



Certificate of Analysis

Human Erythropoietin (EPO), ultrapure

Catalog No: 11965-1

Lot No:

Size: 500 units/vial

Description: Recombinant Erythropoietin (EPO), ultrapure

Volume: ml

Concentration: units/ml

Specific Activity: x 10 units/mg

Buffer: Phosphate buffered saline containing 0.1% bovine serum albumin (BSA)

Endotoxin: < 1 EU/ g

Molecular Weight: 34-36 kDa

Purity: > 95%

Purification Method: Proprietary chromatography method

Source: Human EPO cDNA expressed in CHO cells

Assay Used to Measure Bioactivity: Biological assay in the model of the ex-hypoxic polycythemic mice versus international standard (EPO NIBSC World Health Organization (WHO) #87/684).

Shipping Conditions: Dry Ice

Physical State of Product During Shipping: Frozen

Storage Conditions/Comments: After receipt, the product may be stored at -20°C for short-term use (6 months). For long-term storage, we recommend storing the product at -70°C or below for retention of full activity. When thawing, the contents of the tube should be apportioned in separate tubes so that freezing and thawing is kept to a minimum. Re-freezing should be done on dry ice or in a dry ice/alcohol bath. Further dilution of the product should be in buffers containing protein such as 0.1% bovine serum albumin (BSA). For more information on protein handling, visit our Resource Library at www.pbl assaysci.com.

Authorization

Released by: _____

Date:

Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a PBL Assay Science. For research use only. Not for diagnostic or clinical use in, or administration to, humans. Not for resale in original or any modified form, including inclusion in a kit, for any purpose. Not for use in the preparation of any commercial product.

Rev. 03