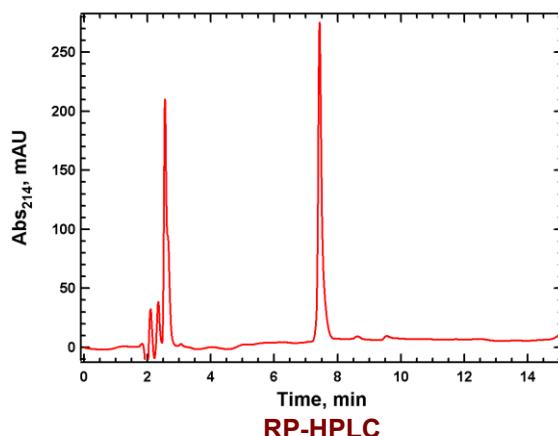


Background	UBE2V2 (also known as MMS2 or UEV-2) is a ubiquitin E2 variant (UEV) protein that forms a heterodimeric complex with the E2 conjugating enzyme UBE2N. This complex functions in the ubiquitination of target substrates alongside E1 and E3 enzymes. The UBE2N/UBE2V2 complex specifically assembles K63-linked polyubiquitin chains, which are utilized in non-degradative ubiquitin signaling pathways, such as DNA repair and signal transduction. UEV proteins are structurally similar to E2 enzymes, displaying the same fold but lacking the catalytic cysteine residue. Notably, the UBE2N/UBE2V2 complex is unique in its ability to form free K63-linked polyubiquitin chains in solution, even in the absence of an E3 ligase.
Alternate Names	Ubiquitin Conjugating Enzyme E2 V2, Enterocyte Differentiation-Associated Factor 1 (EDAF-1), MMS2 Homolog, Enterocyte Differentiation-Promoting Factor 1 (EDPF-1), 1 Alpha,25-Dihydroxyvitamin D3-Inducible DDVit-1
Application	Ubiquitin ligation reactions

Product Specifications

Tag	None
Purity	≥ 95% by RP-HPLC
Molecular Weight	16,363.1 Da by MS (calculated 16,362.8)
Quantity	20 µl or 75 µl of a 40 µM solution (0.8 and 3 nmoles, respectively)
Species	Human, recombinant; Accession No. Q15819
Expression System	<i>E. Coli</i>
Physical State	Liquid
Buffer	25 mM Tris, pH 7.4; 150 mM NaCl; 10 mM DTT; 10% glycerol;
Solubility	> 3 mg/ml
Stability & Storage	1 year at -80° C. Avoid repeated freeze/thaw cycles

Product QC



References

1. Hofmann, R.M. and C.M. Pickart. J Biol Chem, 2001; **276**, 27936-43.
2. VanDemark, A.P., et al., Cell, 2001; **105**, 711-20 (2001).
3. Eddins, M.J., et al., Nat Struct Mol Biol, 2006 **13**, 915-20 (2006).
4. Fletcher AJ., et al., EMBO J. 2015;34(15):2078-95.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2025 LifeSensors, Inc. All Rights Reserved