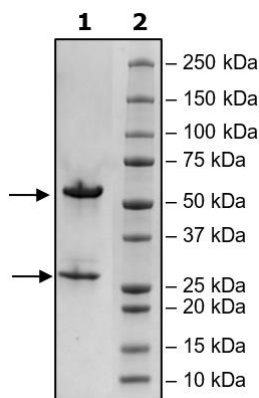


Product Information

Description:	Anti-Claudin-18 isoform 2 (CLDN18.2) IgG1 monoclonal antibody is a purified recombinant antibody which recognizes human CLDN18.2 antigen. This antibody has been tested for specific binding affinity to purified human CLDN18.2 protein.
Concentration:	1.31 mg/ml
Species:	Human
Clonality:	Monoclonal
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:	Total: 150 kDa; Light Chain: 24 kDa; Heavy Chain: 51 kDa
Purity:	≥90%
Assay Conditions:	Functional validation: The antibody was validated by measuring its binding to CLDN18.2 antigen in ELISA. The CLDN18.2 protein was 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) with DDM/CHS detergent (dodecyl-β-D-maltoside (DDM) 0.05%, cholesteryl hemisuccinate (CHS) 0.01%) for 1 hour at room temperature. After removing the blocking buffer, 50 µl/well of purified anti-CLDN18.2 antibody (BPS Bioscience #101564), serially diluted in Blocking Buffer 2 with DDM/CHS detergent, was added for 60 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 Colorimetric HRP substrate (BPS Bioscience #79651). The reaction was stopped, and absorbance was read at 450 nm. The Blank value was subtracted from all values.
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



CLDN18.2: Anti-CLDN18.2 Binding Assay

