

## Erythropoietin human

Erythropoietin (EPO) is a glycoprotein renal hormone responsible for promoting erythropoiesis, formation of red blood cells by bone marrow. Oxygen scarcity or bleeding stimulate release of EPO. It can be used to test for and treat certain blood conditions like anemia, and has been misused by athletes as a performance-enhancer.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1224</b>
<b>Quantity Supplied:</b>	1 vial(s), 10 µg per vial
<b>Reconstitution Buffer:</b>	1 mL

### Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

### Product Information

<b>Molecular Weight:</b>	36 kg/mol
<b>Source:</b>	Recombinant
<b>Purity:</b>	≥ 95% SDS PAGE
<b>Endotoxin Level:</b>	< 1EU/µg
<b>Formulation:</b>	N/A
<b>Storage:</b>	-20°C Please avoid multiple freeze/thaw cycles.

### Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

<b>Stock Concentration:</b>	100 µg/mL
<b>Reconstitution Volume:</b>	0.100 mL
<b>Reconstitution Solvent:</b>	Reconstitution Buffer

***For Research Use Only***

Ordering: +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
Technical support: +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
General product information: [www.discoverx.com](http://www.discoverx.com)