

# Erythropoietin $\alpha$ (human), (recombinant)

**Recombinant human EPO that stimulates red blood cell production and increases oxygen capacity.**

Erythropoietin (EPO) is a 34 kDa glycoprotein hormone produced by the kidney that promotes the formation of red blood cells in the bone marrow, thereby increasing the oxygen carrying capacity of blood in times of oxygen demand.

**Manuals, SDS & CofA**

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name EPO

Application Notes Western blot control.

Formulation Lyophilized. Each milligram contains 5.8mg sodium chloride and 0.06mg citric acid. Reconstitute with sterile 18W-cm water not <100µg/mL

MW Predicted MW of ~21kDa (single chain) and ~30kDa (glycosylated)

Purity ≥98% (SDS-PAGE; RP-HPLC)

Purity Detail Purified by multi-step chromatography.

Source Produced in CHO cells.

UniProt ID P01588

Last modified: October 9, 2025



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)