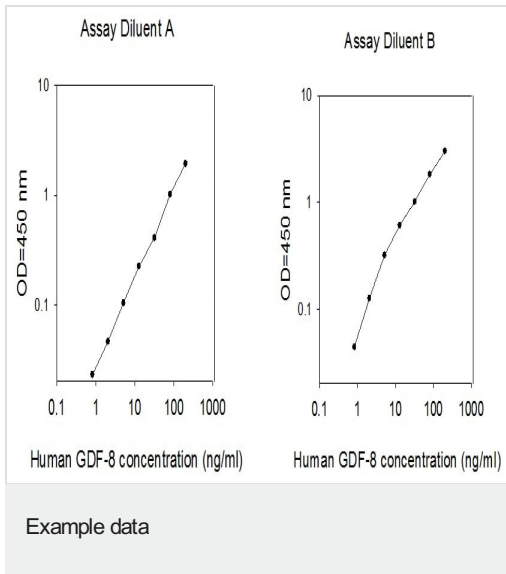
Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Anti-Human Myostatin coated Microplate (12 x 8 wells)	1 unit
Assay Diluent A	1 x 30ml
Biotinylated Anti-Human Myostatin Detection Antibody	2 vials
Human Myostatin Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

Function Acts specifically as a negative regulator of skeletal muscle growth.

Sequence similarities Belongs to the TGF-beta family.

Cellular localization Secreted.

Images



These standard curves are for demonstration only. A standard curve must be run with each assay

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors