

## Data Sheet

### Zebov GP1,2

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

**Catalog #:** 21003

**Lot #:** 141118-D

**Conc:** 1.32 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  $-80^{\circ}\text{C}$ . *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 with a C-terminal His-tag, a.a. 33-650, GenBank Accession No. NP\_066246, MW=68.7 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.

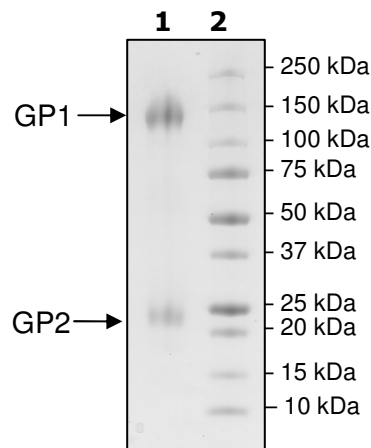
### Quality Assurance

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

**Lane 1:**  
7  $\mu\text{g}$  Zebov GP1,2

**Lane 2:**  
Protein Marker

**MW:** 68.7 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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)