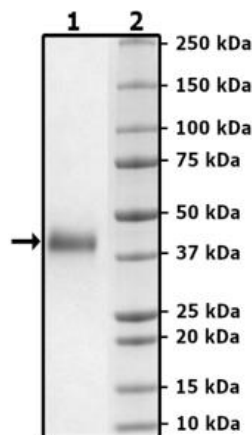


Product Information

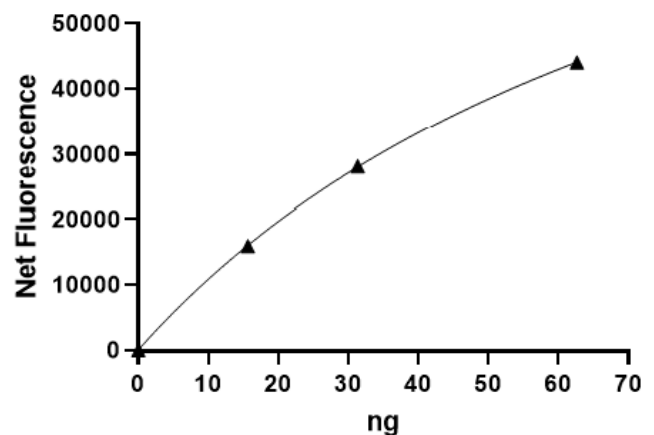
Description:	Recombinant human CD38 (CD38 molecule), encompassing amino acids 43-300(end). This construct contains a C-terminal His-tag (6xHis). This recombinant protein was affinity purified. HiP™ indicates a high purity protein (≥90% pure) and less than 10% aggregation as measured by gel filtration.
Species	Human
Construct:	CD38 (43-300(end)-His)
Concentration:	4.36 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:	31 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation
Aggregation:	<10%
Genbank Accession:	NM_001775
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.
Assay Conditions:	Assay was done according to CD38 Inhibitor Screening Assay Kit (Cyclase Activity) (BPS Bioscience #71275) with CD38 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



CD38 (Cyclase) Activity



Gel Filtration Curve

