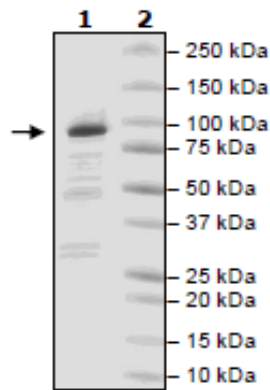


Product Information

Construct:	A20 (FLAG-2-790(end)-His)
Concentration:	0.09 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 3mM DTT and 20% glycerol.
Expression System:	Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_006290
MW:	91 kDa
Purity:	79%
Specific Activity:	0.77 pmol/min/μg
Assay Conditions:	Reaction was performed in 50 mM Tris pH 7.4, 1 mM DTT, 0.5 mM EDTA, 500 nM Ub-AMC, and various amounts of A20 in a volume of 50 μl. Reaction was incubated at 37°C for 30 min and fluorescent signal was measured at excitation of 360 nm, and emission at 460 nm.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

