

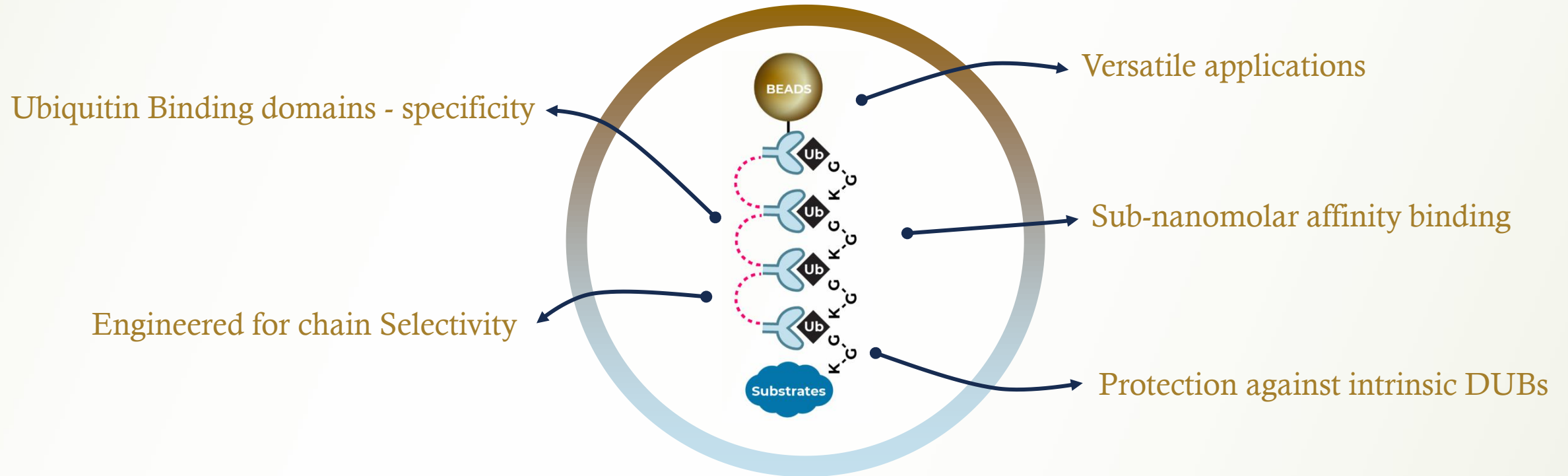
# LifeSensors TUBEs: TANDEM UBIQUITIN BINDING ENTITIES

