

B7-H5 Fc Fusion, Avi-Tag, Biotin-Labeled (Human) Recombinant

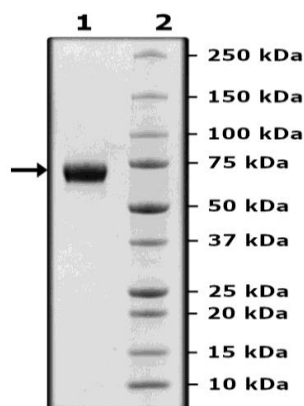
Catalog: 71327
Lot: 230301

Product Information

Description:	Recombinant human secreted B7-H5 Fc fusion protein also known as V-set immunoregulatory receptor (VSIR), encompassing amino acids 33-194. This construct contains a C-terminal Fc immunoglobulin domain followed by Avi-tag™ and biotin label. The recombinant protein was enzymatically biotin-labeled using Avi-tag™ and affinity purified.
Species	Human
Construct:	B7-H5 (33-194-Fc(IgG1)-Avi)-(Biotin)
Concentration:	1.55 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:	47 kDa + glycosylation
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_022153
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.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown

