

His₆-TUBE1

Cat. # UM201

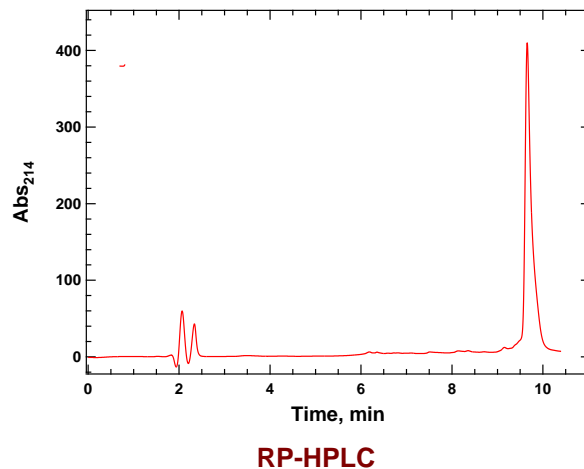
Background: Based on protein domains known to possess an affinity for ubiquitin, Tandem Ubiquitin Binding Entities (TUBEs) have been developed for the isolation and identification of ubiquitylated proteins. TUBEs display up to a 1000-fold increase in affinity for poly-ubiquitin moieties over the single ubiquitin binding associated domain (UBA). In addition, TUBEs display a protective effect on polyubiquitinated proteins, allowing for detection at relatively low abundance. These properties effectively "capture" proteins in their polyubiquitylated state.

The His₆-TUBE can be used in conjunction with Ni-IMAC supports to isolate or pull-down polyubiquitylated proteins.

- Application:**
- Protection of polyubiquitylated proteins from deubiquitylation and degradation by the proteasome
 - Pull down of polyubiquitylated proteins from cell lines, tissues and organs

Product Information

Purity:	≥ 90% by RP-HPLC
Molecular Weight:	37,696.9 Da
Tag:	His ₆ -SUMO
Physical State:	liquid
Quantity:	200 µg
Storage:	-80° C. Avoid repeated freeze/thaw cycles



References

Hjerpe, R, Aillet, F, Lopitz-Otsoa, F, Lang, V, England, P, and Rodriguez, MS., [Efficient protection and isolation of ubiquitylated proteins using tandem ubiquitin-binding entities.](#) *EMBO Rep.* **10**,1250-1258 (2009).

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