POLYUBIQUITIN BINDING DOMAINS

# LifeSensors Inc. Mission

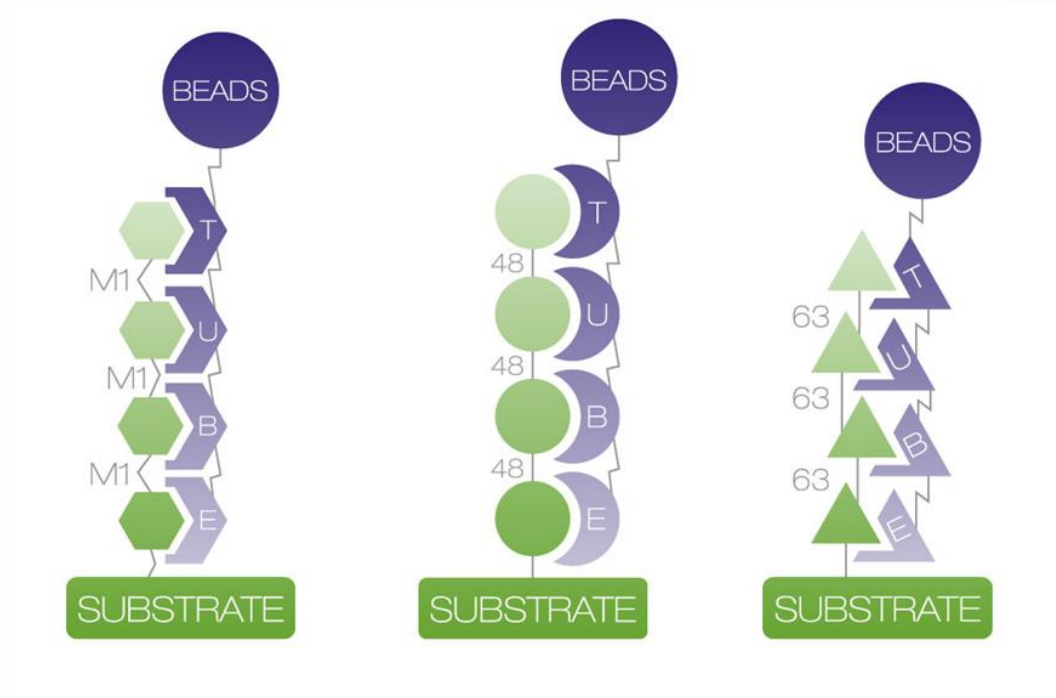
- Leadership in UPS, [PROTAC](#), [DUBTAC](#), and [Molecular Glues](#)
- Drug Discovery, [UPS Enzymes](#), [DUBs](#), [PROTAC Screening Services](#)
- Biomarker Development and Collaborative Research
- ~500 Products, [DUBs](#), [E3 ligases](#), [Ubiquitin Affinity Matrices \(TUBEs\)](#), [Assay Kits](#) and Proprietary [Protein Expression Systems \(SUMO\)](#)
- Profiling Compounds Against [Ubiquitin Ligases](#) and [DUBs](#)

# TUBEs: TANDEM UBIQUITIN BINDING ENTITIES



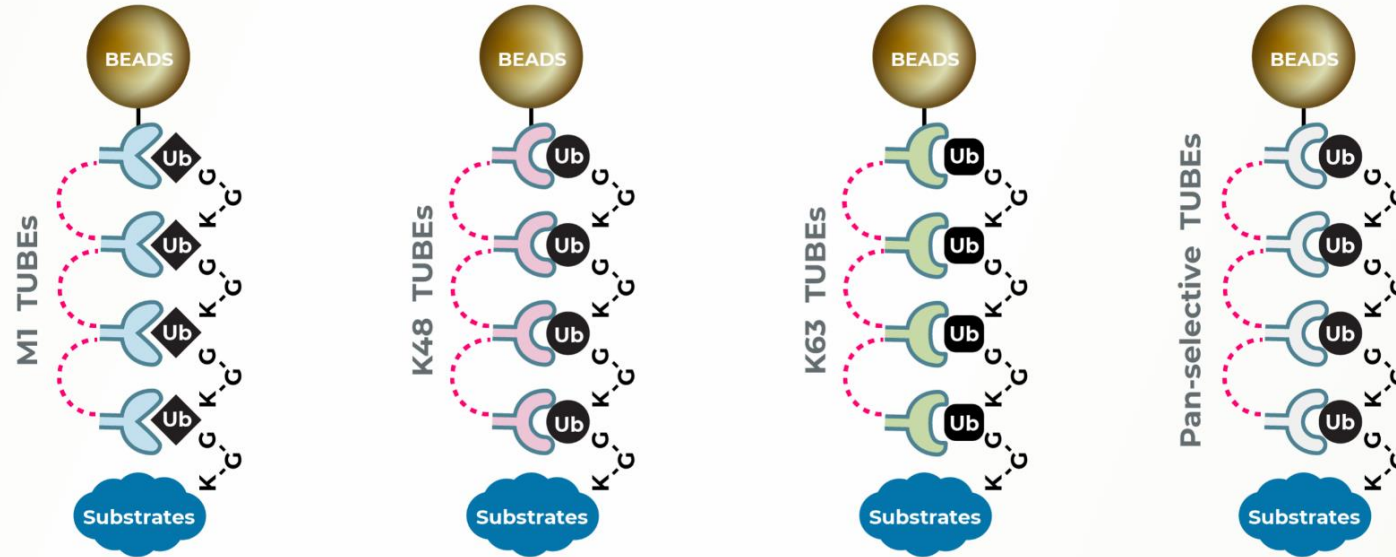
# What are TUBEs?

