



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

BCA1 (CXCL13)

Catalog #: 90104-B (20 µg)
Lot #: 150723

Description: Recombinant human BCA1 (B cell-attracting chemokine 1), also known as CXCL13 or BLC, is a disulfide-linked monomeric protein consisting of 88 amino acids and migrates as an approximately 10 kDa protein under reducing and non-reducing conditions. Human CXCL13, generated by the proteolytic removal of the signal peptide and propeptide, has a calculated molecular mass of ~10 kDa.

Source: Optimized DNA sequence encoding human BCA1 mature domain was expressed in *E. coli*.

Formulation: Lyophilized from a 0.2 µm filtered solution containing 20 mM phosphate buffer, pH 7.5, 100 mM NaCl.

Reconstitution: Reconstitute at 0.1-1.0 mg/mL in distilled water. This solution can then be diluted into other buffers. To maximize product collection from vial surface, vortex briefly and then spin down to recollect the liquid.

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: >98%, as determined by SDS-PAGE and HPLC.

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

Biological Activity: Determined by its ability to chemoattract human CXCR5 transfected mouse BaF3 B cells with ED₅₀ >20 ng/ml.

Protein Sequence:

VLEVYYTSLRCRCVQESSVFIPRRFID
RIQILPRGNGCPRKEIIVWKKKNKSIVCV
DPQAEWIQRMMEVLRKRSSSTLPVPV
FKRKIP

References:

1. *Neurology*, Dec 2009; **73**: 2003 - 2010.
2. *Nephrol. Dial. Transplant.*, Dec 2009; **24**: 3708 - 3712.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com