

E3 Ligase Profiling & Screening Services

LifeSensors

271 Great Valley Parkway

Malvern PA 19355

Phone: 610-644-8845

bd@lifesensors.com

www.lifesensors.com

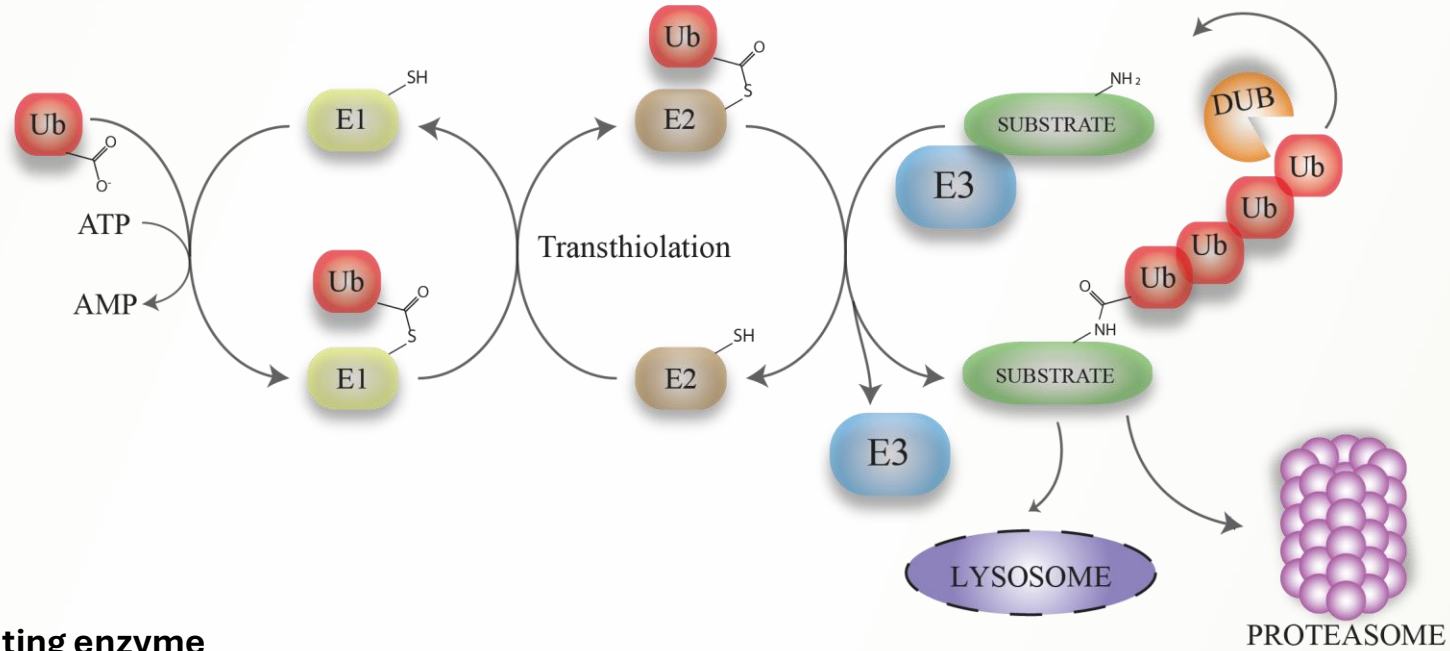
LifeSensors

- Leading Biotech in Ubiquitin Proteasome System (UPS) Drug Discovery
- ~500 Products: Proteins, Ubiquitin Affinity Reagents (TUBEs), Inhibitors, Assays, Kits and Proprietary SUMO Protein Expression Systems
- Drug Discovery, UPS and PROTAC Screening Services
- Profiling Compounds Against Ubiquitin Ligases and De-Ubiquitinases (DUBs)
- Custom Assay Development and Collaborative Research

E3 Ligase Drug Discovery Capabilities

- Expressed and purified ~40 biologically active E3 ligases
- Developed 30 different assays for E3 ligases (auto and substrate ubiquitylation)
- Custom E3 Assay Development and HTS Validation
 - Ability to screen ~650,000 compounds
- E3 and DUB Enzyme selectivity panels for compound profiling
- Determine compound MOA, cellular and target tissue PD markers
 - Enabling technologies based on TUBE applications