TUBEs are high affinity 'ubiquitin traps' that capture poly-ubiquitinated proteins with low nanomolar affinity.

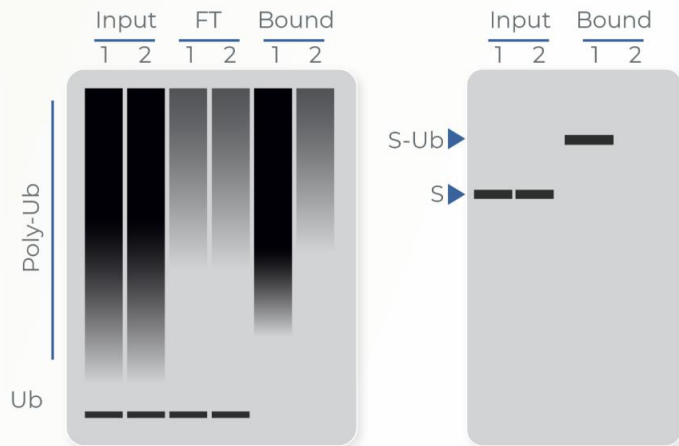


TUBEs can be used for capture or detection of total polyubiquitin or enrichment of linkage specific proteins K63-linked polyubiquitin K48-linked polyubiquitin, and M1-linked polyubiquitin.

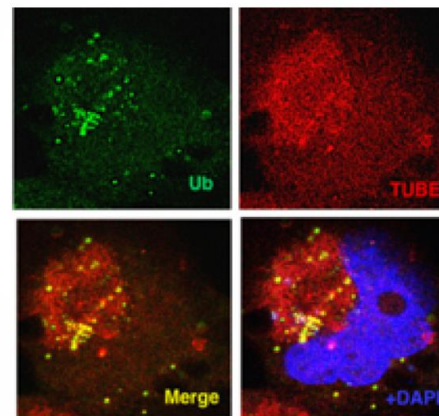
# TANDEM UBIQUITIN BINDING ENTITIES (TUBE<sub>s</sub>)



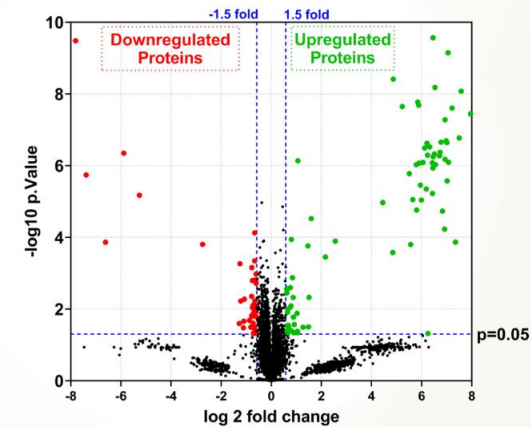
Poly-Ubiquitin Substrate Enrichment & Western Blotting



