

## Data Sheet

### **MICA, Fc fusion, Avi-Tag, Biotin-labeled**

Human, recombinant, C-terminal Fc-Avi-Tag,  
Biotin-Labeled

**Catalog #:** 100300

**Lot #:** 190307 **Conc.:** 2.53 mg/ml

**Formulation:** 8 mM Phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

**Stability:** Stable for at least 12 months at -80°C. Avoid freeze/thaw cycles.

**Description:** Human MICA, also known as, MHC Class I Polypeptide-Related Sequence A, PERB11.1, and Stress Inducible Class I Homolog, GenBank Accession No. NM\_001177519, a.a. 23-308, fused at the C-terminus to the Fc portion of human IgG1 followed by a C-terminal Avi-Tag™ and expressed in a HEK293 cell expression system. MW=62 kDa. This protein is enzymatically biotinylated using Avi-Tag technology.

### **References:**

1. Romphruk, A.V., *et al.*, *HLA*. 2001. Aug; **58(2)**:83-89.
2. Gaudieri, S., *et al.*, *Immuno-genetics*. 1997. Jan; **45(3)**: 209-216

---

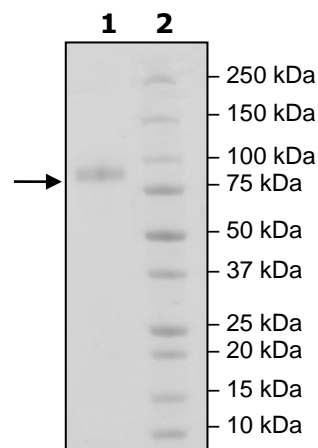
## **Quality Assurance**

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

**Lane 1:**  
4µg MICA

**Lane 2:**  
Protein Marker

**MW:** 62 kDa  
**Purity:** ≥90%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** 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)