

E6-AP (Human papillomavirus E6-associated protein)

Cat. # UB304

Background: E6-AP (E6-associated protein) is an E3 (ubiquitin protein ligase) that is involved in the conjugation of ubiquitin to target substrates along with E1 and E2 enzymes. It is a HECT domain E3 containing an active site cysteine which accepts a charged ubiquitin from the ubiquitin-E2 thiolester complex. E6-AP has been linked to a neurodevelopmental disorder, Angelman syndrome. In addition the oncoprotein E6 of human papillomaviruses (HPV) utilizes E6-AP to target numerous cellular proteins for degradation including the tumor suppressor protein p53.

Alternate names: Ubiquitin-protein ligase E3A

Product Information

Purity:	≥ 85% by SDS-PAGE
Molecular Weight:	101kDa
Quantity:	25 µg
Physical State:	Liquid
Source:	Human Recombinant
Tag:	None
Activity:	Typical enzyme concentration of 100 nM - 5 µM is used for in vitro conjugation, depending on conditions.
Storage:	-80° C. Avoid repeated freeze/thaw cycles

References

- 1) Huibregtse, J.M., M. Scheffner, and P.M. Howley, Cloning and expression of the cDNA for E6-AP, a protein that mediates the interaction of the human papillomavirus E6 oncoprotein with p53. *Mol Cell Biol*, 1993. **13**:775-84.
- 2) Huibregtse, J.M., M. Scheffner, and P.M. Howley, Localization of the E6-AP regions that direct human papillomavirus E6 binding, association with p53, and ubiquitination of associated proteins. *Mol Cell Biol*, 1993. **13**:4918-27.
- 3) Scheffner, M., et al., The HPV-16 E6 and E6-AP complex functions as a ubiquitin-protein ligase in the ubiquitination of p53. *Cell*, 1993. **75**:495-505.
- 4) Huibregtse, J.M., et al., A family of proteins structurally and functionally related to the E6-AP ubiquitin-protein ligase. *Proc Natl Acad Sci USA*, 1995. **92**:5249.
- 5) Huang, L., et al., Structure of an E6AP-UbcH7 complex: insights into ubiquitination by the E2-E3 enzyme cascade. *Science*, 1999. **286**:1321-1326.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2009 LifeSensors, Inc. All Rights Reserved