

Data Sheet

EPOR, Fc fusion, Avi-Tag,

Biotin-labeled

Human, Recombinant, C-terminal Fc-Avi-tag,
Biotin-labeled.

Catalog #: 100612

Lot #: 200130

Conc.: 1.03 mg/ml

Formulated in: 8 mM Phosphate, pH 7.4,
110 mM NaCl, 2.2 mM KCl, and 20%
glycerol.

Stability: At least 6 months at -80°C . *Avoid
freeze/thaw cycles.*

References:

1. Pinto, J.P., *et al. Blood*. 2008; **111(12)**:
5727-5733.
2. Lindern, M.V, *et al., Oncogene*. 2001;
20: 3651-3664.

Description: Human Erythropoietin Receptor, also known as EPOR, GenBank Accession No. NM_000121, a.a 25-250, fused at the C-terminus of the Fc portion of human IgG1, with C-terminal Avi-tag, expressed in a HEK293 cell expression system and enzymatically biotin-labeled using Avi-tag™ technology. Biotinylation is confirmed to be $\geq 90\%$. MW=54 kDa.

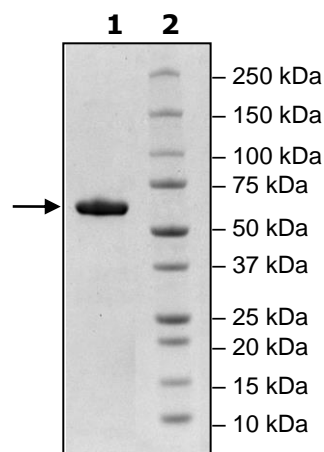
Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
4 μg EPOR
Lane 2:
Protein Marker

MW: 54 kDa

Purity: $\geq 90\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

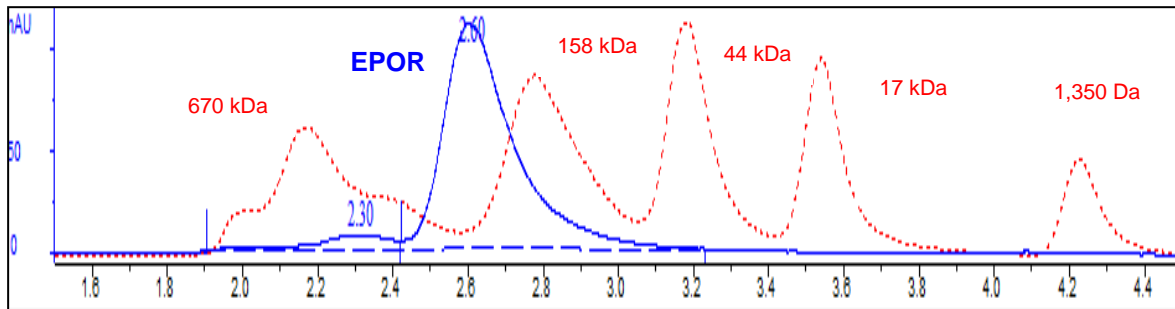
To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com



<10% Aggregation

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com