

Data Sheet

Zebov GP1

Zaire Ebola Virus, Recombinant GP1,
C-terminal His-tag

Catalog #: 21002

Lot #: 141118-D

Conc: 1.11 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. Storing diluted protein is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

References:

1. Lee, J.E., *et al.*, *Nature*. 2008; **454(7201)**: 177-182.
2. Dube, D., *et al.*, *J. Virol.* 2009; **83(7)**: 2883-2891.

Description: Zaire Ebola Virus (Zebov) glycoprotein 1 (GP1), also known as envelope glycoprotein1, with a C-terminal His-tag, GenBank Accession No. NP_066246, MW=57.1 kDa, expressed in a HEK293 cell expression system. This protein runs at a higher apparent molecular weight due to glycosylation.

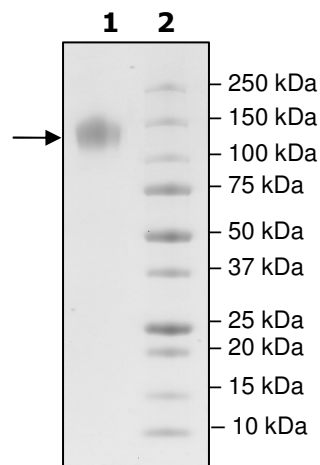
Application: Useful for protein binding studies, screening inhibitors of protein-protein interactions, and studying protein function.

Quality Assurance

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

Lane 1:
4 μg Zebov GP1
Lane 2:
Protein Marker

MW: 57.1 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 **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com