

## EPO

Recombinant human protein expressed CHO cells

**Catalog # E860-40N**

Lot # G3211-28

### Product Description

Recombinant Human EPO was expressed in CHO cells. The protein accession number is [P01588](#)

### Gene Aliases

Erythropoietin  
Recombinant Human Erythropoietin-Alpha (EPO)

### Endotoxin Level

<1.0 EU/µg of recombinant protein as determined by the LAL method

### Formulation

Lyophilized from a 0.2 µm filtered solution in PBS (pH 7)

### Reconstitution Protocol

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.

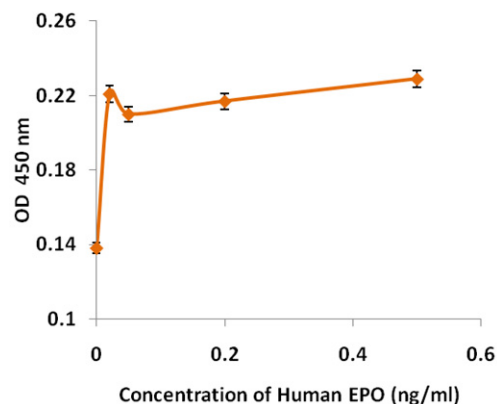
### Storage and Stability

The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

### Scientific Background

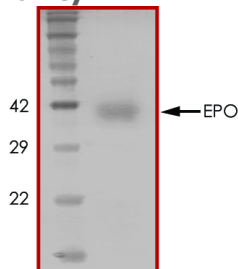
EPO is predominantly synthesized and secreted by tubular and juxtatubular capillary, endothelial, and interstitial cells of the kidney. Approximately 10-15% of the total amount of EPO comes from extrarenal sources and is predominantly produced by hepatocytes and Kupffer cells of the liver. Approximately 40% of the molecular mass of EPO is due to its glycosylation. Glycosylation is an important factor determining the pharmacokinetic behaviour of EPO in vivo. Non-glycosylated Epo has an extremely short biological half life. Recombinant Human EPO is a glycosylated protein that runs at approximately 35 kDa owing to its glycosylation.

### Specific Activity



Activity was determined by the dose-dependent proliferation assay using a factor-dependent human erythroleukemic cell line TF-1 and was found to be less than 0.2ng/ml

### Purity



Recombinant Human EPO resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **20.0 kDa**

## EPO

Recombinant human protein expressed in CHO cells

Catalog Number	E860-40N
Specific Lot Number	G3211-28
Purity	95%
Stability	6 mos. at -20°C from date of shipment
Storage & Shipping	The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**

# MATERIAL SAFETY DATA SHEET

---

## Article 1 - Product Identification and Use

---

**Product Name: Recombinant Human EPO**

**Catalog # E860-40N**

*This product is sold only for research use by qualified laboratory personnel, and is not to be used as a drug, medical device, food additive, cosmetic, nor household chemical. It is not to be used in diagnostic, therapeutic, consumer, agricultural, nor pesticidal applications.*

Manufacturer's Name:	SignalChem Pharmaceuticals Inc.
Street Address:	110-13120 Vanier Place
City, Prov. Postal Code:	Richmond, BC, V6V 2J2
Fax:	604-232-4601
EMERGENCY PHONE:	604-232-4600

---

## Article 2 - Hazardous Ingredients

---

NOT AVAILABLE. We are not aware of any hazards associated with this product or its ingredients, but the chemical, physical, and toxicological properties of this product have not been investigated thoroughly. Observe normal laboratory precautions.

---

## Article 3 - Physical Data

---

This product consists of purified protein in PBS buffer shipped on dry ice. The physical properties of this product have not been investigated thoroughly.

---

## Article 4 - Fire and Explosion Hazard

---

NOT APPLICABLE

---

## Article 5 - Reactivity Data

---

NOT APPLICABLE

---

## Article 6 - Toxicologically Data

---

May be harmful by inhalation, ingestion, or skin absorption. The toxicological properties of this product have not been investigated thoroughly. Exercise due caution.

---

## Article 7 - Preventative Measures

---

Wear chemical safety goggles and compatible chemical-resistant gloves. Avoid inhalation, contact with eyes, skin or clothing.

\*\*\*\*\*MULTIPLE COMPONENT SPILL OR LEAK PROCEDURES\*\*\*\*\*

- Wear protective equipment.
- Absorb on sand or vermiculite and place in closed containers for disposal.
- Observe all federal, state and local environmental regulations.

---

## Article 8 - First Aid Measures

---

- If swallowed, wash out mouth with water, provided person is conscious. Call a physician.
- In case of skin contact, flush with copious amounts of water for at least 15 minutes. Remove contaminated clothing and shoes. If a rash or other irritation develops, call a physician.
- If inhaled, remove to fresh air. If breathing becomes difficult, call a physician.
- In case of eye contact, flush with copious amounts of water for at least 15 minutes while separating the eyelids with fingers. Call a physician.

---

## Article 9 - Preparation

---

Prepared by: Jun Yan

Phone#: 1-866-954-6273

The above information is believed to be correct but does not purport to be all-inclusive and shall be used only as a guide. SignalChem shall not be held liable for any damage resulting from handling or from contact with the above product. See the Technical Specification, Packing Slip, Invoice, and Product Catalogue for additional terms and conditions of sale.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**