

Human Erythropoietin (EPO), ultrapure

Ordering Information

[Order Online »](#)

PBL-11965-1

500U

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	EPO
Endotoxin Content	< 1 EU/μg
Formulation	Liquid. In PBS containing 0.1% BSA.
MW	~34.0-36.0kDa
Purity	≥95%
Purity Detail	Proprietary chromatography methods
Quantity	0.1ml
Source	Produced in CHO cells.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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