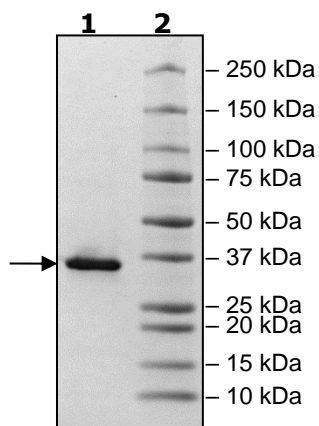


Product Information

Construct:	CD38 (44-308(end)-His) (dog)
Concentration:	0.39 mg/ml
Species:	Dog
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_001003143
MW:	31 kDa
Purity:	>90%
Assay Conditions:	Assay was done according to Dog CD38 Inhibitor Screening Assay Kit (Hydrolase Activity) (BPS Bioscience #78108) with CD38 titrated from 0.5 ng/μl (10 ng/rxn) to 0 ng/μl.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



CD38 (Dog) Hydrolase Activity

