

Data Sheet

MICA, Fc fusion, Avi-Tag

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

Catalog #: 100295

Lot #: 190306

Conc.: 1.3 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.

References:

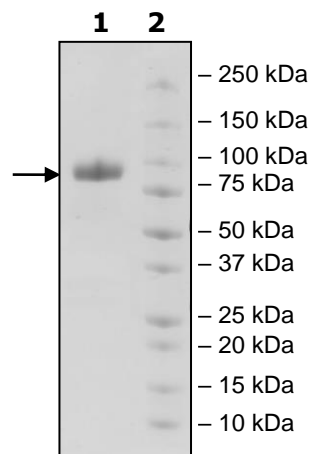
1. Romphruk, A.V., *et al.*, *HLA*. 2001. Aug; **58(2)**:83-89.
2. Gaudieri, S., *et al.*, *Immuno-genetics*. 1997. Jan; **45(3)**: 209-216

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
Please visit our website at: www.bpsbioscience.com