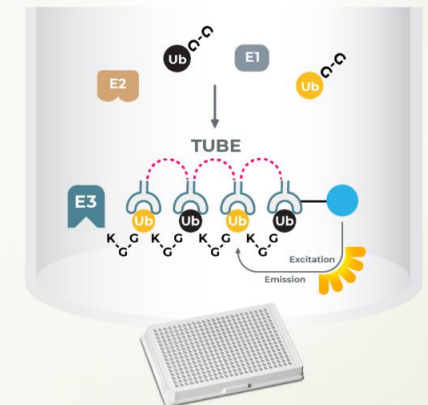
Fluorescent Microscopy



Mass Spectrometry Ubiquitomics



High-Throughput Screening





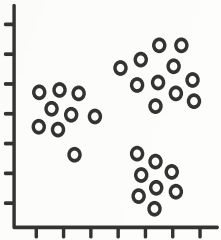
# TUBEs address critical TPD challenges



**Strong binders / better  $K_D$  does not guarantee better degradation**



**Ternary complex rigidity**



**Ligand dependent ubiquitination, true functional HTS**

# The Power of **TUBE** Applications

**Magnetic  
TUBEs**

**Pull-down of  
polyubiquitinated  
proteins**

**Biotin  
TUBEs**

**Histochemistry/  
Cytochemistry**

**Agarose  
TUBEs**

**GST  
TUBEs**

**Biomarkers**

**His6  
TUBEs**

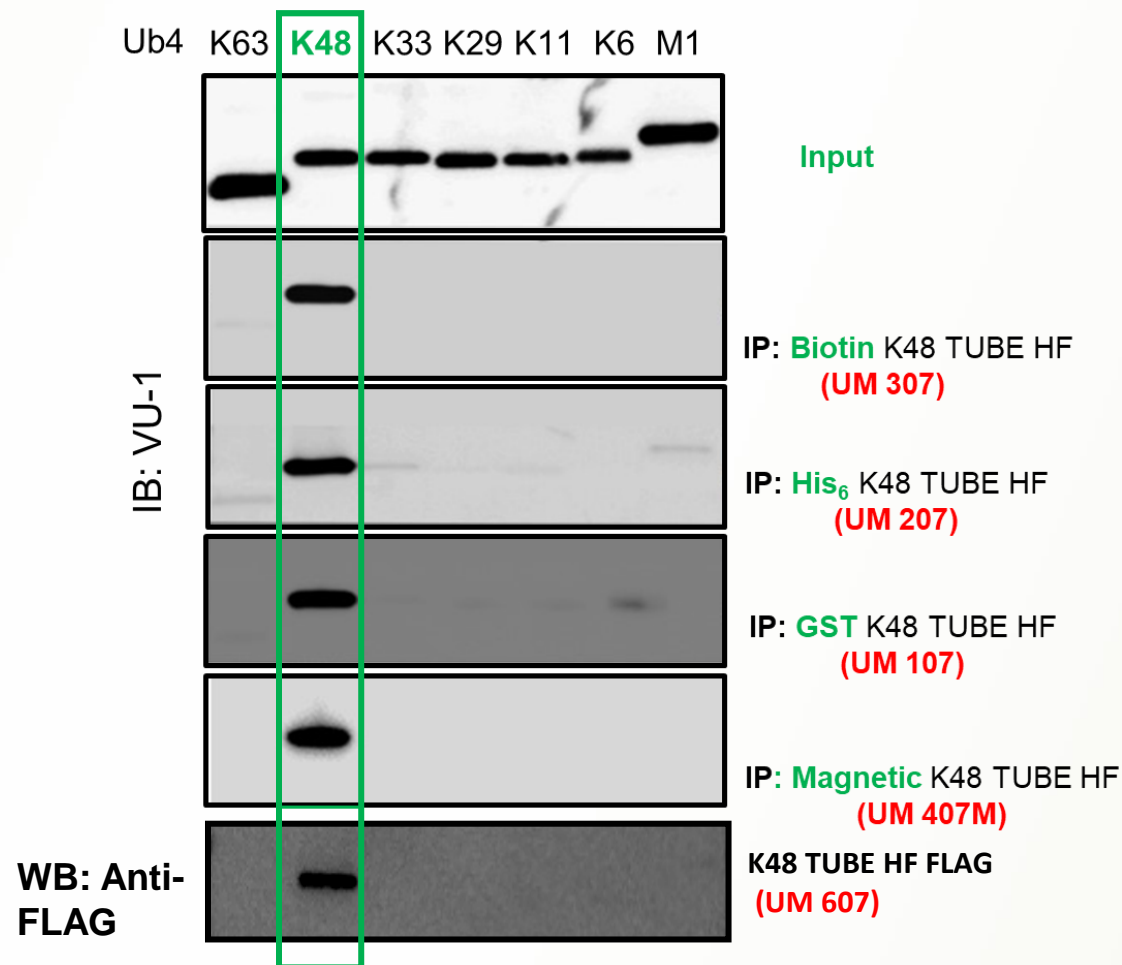
**‘Far Western’  
blot**

**TAMRA-  
TUBEs**

**Mass Spec  
&  
Proteomics**

