

MMS2 (yeast)

Cat. # UB206

Background

MMS2 is a ubiquitin E2 variant (UEV) protein that acts in concert with the E2 enzyme Ubc13. This heterodimeric complex assembles K63-linked polyubiquitin chains for non-degradative ubiquitin signaling pathways (DNA repair, signal transduction). The UEV proteins are structurally similar to E2 enzymes; they display the same fold, yet they lack the active site cysteine indicative of E2 enzymes. The Ubc13/MMS2 complex is unique in that it has the ability to form free K63-linked polyubiquitin chains in solution in the absence of an E3.

Alternate Names

Methyl Methanesulfonate Sensitive 2, Ubiquitin-Conjugating Enzyme E2 Variant 2 (UBE2V2), UEV protein, and Mms2p

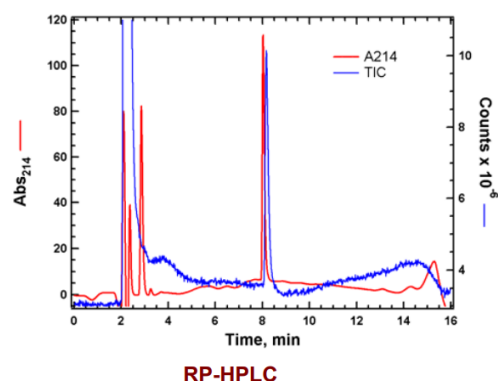
Application(s)

Ubiquitin ligation reactions

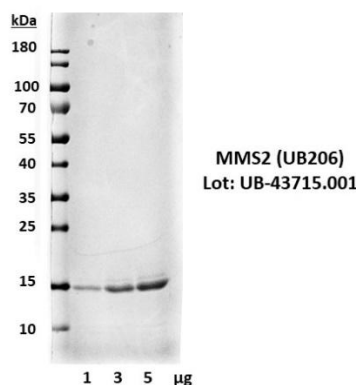
Product Specifications

Affinity tag	None
Purity	> 95% by RP-HPLC
Molecular Weight	15,544.9 Da by MS (calculated 15,544.7)
Quantity	20 μ l or 75 μ l of a 40 μ M solution (0.8 and 3 nmoles, respectively)
Species	<i>S. cerevisiae</i>
Expression System	<i>E. coli</i>
Physical State	Liquid
Buffer	50 mM Tris, pH 7.5; 150 mM NaCl; 10 mM DTT; 10% glycerol
Solubility	> 3 mg/ml
Stability & Storage	Over 1 year at -80°C. Avoid repeated freeze/thaw cycles

Product QC



MMS2 purity gel



References

1. Andersen, P. L., et al., The Journal of cell biology, 2005; 170(5), 745–755.
2. VanDemark, A.P., et al., Cell, 2001; 105:711-20.
3. Eddins, M.J., et al., Nat Struct Mol Biol, 2006; 13:915-20.

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