

Data Sheet

FcRn (FCGRT/B2M), His-Tag (Mouse)

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

Catalog #: 11349

Lot #: 180517

Conc.: 3.0 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. 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

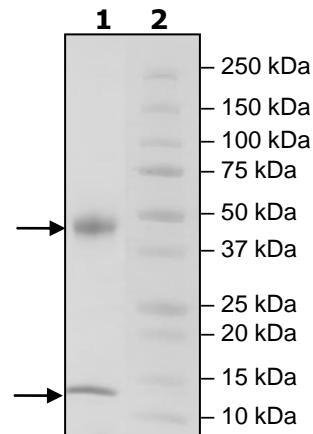
4-20% SDS-PAGE Coomassie staining

Lane 1:
6 μg FcRn/B2M

Lane 2:
Protein Marker

MW:
FcRn: 34 kDa
B2M: ~14 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**

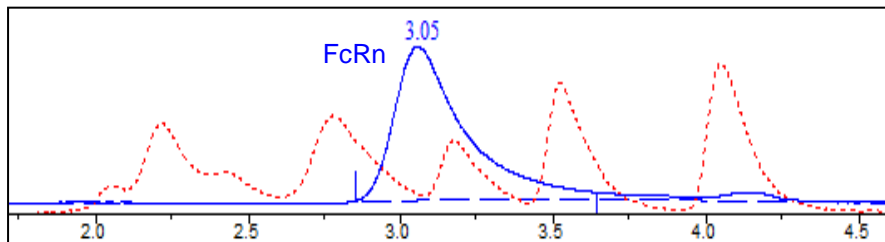
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Gel Filtration Trace



<10% Aggregation

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