Ubiquitin Proteasome System



E1 – Ubiquitin activating enzyme

Requires ATP to attach Ub to E1

E2 – Ubiquitin conjugating enzyme

Transfers Ub from E1 to E3

E3 – Ubiquitin ligases

Transfers Ub to self or substrate

Forms mono-Ub or poly-Ub chains

DUB – Deubiquitinase

Removes mono-Ub or poly-Ub chains

Proteasome – Degrades ubiquitylated proteins

E3 Ligase Drug Screening Overview

➤ **Step One: Assay development, optimization and HTS**

TR-FRET E3 Assay

E3 ELISA Assay

➤ **Step Two: Hit-to-lead optimization**

Working with medicinal chemistry team

Selectivity panel, compound profiling

➤ **Step Three: Validate hits in cellular assays**

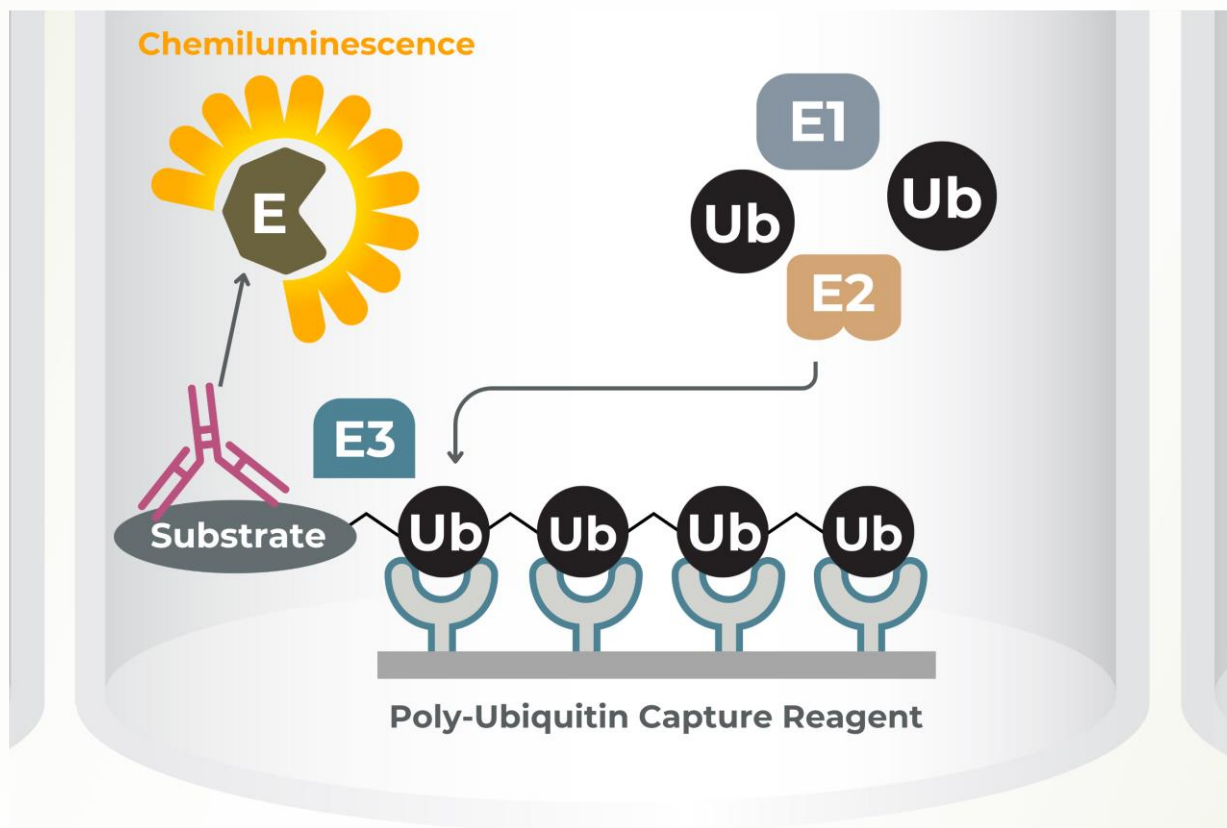
UbiQuant S assay (ELISA / AlphaLISA)

UbiTest (Immunoblot-based assay)

