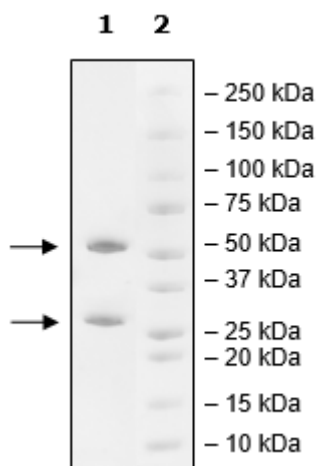


Product Information

Description:	This purified recombinant antibody is a human anti-CD38 (cluster of differentiation 38) antibody. This antibody recognizes human CD38.
Background:	CD38 (cyclic ADP-ribose hydrolase 1, ADPRC1) is a glycoprotein and ectoenzyme which plays an important role in regulating intracellular calcium. CD38 is a highly attractive target antigen for immunotherapy because it is highly expressed on multiple myeloma cells, and at relatively low levels on normal lymphoid and myeloid cells. Expression of CD38 has also been associated with HIV infection, leukemia, and type II diabetes mellitus. In 2015, the FDA approved Daratumumab (Darzalex), a breakthrough therapy drug targeting CD38, for the treatment of multiple myeloma.
Species:	Human
Clonality:	Monoclonal
Concentration:	1.37 mg/ml
Expression System:	HEK293
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	8 mM phosphate, 110 mM NaCl, 2.2 mM KCl, pH 7.4, and 20% glycerol
MW:	Heavy Chain: 49 kDa; Light Chain: 23 kDa
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:	The antibody was validated by measuring anti-CD38 binding to a CD38 antigen in an ELISA assay. The CD38 protein (BPS Bioscience #71277) was coated onto a 96-well plate overnight at 4°C (50 µl/well at a concentration of 4 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 non-biotinylated anti-CD38 antibody (BPS Bioscience #79120), 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 CD38 in ELISA and in cellular assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



CD38:Anti-CD38 Binding Assay

