

Data Sheet

EPOR-(GSSG)₆- JAK2(V617F, Y1007F, Y1008F)

Human, Recombinant

Catalog #: 100713**Lot #:** 200623-GA**Conc.:** 2.22 mg/ml**Formulated in:** 50 mM Tris-HCl pH 8.0, 500 mM NaCl, 10% glycerol, 3 mM DTT.**Stability:** At least 6 months at -80°C. Avoid freeze/thaw cycles.**References:**

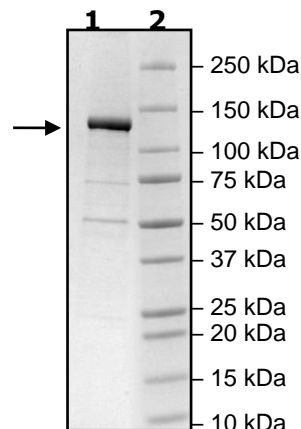
1. Tono, C., *et al.*, *Leukemia*. 2005; **19(10)**: 1843-1844.
2. Goertler, P.S., *et al.*, *Blood*. 2005; **106(8)**: 2862-2864.

Description: Complex of Human **EPOR**, also known as Erythropoietin receptor, GenBank Accession No. NM_000121, a.a. 273-338 with N-terminal His-tag, MW=9 kDa, and **JAK2**, also known as Janus Kinase 2, with V617F, Y1007F and Y1008F mutations, GenBank Accession No. NM_004972, a.a. 36-1132(end), with N-terminal FLAG-tag, MW=128 kDa. The proteins were expressed as single fusion protein with a (GSSG)₆ linker in a Baculovirus infected Sf9 cell expression system.

Quality Assurance

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

Lane 1:
2 µg EPOR-
(GSSG)₆-JAK2
Lane 2:
Protein Marker

MW: 139 kDa**Purity:** 89%

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.comPlease visit our website at: www.bpsbioscience.com