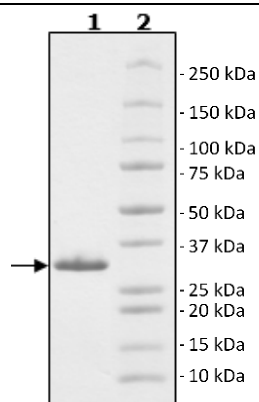


Product Information

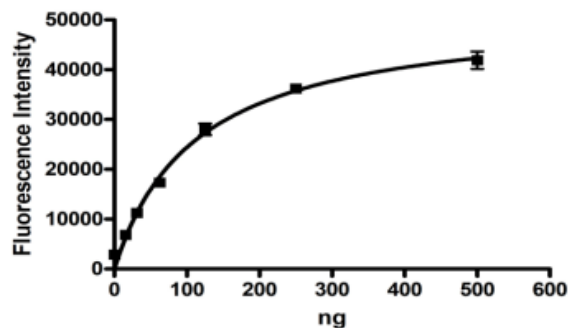
Construct:	CD38 (Pig) (54-286(end)-FLAG)
Concentration:	0.97 mg/ml
Species:	Pig
Formulated In:	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_001243883
MW:	28 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	≥90%
Aggregation:	<10%
Assay Conditions:	Assay was ran according to the CD38 (pig) Hydrolase Activity Assay Kit (BPS Bioscience #78178). Various amounts of enzyme incubated with etheno-NAD and incubated overnight. Fluorescence intensity measured at ex360 nm, em460 nm.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Pig CD38 Hydrolase Activity



Gel Filtration

