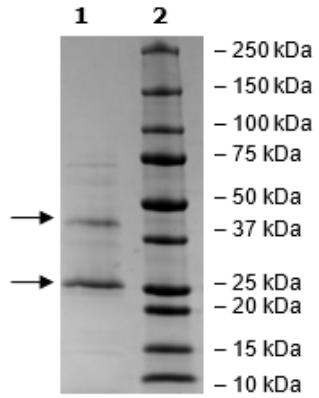


Product Information

Description:	Recombinant human Claudin-18, Isoform 2 (CLDN 18), full length, encompassing amino acids 1-261(end). This construct contains a C-terminal Flag-tag. The recombinant protein was affinity purified.
Species:	Human
Construct:	Claudin-18 Isoform 2 (1-261(end)-FLAG)
Concentration:	0.31 mg/ml
Expression System:	HEK293
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	50 mM HEPES, pH 8, 500 mM NaCl, 10% glycerol, 0.1% DDM, 0.01% CHS, and 100 µg/ml FLAG peptide
MW:	29 kDa
Genbank Accession:	NM_001002026
Stability:	At least 6 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:	The protein was validated by measuring its binding to Anti-Claudin-18 Isoform 2 IgG Antibody (BPS Bioscience #101564) in biochemical binding (ELISA) assay. The Claudin-18 Isoform 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-Claudin-18 Isoform 2 IgG 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 HRP Colorimetric Substrate (BPS Bioscience #79651). The reaction was stopped, and absorbance was read at 450 nm. The Blank value was subtracted from all values.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



CLDN 18.2: CLDN 18.2 Antibody Binding Assay

