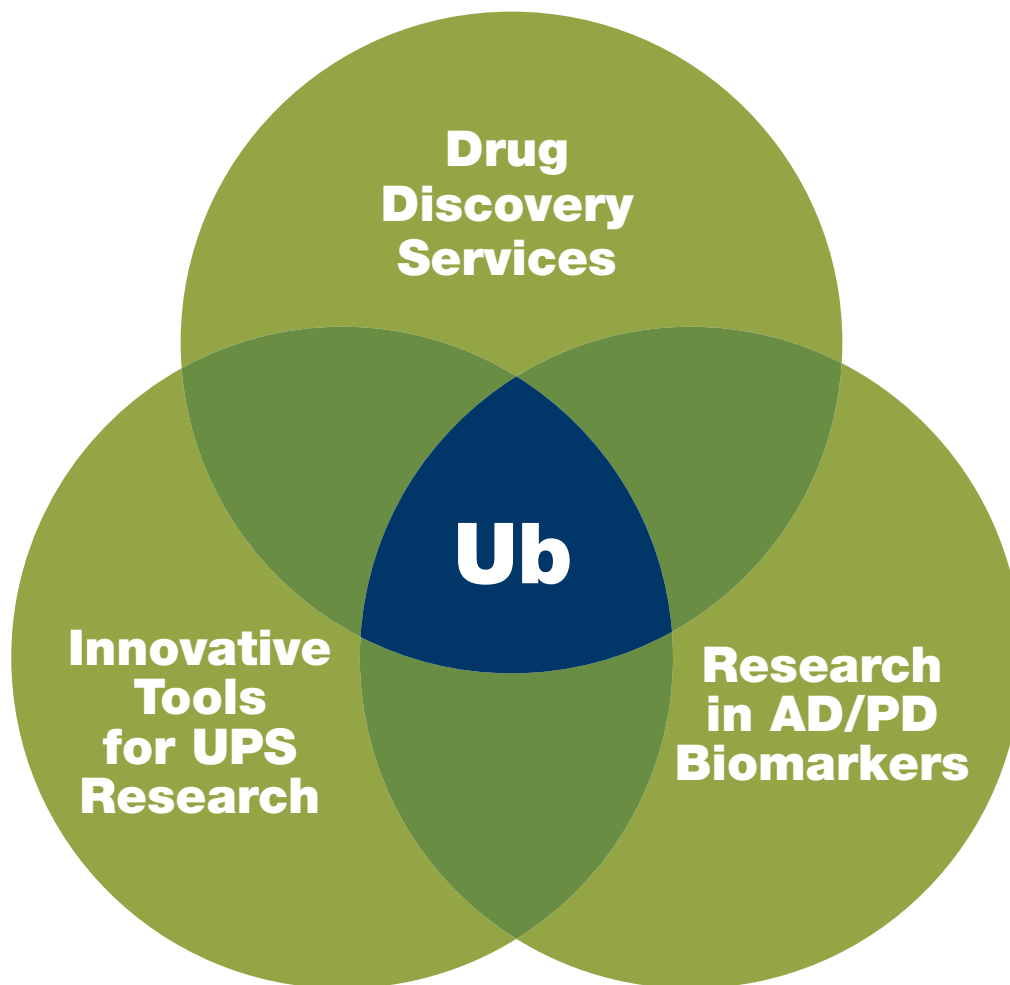


LifeSensors



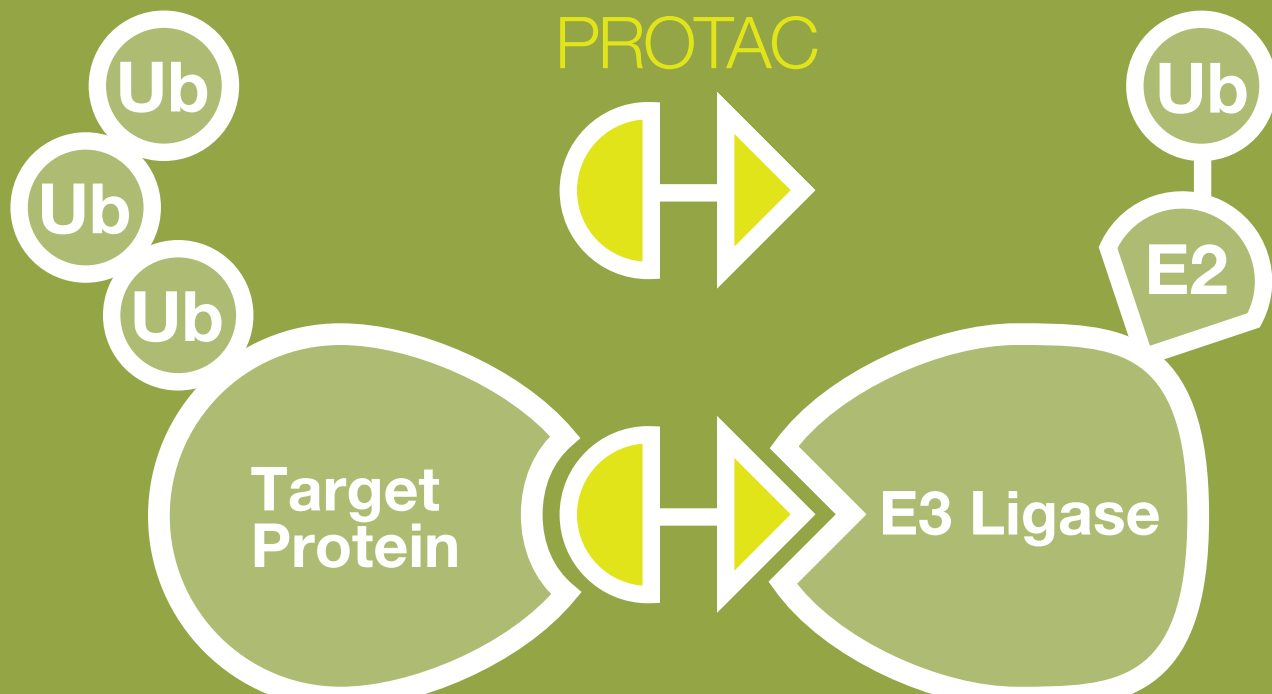
UBIQUITIN

Always our Central Focus



As the ubiquitin leaders for over 20 years,
LifeSensors is a proud sponsor of this conference.

PROTAC Drug Discovery



Let us help you with PROTAC biology.

Ubiquitin Ligase Function and Target Degradation Service

PROTAC compound potency assessment:

PROTAC Compound Profiling • Building SARs • Rank Order Potency

**UBiTest, the most sensitive assay to show
in vivo target ubiquitylation:**

Measure PROTAC activity on 96 or 384 well plate format.

Drug Biomarkers in Cell Culture and Animal Tissue

Assess the state of ubiquitylation in relation to drug efficacy.

NEW!

TUBE-Based Proteomics

Advantages

No need to purify your protein

High quality reproducible data

Complete service starting with your cells or tissue

Excellent support from experiment design to data analysis

Applications

Uncover new substrates for your E3 ligase

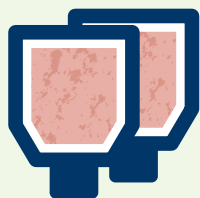
Discover targets for your PROTAC drug

Capture complexed binding partners for your protein

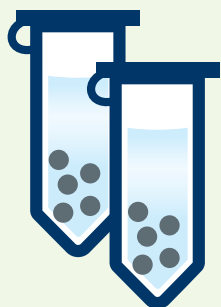
