

Catalog #	Aliquot Size
M838-40N-20	20 µg
M838-40N-100	100 µg
M838-40N-1000	1 mg

Recombinant Human Midkine (NEGF2)

Recombinant human protein expressed in E. coli cells

Catalog # M838-40N

Lot # G3211-94

Product Description

Recombinant Human Midkine (NEGF2) was expressed in E. coli cells. The protein accession number is [P21741](#)

Gene Aliases

MK, NEGF-2

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 0.1 mg/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.

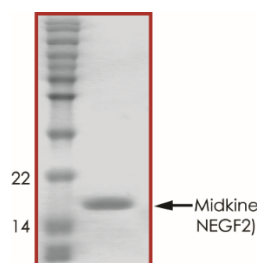
Scientific Background

Midkine and its functionally-related protein Pleiotrophin (PTN), are heparin-binding neurotrophic factors that signal through the same receptor. Midkine is involved in a number of biological processes including the migration of neutrophils, inflammatory leukocytes, macrophages, neuronal cells, and osteoclasts. Midkine is a basic, non-glycosylated polypeptide that contains five intra-chain disulfide bonds. There is 87% identity between the human and murine Midkine and approximately 50% identity between human Midkine and human PTN with the conservation of all 10 cysteines.

Specific Activity

The ED (50) was determined by the dose-dependent proliferation of AsPC-1 cells and was found to be <0.5ng/mL.

Purity



Recombinant Human Midkine (NEGF2) resolved on a 17% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **14.0 kDa**

Recombinant Human Midkine (NEGF2)

Recombinant human protein expressed in E. coli cells

Catalog Number	M838-40N
Specific Lot Number	G3211-94
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.