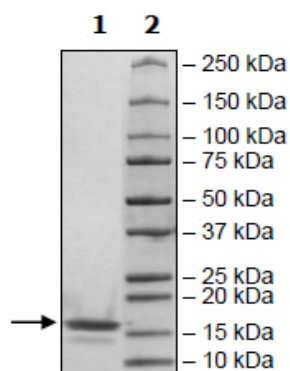


Product Information

Construct:	Ubch5a (UBE2D1) (His-2-148(end))
Concentration:	1.68 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, 20% Glycerol and 250 mM Imidazole.
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_003338
MW:	17 kDa
Purity:	90%
Assay Conditions:	Reaction was performed according to the UBCH5a TR-FRET Assay Kit (BPS Bioscience #79900) with UBCH5a titrated from 80 ng/rxn – 1.25 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



UBCH5a Activity

