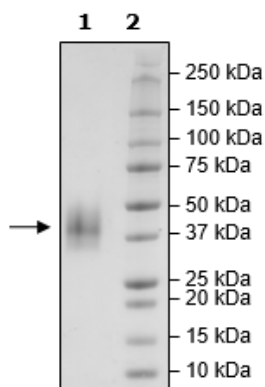


Product Information

Construct:	CD47 (19-139-His)
Concentration:	0.55 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:	NM_001777
MW:	15 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	≥90%
Assay Conditions:	Assay was performed according to CD47:SIRP-γ[Biotinylated] Inhibitor Screening Assay Kit (Human) (BPS Bioscience #72059) with SIRP-γ [Biotinylated] (BPS Bioscience #71171) titrated at various concentrations.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



CD47:SIRP- γ [Biotinylated] Activity

