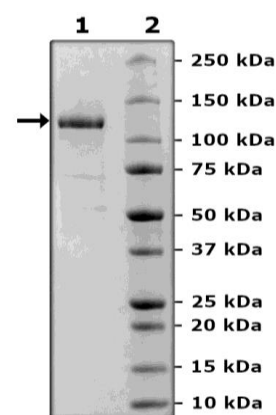


Product Information

Description:	Human USP20, also known as ubiquitin specific peptidase 20 and Deubiquitinating Enzyme 20, GenBank Accession No. NM_006676, a.a. 2–915 (end) with N-terminal His-Tag and C-terminal FLAG-Tag, expressed in a HEK293 expression system. MW = 104 kDa.
Species	Human
Construct:	USP20 (His-2-915(end)-FLAG)
Concentration:	0.31 mg/ml
Expression System:	HEK293
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 100 µg/ml FLAG peptide
MW:	104 kDa
Genbank Accession:	NM_006676
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.
Specific Activity:	10 pmol/min/µg
Assay Conditions:	50 mM Tris-HCl, 0.5 mM EDTA, 0.05% Tween 20, 1 mM DTT, 500 nM Ubiquitin-AMC. Incubate at room temperature for 30 minutes. Fluorescence intensity is measured at exc/em of 360nm/460nm using Tecan M1000 plate reader.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



USP20 Activity

