

Recombinant Human Thymosin-Beta 4 (TMSB4X)

Recombinant human protein expressed in E. coli cells

Catalog # T897-40N

Lot # G3223-17

Product Description

Recombinant Human Thymosin-Beta 4 (TMSB4X) was expressed in E. coli cells. The protein accession number is [P62328](#)

Gene Aliases

T beta-4, Hematopoietic system regulatory peptide, Seraspenide, TMSB4X, TB4X, TMSB4

Endotoxin Level

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

Formulation

Lyophilized from a 0.2 µm filtered solution in PBS

Reconstitution Protocol

Reconstitute at 100 µg/mL in sterile PBS

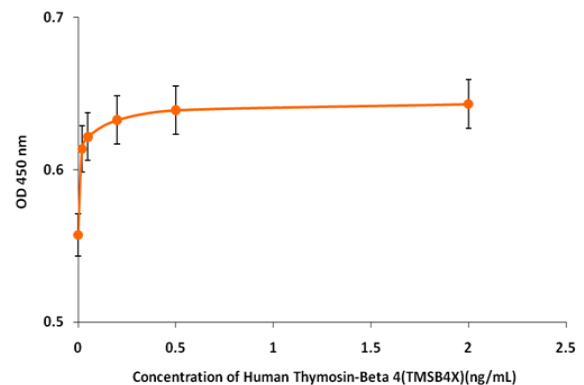
Storage and Stability

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

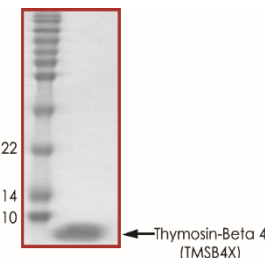
The primary function of the T beta-4 member of the Thymosin-beta family is to maintain the organization of the cell's cytoskeletal structure. Produced by myoblasts and fibroblasts, TMSB4X acts as a modulator for the production and digestion of actin and to inhibit filamentous actin polymerization. TMSB4X has also been found to induce phenotypic changes in T cell lines during early host defence mechanisms and inhibit the progression of hematopoietic pluripotent stem cells into the s-phase.

Specific Activity



The ED (50) was determined by the dose-dependent proliferation of immortalized mouse brain endothelial cells and was found to be <0.5ng/mL.

Purity



Recombinant Human Thymosin-Beta 4 (TMSB4X) resolved on a 17% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **5.2 kDa**

Recombinant Human Thymosin-Beta 4 (TMSB4X)

Recombinant human protein expressed in E. coli cells

Catalog Number	T897-40N
Specific Lot Number	G3223-17
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.