**Fluorescein  
TUBEs**

# Lysine 48 Poly-ubiquitin TUBES are Highly Selective

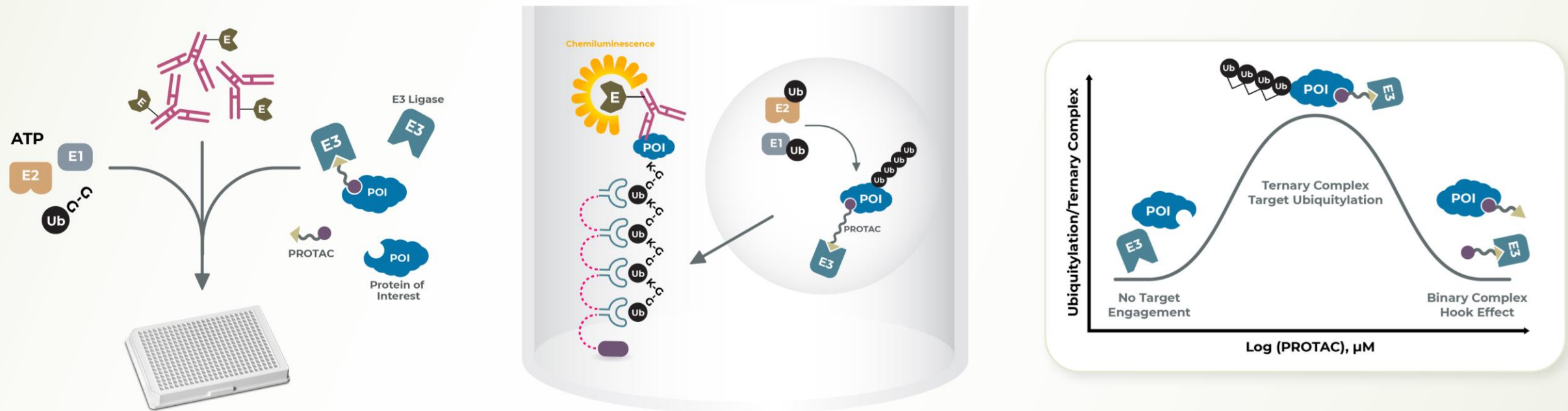


K48 TUBE HF FLAG detects only  
K48-linked ubiquitin chains in  
Western Blots



# TUBE Application for HTS- In Vitro Biochemical Assay

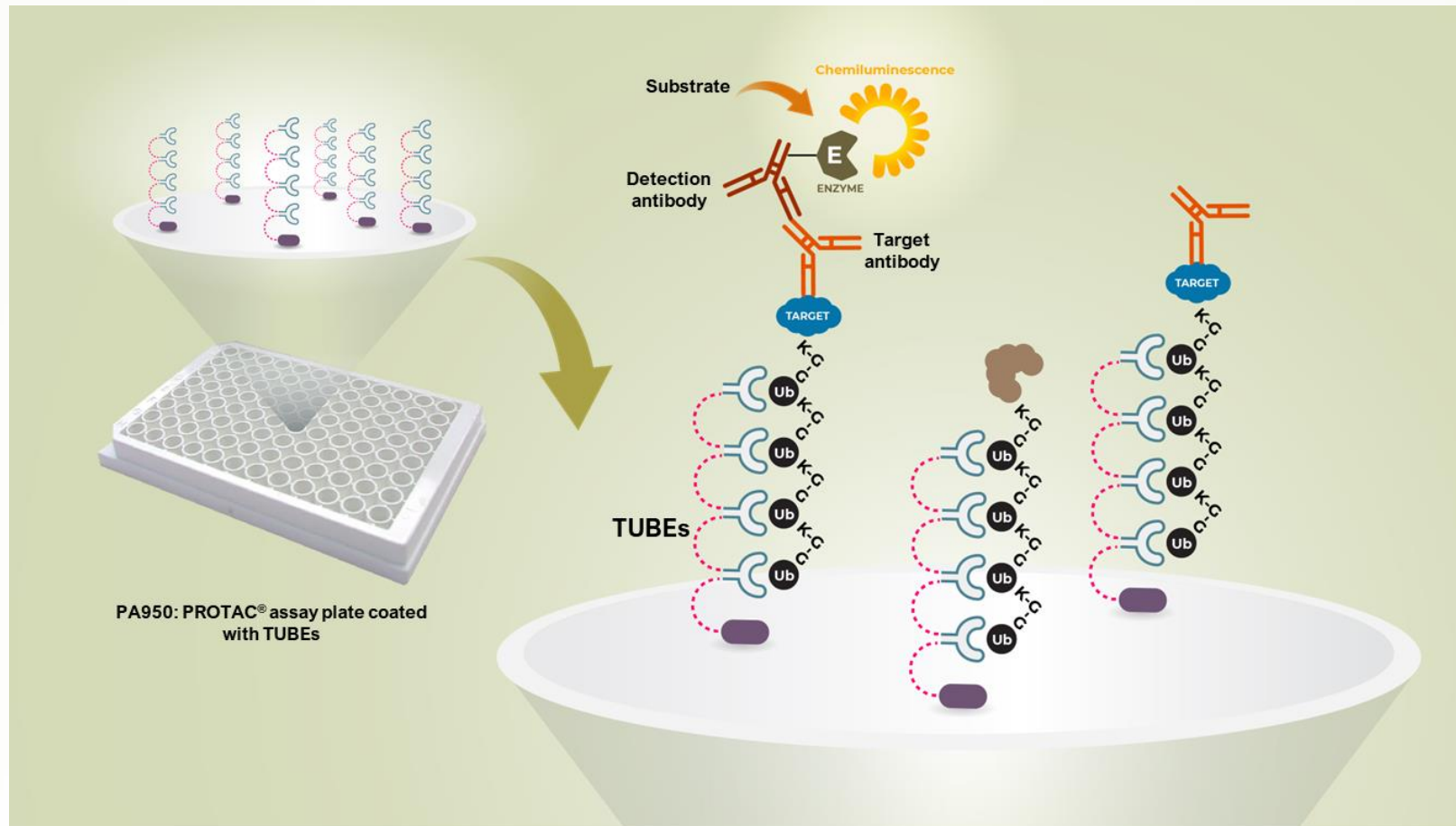
To study functional ternary complex and 'PROTACability'



