

Tandem Ubiquitin Binding Entity

ISOLATION | DETECTION | PROTECTION

Up to 1000X higher affinity than a single UBA

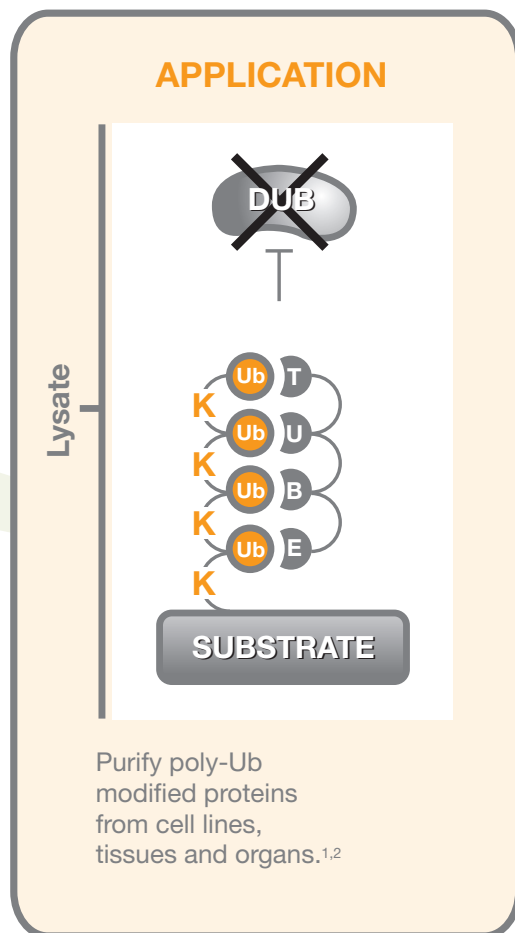
Pull-down Ubiquitylated proteins without overexpression or addition of DUB or Proteasome Inhibitors

Protect ubiquitylated proteins from DUBs

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; capture total ubiquitylated proteins from cell cultures, tissues extracts and for other applications.

TUBEs Mechanism of Action



Ref: Hjerpe et al. EMBO Reports. 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

* TUBE1 is derived from ubiquitin UBA
TUBE2 is derived from human Rad23A
(UBA domain 1)

Immobilized TUBEs**

Agarose-TUBE1
Agarose-TUBE2

** For additional protection from
DUBs, use PR619 (Cat.no. SI619),
a reversible pan DUB inhibitor, during
purification with Agarose-TUBEs

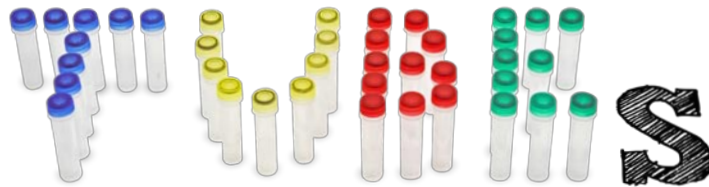


**Polyubiquitylated Proteins:
Improved Recovery and Detection
with Our Versatile TUBE Collection**

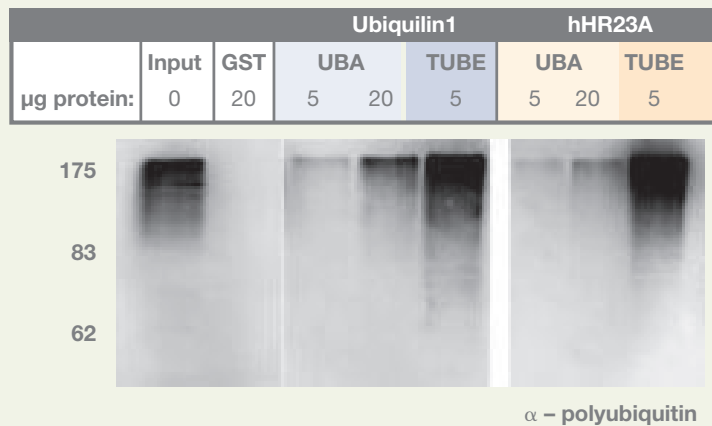
When it comes to polyubiquitylated proteins, TUBEs rapidly eclipsed conventional technologies and are a great alternative to ubiquitin antibodies. Now with new variants currently available, TUBEs truly let you do for more for less.

- High Binding Capacity for Pull-Downs
- Higher Specificity than Antibodies and UBAs
- Lower Cost per Experiment

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 additional flexibility for the pull-down and recovery of polyubiquitylated proteins. For one step pull-down, each TUBE is available in immobilized form (Agarose-TUBEs). Select from several affinity-tagged TUBEs (His₆- or GST) for IMAC purification or use Biotin-TUBEs for the Western blot analysis. No matter what your experimental needs, TUBEs afford high affinity binding and protection from deubiquitylation for your poly-ubiquitylated protein.



**Purification of Poly-Ubiquitylated proteins
using TUBE Technology**



Hjerpe et al. EMBO Reports, 2009