

Detection Reagent 1 for UC101

Cat. # DR101

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.

- Application:**
- Detection of polyubiquitylated proteins by ligand blotting
 - Pull down of polyubiquitylated proteins from cell lines, tissues and organs using a variety of readily available avidin supports
 - *In situ* labeling for detection of polyubiquitin by histochemistry
-

Product Information

Purity:	>95% by SDS-PAGE
Physical State:	Liquid
Quantity:	35 µL
Storage:	-80° C. Avoid repeated freeze/thaw cycles

References

- Donaghy, Ryan, et al. “The BRISC deubiquitinating enzyme complex limits hematopoietic stem cell expansion by regulating JAK2 K63-ubiquitination.” *Blood*, vol. 133, no. 14, 2019, pp. 1560-1571.
- Lv, Kaosheng, et al. “CBL family E3 ubiquitin ligases control JAK2 ubiquitination and stability in hematopoietic stem cells and myeloid malignancies.” *Genes & Development*, vol. 31, 2017, pp. 1007-1031.

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