

Product Information

Description:	Recombinant human USP11 (ubiquitin specific peptidase 11), encompassing amino acids 2-920. This construct contains a N-terminal His-tag (6xHis) and a C-terminal FLAG-tag. This protein was affinity purified.
Background:	USP1 (ubiquitin specific protease 11), or ubiquitin carboxyl-terminal hydrolase 11, belongs to the USP family of proteins. USP11 is a DUB (deubiquitinating enzymes) and is involved in removing K-48 and K-63 ubiquitin chains from several substrates. It participates in DNA repair and inflammation, can regulate immune cell differentiation via the NF- κ B and TNF (tumor necrosis factor) pathways. It has been linked to the development and progression of several cancer types, such as breast, hepatocellular and ovarian cancer. The development of USP11 specific inhibitors will bring therapeutic benefits in cancer therapy.
Species:	Human
Construct:	USP11 (His-2-920(end)-FLAG)
Concentration:	0.11 mg/ml
Expression System:	HEK293
Purity:	80%
Format:	Aqueous buffer solution.
Formulated In:	45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 0.045% Tween-20, and 10% glycerol
MW:	107 kDa
Genbank Accession:	NM_001371072.1
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.
Applications:	Useful for SDS-PAGE.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

