

Recombinant Human MSTN

Recombinant human protein expressed in E. coli cells

Catalog # **G818-40N**

Lot # G3211-99

Product Description

Recombinant Human MSTN was expressed in E. coli cells. The protein accession number is [O14793](#)

Gene Aliases

MSTN, Growth differentiation factor 8, GDF-8, Recombinant Human Myostatin (MSTN)

Endotoxin Level

<1.0 EU/µg of recombinant protein as determined by the LAL method.

Formulation

Recombinant Myostatin was lyophilized from a 0.2 µm filtered Tris solution pH 8.0.

Reconstitution Protocol

A quick spin of the vial followed by reconstitution in distilled water to a concentration not less than 0.1 mg/mL. This solution can then be diluted into other buffers.

Storage and Stability

The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

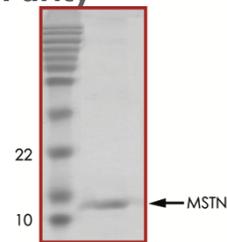
Scientific Background

Growth differentiation Factor 8 (GDF-8), also known as myostatin, is a secreted protein that is expressed specifically in developing and adult skeletal muscle. It controls myoblast proliferation and is a potent negative regulator of skeletal muscle mass. GDF-8 belongs to the transforming growth factor beta (TGF-beta) superfamily, which includes the TGF-betas, bone morphogenetic proteins (BMPs), growth differentiation factors (GDFs), activins, inhibins, leftys, nodal, Mullerian inhibitory substance (MIS) and the glial cell line-derived neurotrophic factors (GDNFs). All TGF-β superfamily members are synthesized and secreted as a homodimeric prepropeptide that is cleaved by proprotein convertases such as furin to generate the dimeric N-terminal propeptide and the dimeric C-terminal mature active protein.

Specific Activity

The ED(50) was determined by the dose-dependent proliferation inhibition of human MPC-11 cells was found to be in the range of 20-40 ng/mL.

Purity



Recombinant Human MSTN resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **13.0 kDa**

Recombinant Human MSTN

Recombinant human protein expressed in E. coli cells

Catalog Number	G818-40N
Specific Lot Number	G3211-99
Purity	95%
Stability	6 mos. at -20°C from date of shipment
Storage & Shipping	The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.