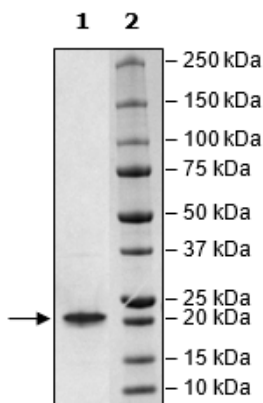


Product Information

Construct:	UEV1A (FLAG-2-170(end))
Concentration:	0.20 mg/ml
Species:	Human
Formulated In:	45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 10% glycerol, 3mM DTT, 100 µg/ml FLAG peptide.
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_021988
MW:	20 kDa
Purity:	≥90%
Assay Conditions:	Reaction was performed according to the UBCH13 TR-FRET Assay Kit (BPS Bioscience #79741) with various amounts of UEV1A.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



UEV1A Activity

