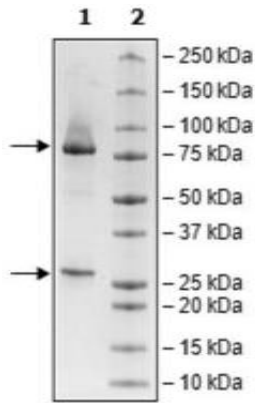


Product Information

Description:	Anti-FcGR3A-anti-EGFR bispecific IgG antibody, which recognizes FcGR3A (also known as CD16A) and EGFR antigen binding proteins. The amino acid sequences correspond to the sequences from the bispecific EGFR/FcGR3 antigen-binding protein patent WO2019175368A1. This bispecific antibody has been tested for specific binding affinity to purified human FcGR3A and EGFR recombinant protein.
Species:	Human
Clonality:	Comprised of two monoclonal antibodies
Concentration:	3.14 mg/ml
Expression System:	HEK293
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	Heavy Chain: 76 kDa; Light Chain: 23 kDa + glycans
Glycosylation:	This antibody runs at a higher MW by SDS-PAGE due to glycosylation.
Stability:	At least 12 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Assay Conditions:	Functional validation: The antibody was validated by measuring anti-FcGR3A-anti-EGFR binding to FcGR3A and EGFR antigens in ELISA assay. The FcGR3A protein (BPS Bioscience #79013) and EGFR protein were coated onto a 96-well plate overnight at 4°C (50 µl/well at a concentration of 2 µg/ml in PBS). The plate was washed 3 times with Immuno Buffer 1 (BPS Bioscience #79311) and blocked using 100 µl of Blocking Buffer 2 (BPS Bioscience #79728) for 1 hour at room temperature. After removing the blocking buffer, 50 µl/well of purified anti-FcGR3A-anti-EGFR antibody (BPS Bioscience #101535), serially diluted in Blocking Buffer 2, was added for 30 minutes at room temperature. After 3 more washes, the plate was incubated with Goat Anti-Human IgG Fc (HRP) (Abcam #ab97225), washed, and incubated with the Colorimetric HRP substrate. The reaction was stopped, and absorbance was read at 450 nm. The Blank value was subtracted from all values.
Applications:	Useful for studying the binding of FcGR3A and EGFR in ELISA and in cellular assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Binding Assay

