

Catalog #	Aliquot Size
S801-40N-10	10 µg
S801-40N-50	50 µg
S801-40N-1000	1 mg

Recombinant Human S100A4

Recombinant human protein expressed in E. coli cells

Catalog # S801-40N

Lot # G3223-10

Product Description

Recombinant Human S100A4 was expressed in E. coli cells. The protein accession number is [P26447](#)

Gene Aliases

Protein S100-A4, Calvasculin, Metastasin, Placental calcium-binding protein, Protein Mts1, S100 calcium-binding protein A4, CAPL, MTS1, Recombinant Human S100 Calcium Binding Protein A4 (S100A4)

Endotoxin Level

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

Formulation

Recombinant S100-A4 was lyophilized from a 0.2 µm filtered PBS solution and 1.0 mM DTT.

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.

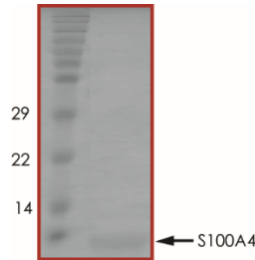
Storage and Stability

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

S100A4 is a calcium-binding protein related to the S100 family of proteins. It is known also as a growth-related protein, CAPL. It is expressed at high levels in different mouse metastatic cells and in normal tissues (thymus, spleen) and activated macrophages, T cells, neutrophils, monocytes, macrophages.

Purity



Recombinant Human S100A4 resolved on a 17% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **12.0 kDa**

Recombinant Human S100A4

Recombinant human protein expressed in E. coli cells

Catalog Number	S801-40N
Specific Lot Number	G3223-10
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.