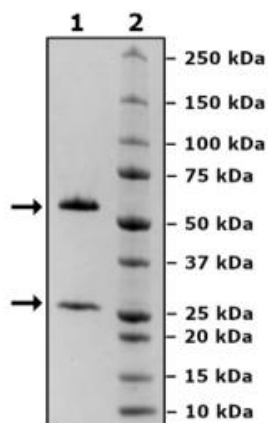


Product Information

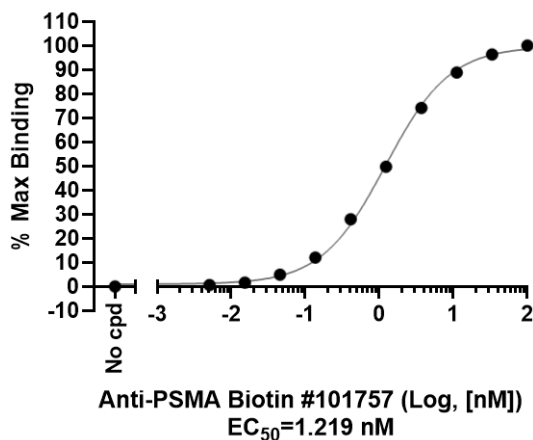
Description:	Recombinant human anti-PSMA antibody recognizing human PSMA protein. This anti-PSMA antibody is a purified recombinant antibody, which contains an Avi-Tag™ and is biotinylated at the C-terminus of the heavy chain unit.
Background:	PSMA (prostate-specific membrane antigen, also known as Folate hydrolase 1, FOLH1), is highly expressed in prostate cells and is used as a diagnostic and prognostic indicator of prostate cancer. The enzyme has folate hydrolase and peptidase activity. It plays a role in prostate cancer progression through the PI3K-Akt and MAPK-ERK1/2 pathways. PSMA/FOLH1 is used in the clinic in PET (positron emission tomography) imaging of prostate cancer whereas radioactive analogs are used for the treatment of prostate cancer. PSMA is also expressed in other tumor types and in a few normal tissues, such as kidney. PSMA/FOLH1 is a target of CAR-T cells and bi-specific antibodies are currently under development for cancer therapy.
Species:	Human
Isotype:	IgG1
Clonality:	Monoclonal
Concentration:	2.63 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: 52 kDa; Light Chain: 23 kDa + glycans
Glycosylation:	This antibody runs at a higher MW by SDS-PAGE due to glycosylation.
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be ≥90%.
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 antibody was validated by measuring Anti-PSMA binding to PSMA antigen in ELISA assay. The PSMA protein (BPS Bioscience #100463) 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 biotinylated anti-PSMA antibody (BPS Bioscience #101757), serially diluted in Blocking Buffer 2, was added for 30 minutes at room temperature. The plate was washed, incubated with Streptavidin-HRP, washed again, 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 PSMA in ELISA and in cellular assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



PSMA: anti-PSMA-Biotin Binding Assay



Biotin-Avidin Pulldown

