

CD28, Fc Fusion, Avi-Tag, Biotin Labeled Recombinant

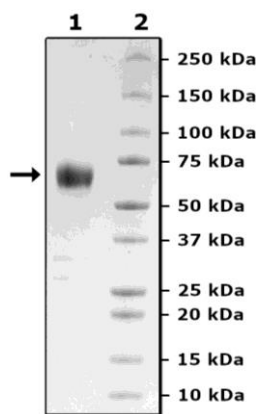
Catalog: 101580
Lot: 221026

Product Information

Construct:	CD28 (19-152-Fc(IgG1)-Avi)-(Biotin)
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
Concentration:	1.14 mg/ml
Species:	Human
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
Expression System:	HEK293
Format:	Aqueous buffer solution.
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession No.:	NM_006139
MW:	44 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	≥90%

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown

