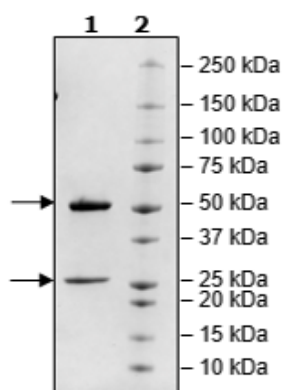


Product Information

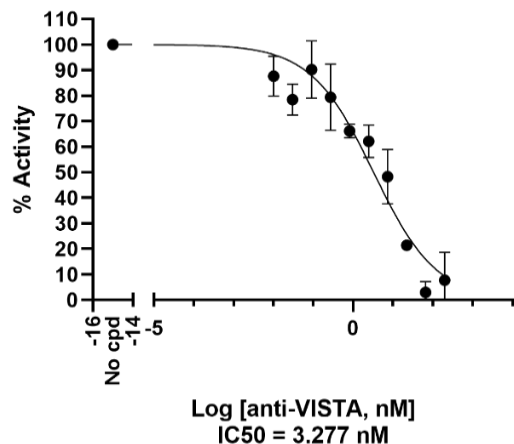
Description:	This human monoclonal antibody recognizes human V-domain immunoglobulin (Ig) suppressor of T cell activation (VISTA), and neutralizes its interaction with VSIG3 (also called IGSF11). VISTA is also known as B7-H5 or VSIR (GenBank Accession No. NM_022153). VSIG3 is a ligand involved in cell adhesion as a VISTA binding partner. This neutralizing antibody has been functionally tested using the "VSIG-3:VISTA [Biotinylated] Inhibitor Screening Assay Kit" (BPS Bioscience #79782).
Concentration:	2.15 mg/ml
Specificity:	The antibody recognizes a unique epitope on the surface of human VISTA protein. It has not been tested on other species.
Isotype	Human IgG1
Secondary detection:	Anti-human secondary
Formulated In:	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
Purification:	Protein A affinity chromatography
Format:	Aqueous buffer solution
Storage:	Store in aliquots at -80°C. Stable for at least 12 months from date of receipt. Avoid freeze/thaw cycles.
MW:	~150 kDa
Assay Conditions:	The functional activity of the antibody was measured using "VSIG-3:VISTA [Biotinylated] Inhibitor Screening Assay Kit" (BPS Bioscience #79782).
Applications:	This product is for research use only. It is not suitable for human, diagnostic or therapeutic use. The human monoclonal neutralizing IgG can be used for functional assays to block VISTA and VSIG3 binding.

Quality Control Data

4-20% SDS-Page Coomassie Staining



VSIG-3:VISTA-Biotin Activity



Anti-VISTA neutralizing antibody competes with and blocks the binding of VISTA and VSIG3. The IC₅₀ was calculated to be 3.28 nM using the “VSIG-3:VISTA [Biotinylated] Inhibitor Screening Assay Kit” (BPS Bioscience #79782). The assay was performed following the recommended kit protocol. Briefly, VSIG3 (BPS Bioscience #79491) was coated onto the surface of a 96-well microtiter plate overnight. The following day, VSIG3 was incubated with serial dilutions of anti-VISTA (0 to 200 nM, three-fold dilutions) and with 12 ng/μl (approximately 250 nM) Biotin-labeled VISTA (BPS Bioscience #71327). Next, the unbound proteins are washed off and Streptavidin-HRP (BPS Bioscience #79742) was added to the plate. Detection was accomplished by the addition of ELISA ECL Substrates A and B (BPS Bioscience #79670), prior to reading the signal using a luminescence plate reader.

Related Products

Products	Catalog #	Size
VSIG-3:VISTA [Biotinylated] Inhibitor Screening Assay Kit	79782	96 reactions
VSIG3, Avi-His-Tag	100349	100 μg
VSIG-3, Fc-fusion (IgG1), Avi-Tag	79491	100 μg
VSIG3, Fc-fusion (IgG1), Avi-Tag, Biotin-Labeled	100044	25 μg/50 μg
B7-H5, Avi-His-Tag	100358	100 μg
B7-H5 (VISTA), Fc fusion (Human) HiP™	71148	100 μg
B7-H5 Fc fusion, Avi-Tag, Biotin-labeled (Human)(VISTA-biotin)	71327	25 μg/50 μg