

Recombinant Human EGF

Recombinant human protein expressed in E. coli cells

Catalog # E801-40N

Lot # G3211-26

Product Description

Recombinant Human EGF was expressed in E. coli cells. The protein accession number is [P01133](#)

Gene Aliases

Urogastrone, EGF, Recombinant Human Epidermal Growth Factor (EGF)

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 buffer with NaCl

Reconstitution Protocol

A quick spin of the vial followed by reconstitution in sterile distilled water to a concentration not less than 0.1 mg/mL. This solution can then be diluted into other buffers.

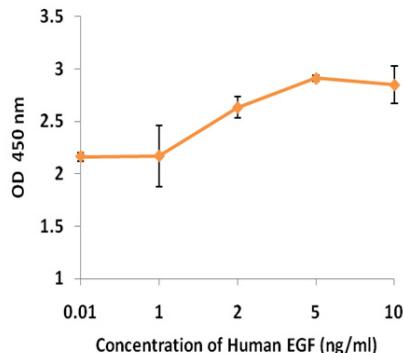
Storage and Stability

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

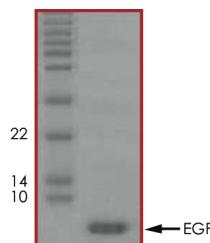
Epidermal growth factor (EGF) is a small mitogenic protein that is thought to be involved in mechanisms such as normal cell growth, inhibition of gastric acid secretion, oncogenesis and wound healing. EGF shares strong sequential and functional homology with the type-alpha transforming growth factor, which is a competitor for EGF receptor sites. EGF signals through a receptor known as c-erbB, which is a class I tyrosine kinase receptor. The EGF receptor is expressed in almost all types of tissues with the exception of the parietal endoderm, mature skeletal muscles and hematopoietic tissue. Recombinant Human EGF is a globular protein that displays the EGF-like domain; characterized by three intra-molecular disulfide bonds that are formed by six similarly spaced conserved cysteine residues. Mature human EGF shares 70% amino acid sequence identity with its counterparts from mouse and rat.

Specific Activity



The ED₅₀ as determined by a cell proliferation assay using 3T3 cells was <1ng/ml.

Purity



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

Approx. MW **6.5 kDa**

Recombinant Human EGF

Recombinant human protein expressed in E. coli cells

Catalog Number E801-40N

Specific Lot Number G3211-26

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.