



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 Erythropoietin-alpha Recombinant

Catalog #: 90257
Lot#: 231011

Description: Recombinant human EPO-a is a glycosylated monomer protein consisting of 166 amino acids and, due to glycosylation migrates as an approximately 35 kDa protein under reducing conditions.

Source: Optimized DNA sequence encoding Human EPO-a mature chain was expressed in CHO cells.

Formulation: Lyophilized from a 0.2 μ m filtered PBS solution pH 7.0.

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: >98%, 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: Determined by the dose-dependent proliferation assay using a factor-dependent human erythroleukemic cell line TF-1, activity was found to be 0.8 x 10⁵ IU/mg.

References:

1. *J. Biol. Chem.*, Sep 2009; **284**: 26988 - 26998.
2. *Invest. Ophthalmol. Vis. Sci.*, Jul 2009; **10.1167**/iovs.09-3544.

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