

Data Sheet

Zebov GP1,2, Biotin-labeled

Zaire Ebola Virus, Recombinant GP1,2, C-terminal

Avi-biotin, His-tag

Catalog #: 21004

Lot #: 141125-BD

Conc: 2.52 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 1,2 (GP1,2), also known as envelope glycoprotein, and GP, a.a. 33-650 with a C-terminal Avi and His-tag, GenBank Accession No. NP_066246, MW = 70.6 kDa, expressed in a HEK293 cell expression system. The protein was enzymatically biotinylated using AviTag™ technology. 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

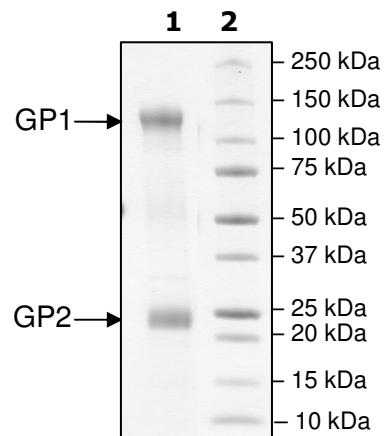
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
3 μg Zebov GP1,2

Lane 2:
Protein Marker

MW: 70.6 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