



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

USP33 Core Domain

Human, recombinant, His₆-SUMO tag

Catalog #: 81102

Lot #: 141217 **Conc.:** 6.5 µM

Formulated in: 25 mM Tris-HCl, pH 7.5, 150 mM NaCl, 5 mM DTT, 10% glycerol.

Stability: At least 12 months at -80°C. *Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

References:

1. De Pitta, C., *et al.*, *Haematologica*. 2005; **90(7)**: 890-898.
2. Li, Z., *et al.*, *J. Biol. Chem.* 2002; **277(7)**: 4656-4662.

Description: Human Ubiquitin-specific protease 33 (USP33) catalytic domain, also known as deubiquitinating enzyme 33, KIA1097, ubiquitin carboxyl-terminal hydrolase 33, VDU1, and VHL-interacting deubiquitinating enzyme 1 with a His6-SUMO tag, GenBank Accession No. NM_015017.4, a.a. 181-712, MW = 61 kDa.

Activity: Activity of this enzyme was confirmed in the Ub-CHOP assay.

Application: This enzyme is useful for the study of enzyme kinetics and in *in vitro* applications of deubiquitination studies.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com