TUBE Capture & PROTAC Mediated Ubiquitination of POI Detection

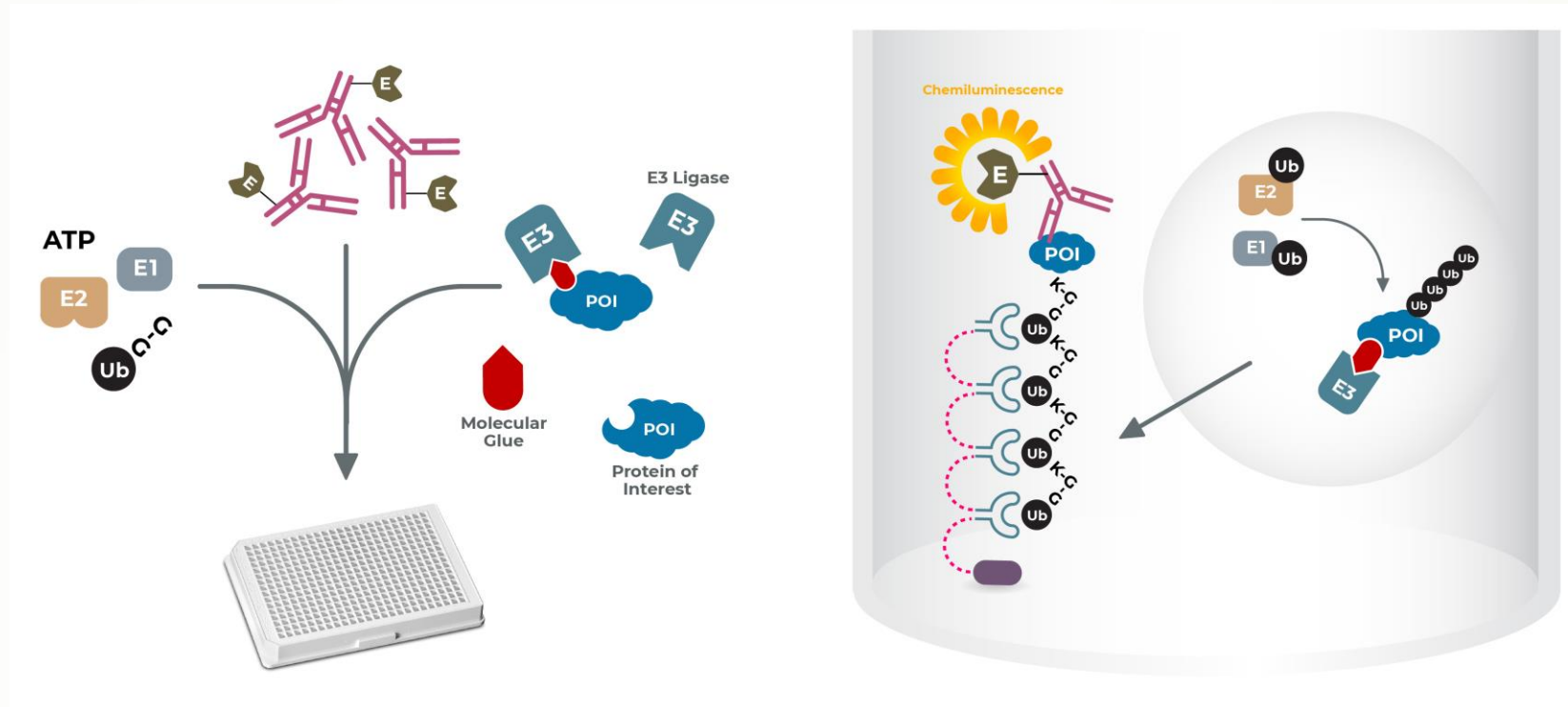
# Cell-Based PROTAC Assay using TUBE-coated plate

Directly monitor PROTAC-mediated ubiquitination and degradation of target protein

