

Tandem Ubiquitin Binding Entity

ISOLATION | DETECTION | PROTECTION

IP Ubiquitylated protein
without overexpression
or Proteasome Inhibitor

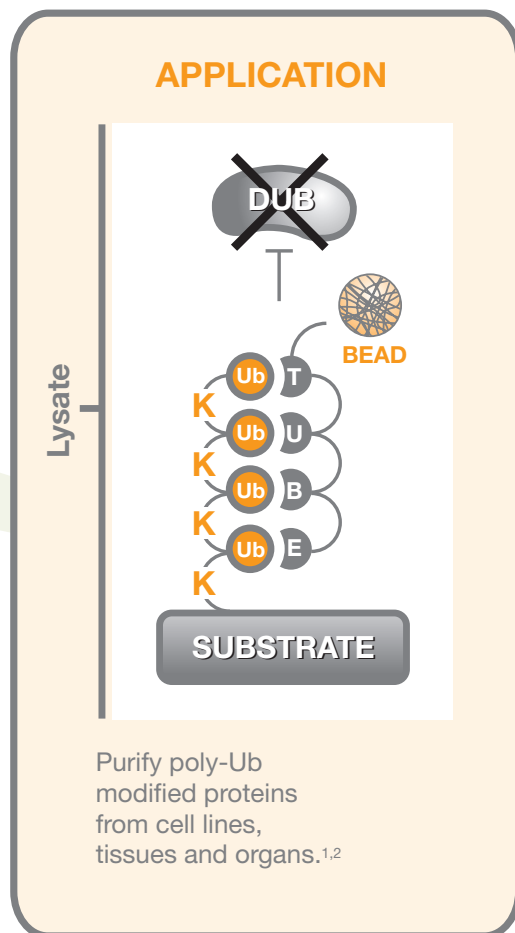
Up to 1000X higher affinity than
a single UBA domain

Protect ubiquitylated proteins
from the degradation

Tandem Ubiquitin Binding Entities (TUBEs)
are high affinity ubiquitin traps that bind to
poly-ubiquitin chains.

TUBEs bind ubiquitin chains with low
nanomolar affinity and can be used for
inhibition of deubiquitylation and protein
degradation; capture of total ubiquitylated
proteins from cell cultures, tissues and
organs; and many other uses.

TUBEs Mechanism of Action



Ref: Hjerpe et al. *EMBO Reports*, Oct 2009



**Contact Us
To Learn More:
info@lifesensors.com**

To order, please call
610-644-8845 or visit
www.lifesensors.com

Proven Results
Improved Versatility
Which TUBE will work for you?

Tagged Tubes

GST-TUBE1
GST-TUBE2

His₆-TUBE1
His₆-TUBE2
His₆-TUBE3

Biotin-TUBE1
Biotin-TUBE2

Immobilized TUBEs

Agarose-TUBE1
Agarose-TUBE2

Poly-Ubiquitylated Proteins: Improve Your Recovery and Detection with Our Expanded TUBEs

When it comes to poly-ubiquitylated protein recovery, TUBEs have rapidly eclipsed conventional technologies and are a great alternative for ubiquitin antibodies. Now, with 10 variants currently available, TUBEs truly let you do more for less.

- Higher Binding Capacity for Pull-Downs
- Higher Specificity than Antibodies/UBAs
- Detect Poly-Ubiquitylation by Western Without Boiling
- Lower Cost per Experiment

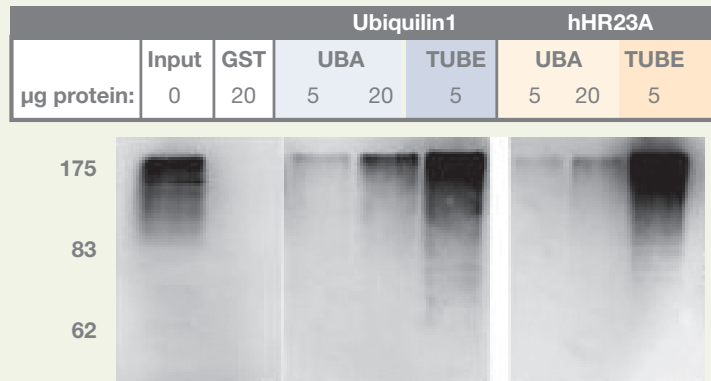


Tandem Ubiquitin Binding Entities (TUBEs) are based on known UBAs, yet display up to 1000-fold higher affinity than the monomer form. An array of TUBE technologies now provides added flexibility for pull-downs and recovery of your poly-ubiquitylated proteins. Solution phase TUBEs have been expanded to offer your choice of affinity tag, including biotinylated TUBEs for detection of poly-ubiquitin by western blotting. And for the added ease and convenience of one-step pull-downs, each TUBE is now available in immobilized form. No matter what your experimental needs, TUBEs afford high affinity binding and protection from degradation for your poly-ubiquitylated protein.

Suggested Applications Include:

- Pulldown and analysis of your poly-ubiquitylated protein
- Detection of poly-ubiquitylated proteins by “far western”
- In situ detection of poly-ubiquitylated by “far western”
- Protection of poly-ubiquitylated proteins from UPP degradation

Purification of Poly-Ubiquitylated proteins using TUBE Technology



α – polyubiquitin

Ref: Hjerpe et al. *EMBO Reports*, Oct 2009