Step One: Assay Development, Optimization and HTS

E3 Ligase ELISA Assays

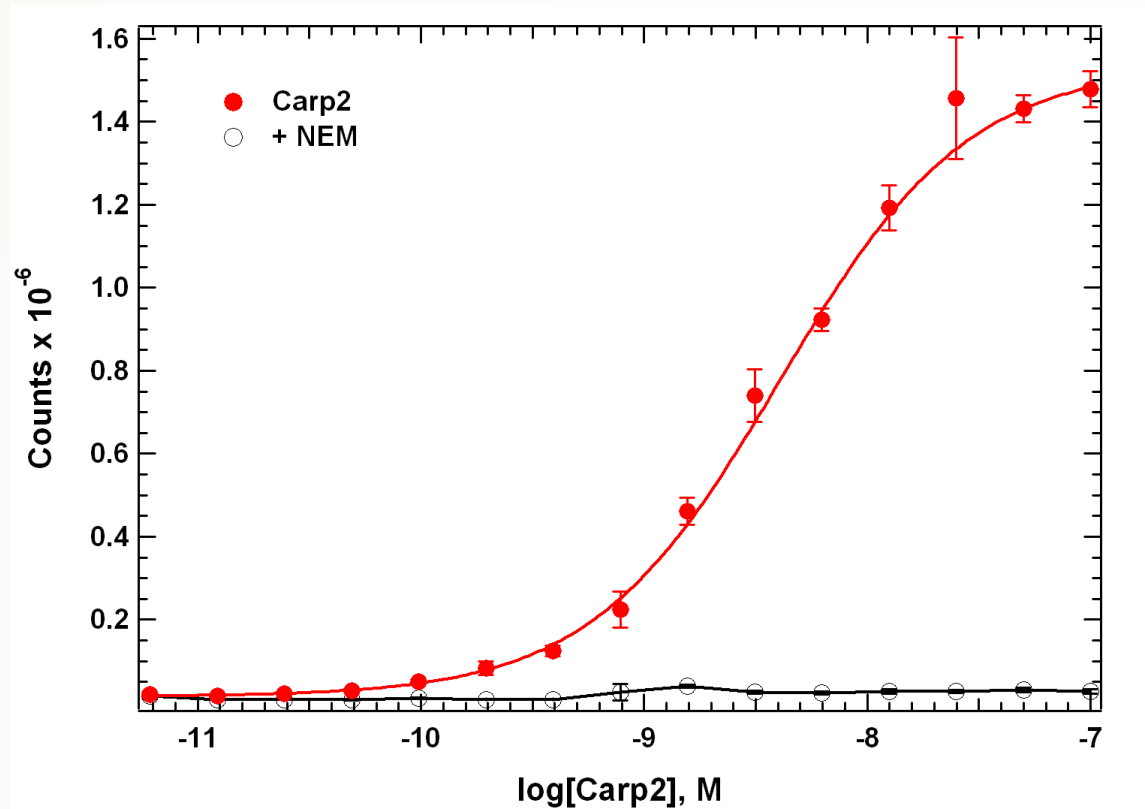
Quantification of E3 ubiquitin ligase activity, employs a proprietary TUBE reagent to capture polyubiquitin chains formed in an E3 ligase dependent manner



- The polyubiquitylated E3 is detected using HRP-conjugated TUBEs.
- Polyubiquitinated substrate is detected using specific antibodies.
- The chemiluminescent signal can be followed over time in a homogenous format
- High-throughput format, ideal for small-molecule screening.