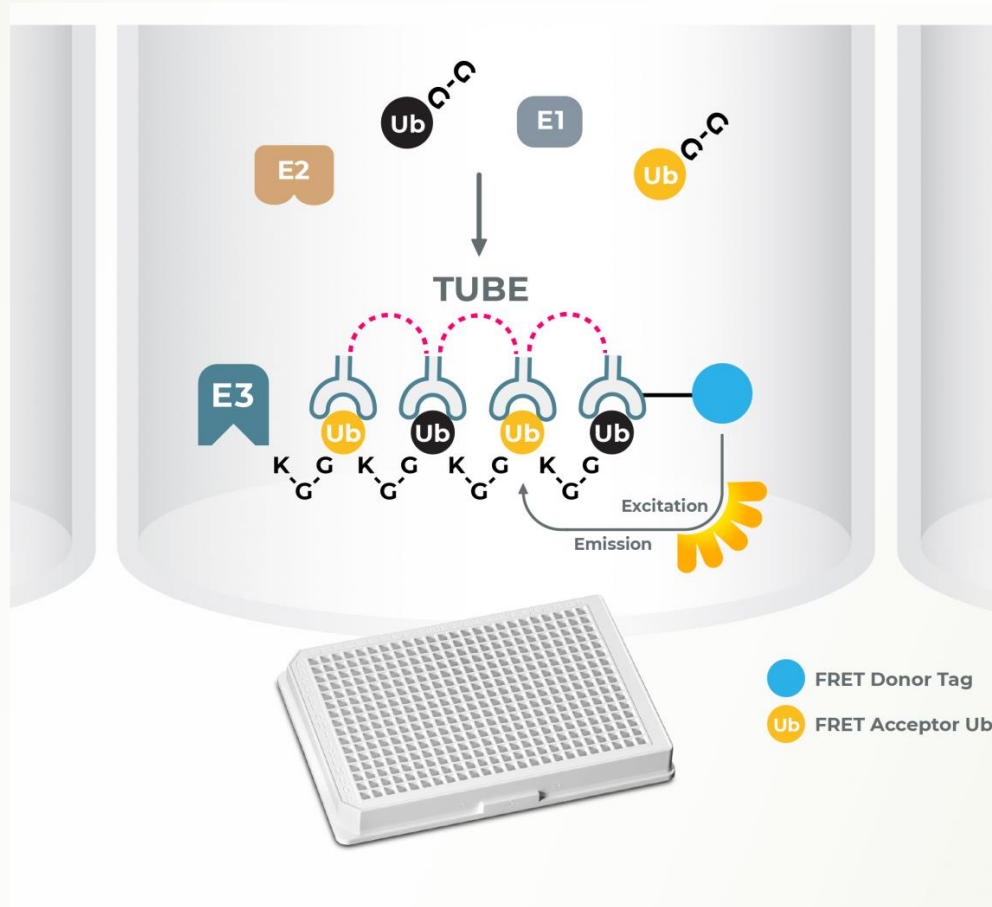


# TUBEs' Role in Molecular Glue Discovery

Study Molecular Glue mediated ubiquitination and degradation simultaneously



# TUBEs Facilitate Discovery of Novel E3 Ligase for KRAS



- ✓ High throughput screening for E3 ligase activators
- ✓ Homogenous assay for library screening
- ✓ Identification of novel E3 ligands
- ✓ SPR / TSA based confirmation and PROTACability



# Novel Application of TUBES, in addition to Ubiquitination Studies

- ✓ Remarkable tool for monitoring PROTAC and molecular glue function
- ✓ HTS of in vivo ubiquitylated proteins
- ✓ Isolation of PROTAC/Mol Glue mediated ubiquitylated substrates from cell lysates
- ✓ Ubiquitin mass spec proteomics bypassing SILAC
- ✓ Perform E3 ligase assays using TR-FRET assays
- ✓ Superior to antibodies, detection by Western blot
- ✓ Imaging tools for In situ detection with fluorescence



# Thank You

We are your partner in UPS, TUBEs, DUBs, E3s, PROTAC, Mol Glue, Protein Expression, CAR-T/Gene therapy and vaccine development

## Contact Us!

Research & Product Inquiries

R&D

[info@lifesensors.com](mailto:info@lifesensors.com)

610-644-8845 (ext 339)

Custom Service & Assays BD

[bd@lifesensors.com](mailto:bd@lifesensors.com)

610-644-8845 (ext 310)