



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

## ***Human BCA1(CXCL13) Recombinant***

**Catalog #:** 90104-B (20 µg)  
**Lot #:** 221017

**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:** 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:** >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  
RIQILPRGNGCPRKEIIVWKKNKSIKCV  
DPQAEWIQRMMEVLRKRSSSTLPVPV  
FKRKIP

### **References:**

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