

Data Sheet

FcRn (FCGRT/B2M), His-Avi-Tag, Biotin-Labeled

Mouse, Recombinant, C-terminal His-Avi-tag

Catalog #: 71286

Lot #: 161215 **Conc.:** 1.85 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 enzyme is not recommended, if necessary use carrier protein (BSA 0.1 – 0.5%).

References:

1. Roopenian, D., *et al.*, *Nature*. 2007 Sep 7: 715-725
2. Shields, R. *et al.*, *Biological Chemistry*. 2000 Nov; **276**:6591-6604.

Description: Mouse FcRn also known as Neonatal Fc receptor is a heterodimer of FCGRT (Fc Fragment of IgG Receptor and Transporter) and B2M (beta 2 microglobulin), GenBank Accession numbers NM_010189 and NM_009735; a.a. 22-297 and a.a. 1-119 respectively. FCGRT has a C-terminal His-Avi-tag. This protein is enzymatically biotinylated using AviTag™ technology. These proteins are co-expressed in a HEK293 expression system. FcRn MW = 34 kDa, B2M MW = 14 kDa.

Application: Useful as a positive control for western blot analysis.

Quality Assurance

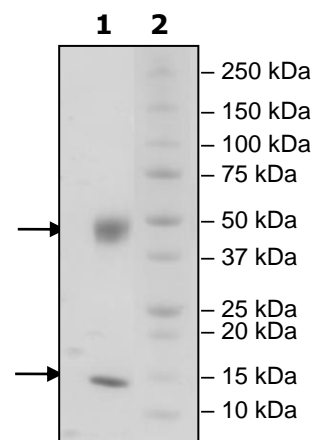
4-20% SDS-PAGE Coomassie staining

Lane 1:
5µg FcRn/B2M

Lane 2:
Protein Marker

MW:
FcRn: 34 kDa
B2M: 14 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 **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com