

## Recombinant Human GMFB

Recombinant human protein expressed in E. coli cells

**Catalog # G842-40N**

Lot # G3211-46

### Product Description

Recombinant Human GMFB was expressed in E. coli cells. The protein accession number is [P60983](#)

### Gene Aliases

GMF-beta, GMFB, Recombinant Human Glia Maturation Factor Beta (GMFB)

### Endotoxin Level

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

### Formulation

Recombinant GMFB was lyophilized from a 0.2 µm filtered 20 mM PB, 130 mM NaCl solution pH 7.5.

### 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.

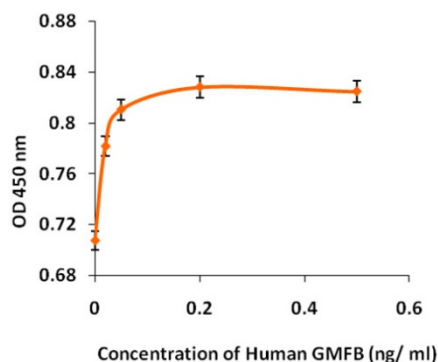
### Storage and Stability

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

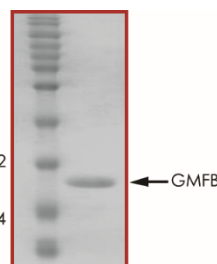
GMF-beta is a 17-kDa brain-specific protein that was isolated from bovine brain homogenate as a substance inducing the maturation of normal neurons as well as glial cells, and at first, it was considered to be a neurotrophic factor. The amino acid sequence of GMFB is highly conserved among many species, suggesting that it plays basic roles across many species. The expression of GMF-b is largely limited to the brain, especially the glial cells and some neurons. Schwann cells of the distal segment of the transected nerve express GMF-b, and this induction of GMF-b coincides with the temporal expression of nerve growth factor receptors in the cell.

### Specific Activity



ED50 not determined.

### Purity



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

Approx. MW **15.0 kDa**

## Recombinant Human GMFB

Recombinant human protein expressed in E. coli cells

Catalog Number	G842-40N
Specific Lot Number	G3211-46
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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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