

Data Sheet

EPOR, Fc fusion, Avi-Tag

Human, Recombinant, C-terminal Fc-Avi-tag

Catalog #: 100611

Lot #: 200115

Conc.: 0.46 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, MW=54 kDa.

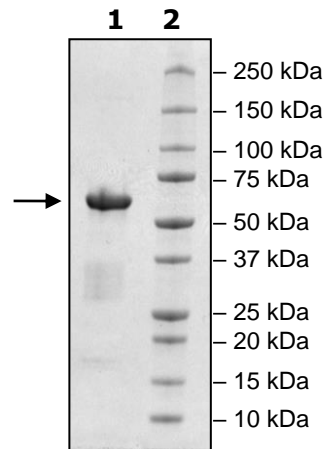
Quality Assurance

4-20% SDS-PAGE **Coomassie staining**

Lane 1:
3µg EPOR
Lane 2:
Protein Marker

MW: 54 kDa

Purity: ≥82%



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