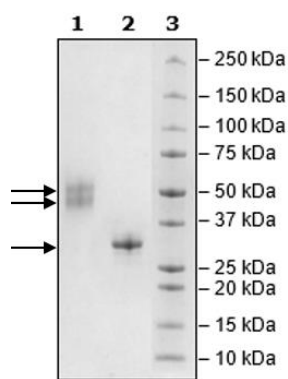


Product Information

Description:	The FcRn (FCGRT/B2M) blocker (Fc-Tag) is a purified human protein corresponding to the Fc region of an antibody which has been engineered to bind to human neonatal Fc Receptor for IgG (FcRn; FCGRT/B2M) with high affinity. This antibody has been tested for specific binding to human FcRn (BPS Bioscience #71283) and it blocks the binding of FcRn with immunoglobulins G.
Concentration:	0.32 mg/ml
Species:	Human
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
Expression System:	HEK293
Purification:	Protein A affinity
Format:	Aqueous buffer solution
Stability:	At least 12 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
MW:	51.2 kDa (homo-dimer); Chain: 25.6 kDa
Purity:	≥90%
Assay Conditions:	Assay 1. A 96-well plate was coated overnight at 4°C with 50 µl/well of the FcRn blocker at a concentration of 2 µg/ml, diluted in PBS. The plate was washed 3 times with Immuno Buffer 1 (BPS Bioscience #79311) and blocked with 100 µl of Blocking Buffer 6 for an hour at room temperature. A serial dilution of biotinylated FcRn protein (BPS Bioscience #71283) prepared in Blocking Buffer 6 was added for 30 minutes. After 3 washes, the plate was incubated with Streptavidin-HRP followed by more washes and addition of a colorimetric HRP substrate. After quenching, color intensity was measured at 450 nm. Assay 2. The effect of FcRn (FCGRT/B2M) Blocker was assessed using the Fc(IgG1):FcRn Inhibitor Screening colorimetric kit (BPS Bioscience #78501) using the following conditions: Fc(IgG1) protein at 2 µg/ml, diluted in PBS, serial dilution of FcRn Blocker prepared in blocking buffer, and 1.5 µg/ml of FcRn-Biotin.
Applications:	This product is for research use only. It is not suitable for human diagnostic or therapeutic use.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

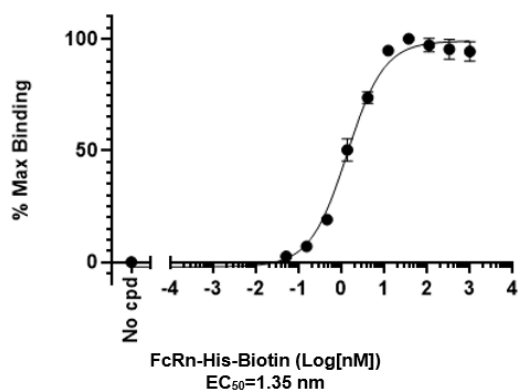


Lane 1: Anti-FcRn (Non-reduced)

Lane 2: Anti-FcRn (Reduced)

Lane 3: Protein Marker

Assay 1. FcRn blocker:FcRn-Biotin Binding Assay



Assay 2. Fc(IgG1):FcRn-Biotin Neutralization

