

E6-AP (Human papillomavirus E6-associated protein)

Cat. # UB304

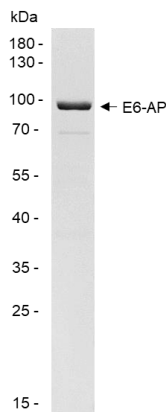
Background

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.

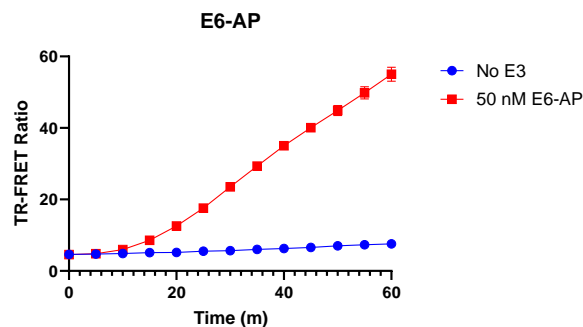
Product Information

Purity	≥ 85% by SDS-PAGE
Molecular Weight	101 kDa
Quantity	25 µg
Physical State	Liquid, 20 mM Hepes pH 7.5, 150 mM NaCl, 10% glycerol
Species	Human
Source	<i>E. coli</i>
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

Product QC



SDS-Page Analysis of purified E6-AP. Two µg of the protein was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of E6-AP. 50 nM E6-AP was tested in a TR-FRET assay showing robust E3 ligase activity.

References

- Owais, A., et al., Cancers (Basel)., 2020. 12(8):2108.
- Sailer, C., et al., Nat Commun., 2018. 9(1):4441.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic use.

Copyright © 2009 LifeSensors, Inc. All Rights Reserved