

## Recombinant Human IGF2

Recombinant human protein expressed in E. coli cells

**Catalog # 1812-40N**

Lot # G3211-53

### Product Description

Recombinant Human IGF2 was expressed in E. coli cells. The protein accession number is [P01344](#)

### Gene Aliases

Insulin-like growth factor II, multiplication stimulating activity, MSA, Somatomedin-A, T3M-11-derived growth factor, PP1446

### 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 Tris and NaCl

### Reconstitution Protocol

Reconstitute in sterile PBS at 100 µg/mL

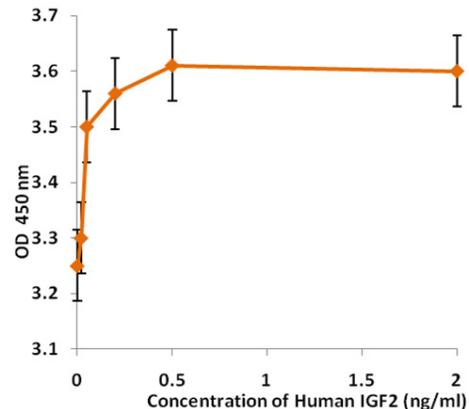
### Storage and Stability

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

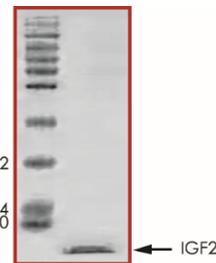
Insulin-like growth factor 2 (IGF2) is one of the mitogenic, polypeptide growth factors that belongs to the insulin family, due to its structural and functional similarity to insulin. IGF2 is mainly produced in the liver and several other tissues at certain times. IGF2 expression can be found in many tissues usually with autocrine, paracrine and endocrine functions. IGF2 mainly acts as a growth promoting hormone during gestation, and does so by binding to the tyrosine kinase type I receptor (IGF1R) as well as the IGF2 receptor cation-independent mannose 6-phosphate receptor. Mature IGF2 is highly conserved between human, bovine, and porcine species. Recombinant Human IGF2 is a single, non-glycosylated, polypeptide chain of 7.8 kDa.

### Specific Activity



The ED50 as determined by dose dependent proliferation of MCF7 cells was <1 ng/ml

### Purity



Recombinant Human IGF2 resolved on a 17% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **7.8 kDa**

## Recombinant Human IGF2

Recombinant human protein expressed in E. coli cells

Catalog Number	1812-40N
Specific Lot Number	G3211-53
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.**