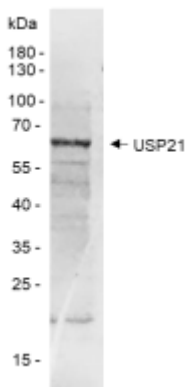


Product Information

| | |
|------------------------------|--|
| Description: | Recombinant human full length USP21 (Ubiquitin Specific Protease 21). This construct contains an N-terminal His-tag followed by a SUMO tag. This protein was affinity purified. |
| Background: | USP21 (Ubiquitin Specific Protease 21) regulates transcription initiation via deubiquitination of histone H2A and exhibits both deubiquitinase and deNEDDylase activities. |
| Species: | Human |
| Construct: | USP21 (His-SUMO-1-565(end)) |
| Concentration: | 1.25 mg/ml |
| Expression System: | <i>E. coli</i> |
| Purity: | ≥85% by SDS-PAGE |
| Format: | Aqueous buffer solution. |
| Formulated In: | 25 mM Tris-HCl, pH 7.4, 150 mM NaCl, 5 mM DTT, 10% glycerol |
| MW: | 74.7 kDa |
| 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. |
| Assay Conditions: | 100 nM USP21 was tested in the Ub-Rh110 assay at several time points. |
| Applications: | Suitable for deubiquitinase and deNEDDylase studies. |

Quality Control Data

SDS-PAGE Coomassie Staining



Activity

