



6405 Mira Mesa Blvd, Suite 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Human Keratinocyte Growth Factor Recombinant

Catalog #: 90205-B
Lot #: 230501

Description: Recombinant human Keratinocyte Growth Factor is a disulfide-linked monomeric protein consisting of 164 amino acid residues. Native human KGF is generated by the proteolytic removal of the signal peptide and propeptide. KGF migrates as an approximately 19 kDa protein under non-reducing conditions and reducing conditions in SDS-PAGE.

Source: Optimized DNA sequence encoding Human Keratinocyte Growth Factor mature chain was expressed in *E. coli*.

Formulation: Lyophilized from a 0.2 μ m filtered 20 mM Phosphate buffer, 1M NaCl, pH 8.0.

Reconstitution: 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: The lyophilized protein is stable for at least 2 years from date of receipt when stored at -20°C. Upon reconstitution, store in working aliquots at 2 - 8° C for up to one month, or at -20°C for up to six months, in the presence of a carrier protein. Avoid repeated freeze/thaw cycles.

Purity: >96%, as determined by SDS-PAGE and HPLC.

Endotoxin Level: Endotoxin level was found to be < 0.1 ng/ μ g (1EU/ μ g), using the LAL gel clot method.

Biological Activity: The ED(50) was determined by the dose-dependent proliferation of BAF3 cells was found to be less than 10 ng/ml.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.
To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: support@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com