And many more...

Surveying the Ubiquitome has NEVER BEEN EASIER.

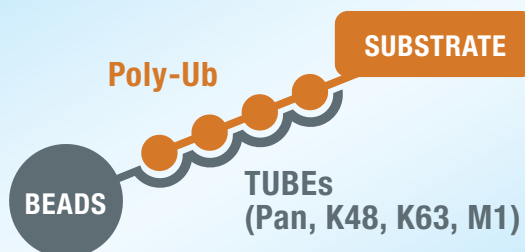
Cell/Tissue Samples



Cell Lysis



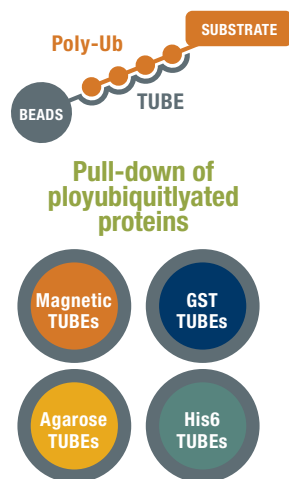
Enrichment of Polyubiquitylated Proteins



Elution & Digestion

TUBE Proteomics

Tandem Ubiquitin Binding Entities



Remarkable TUBE Applications

Biomarkers

Drug Target Validation
DUB Screening
E3 Screening & Profiling

Mass Spectroscopy / Proteomics

TUBE-based M/S
UbiSight Proteomics
Microarrays for Ubiquitomics



"Far Western"
Blot

Pan-Selective
TUBEs (TUBE1)

Linkage Specific
TUBEs (K63, K48, M1)

Histochemistry
Cytochemistry



Ubiquitin Research ONE STOP SHOP

SERVICES

In Vivo Drug Validation

DUB/E3 Screening and Profiling

TUBE-Based Mass Spec

Custom Biomarker Assays

UbiSight Proteomics

Protein Expression

PRODUCTS

TUBEs: Ubiquitin Affinity Matrix

SUMO Fusion Technology

Ubiquitin Enzymes

IQF Diubiquitin for DUB Assays

Ubiquitin Pathway Inhibitors

Ubiquitin Antibodies

Contact LifeSensors to bring your ubiquitin research into focus.

610.644.8845 x339 | info@lifesensors.com

View our full listing of services and products at LifeSensors.com

