



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

## ***Human Thrombopoietin Recombinant***

Catalog #: 90249-B  
Lot#: 221021

**Description:** Recombinant human Thrombopoietin erythropoietin, also known as TPO, is a disulfide-linked monomeric protein consisting of 332 amino acids and migrates as an approximately 35 kDa protein under reducing and non-reducing conditions.

**Source:** Optimized DNA sequence encoding human TPO erythropoietin domain was expressed in Chinese Hamster Ovary Cells.

**Formulation:** Lyophilized from a 0.2  $\mu$ m filtered solution in PBS, pH 7.

**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/ $\mu$ g (1EU/ $\mu$ g), using the LAL gel clot method.

**Biological Activity:** The ED50, was determined by the dose-dependent proliferation of Mo7e cells, was found to be in the range of 1.0-2.0 ng/ml.

### **References:**

1. Blood, Aug 2009; 10.1182/blood-2009-04-219493.
2. J. Biol. Chem., May 2009; 284: 11781-1179