

Data Sheet

FcRn (FCGRT/B2M), His-Tag

Human, Recombinant, C-terminal His-tag

Catalog #: 71285

Lot #: 190702 **Conc.:** 2.26 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.

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: Human complex of FcRn also known as Neonatal Fc receptor, heterodimer of FCGRT (Fc Fragment of IgG Receptor and Transporter), GenBank Accession No. NM_001136019, a.a. 24-297, with a C-terminal His-tag, MW=32 kDa and B2M, also known as beta 2 microglobulin, GenBank Accession No. NM_004048, a.a. 1-119, MW=14 kDa. Individually expressed in a HEK293 expression system.

Quality Assurance

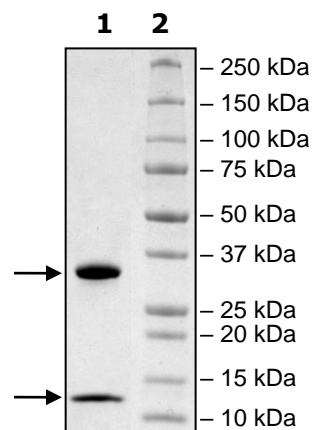
4-20% SDS-PAGE Coomassie staining

Lane 1:
4 μg FcRN/B2M

Lane 2:
Protein Marker

MW:
FcRN: 32 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 **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com