Example of E3 Ligase ELISA Assay

CARP2 E3 Autoubiquitination Assay

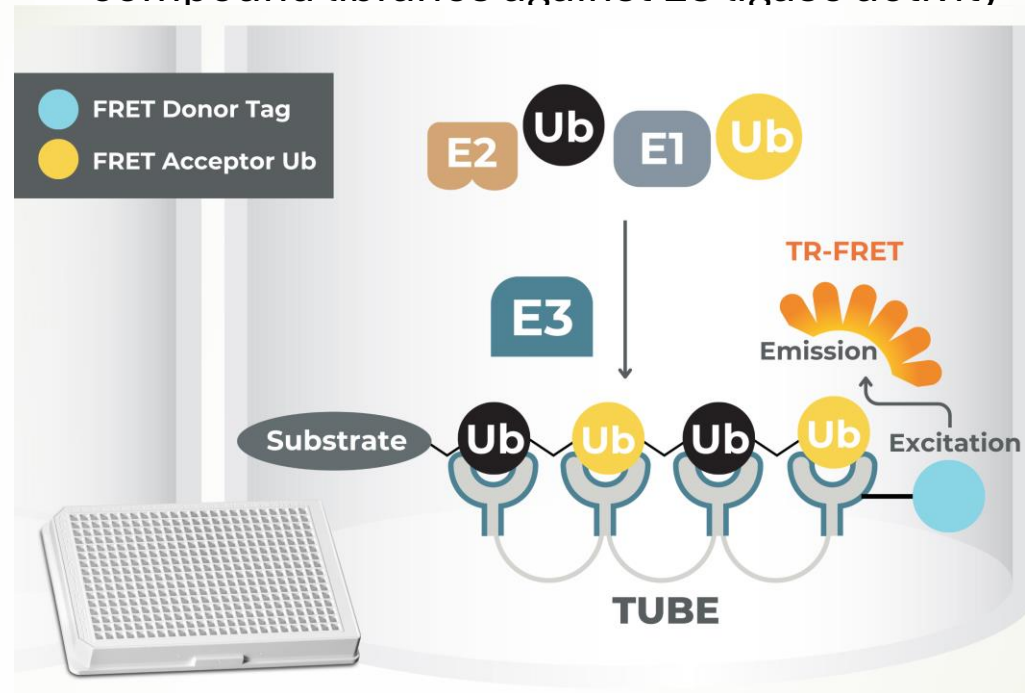


- E3 Dose dependent signal increase
- Robust Assay ($Z' > 0.8$, $S/B > 15$)
- E3 assay inhibited with NEM
- Assay also validated with TAK-243, an E1 inhibitor as positive control for inhibition

Step One: Assay Development, Optimization and HTS

TR-FRET E3 Ligase Assay

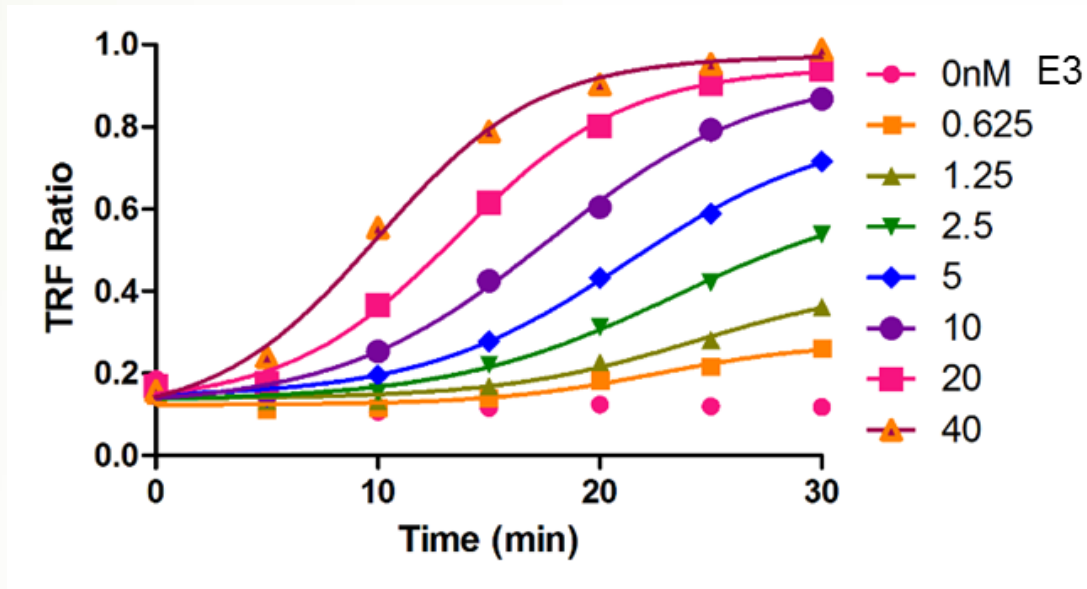
Fluorescence-based high-throughput assay system for screening compound libraries against E3 ligase activity



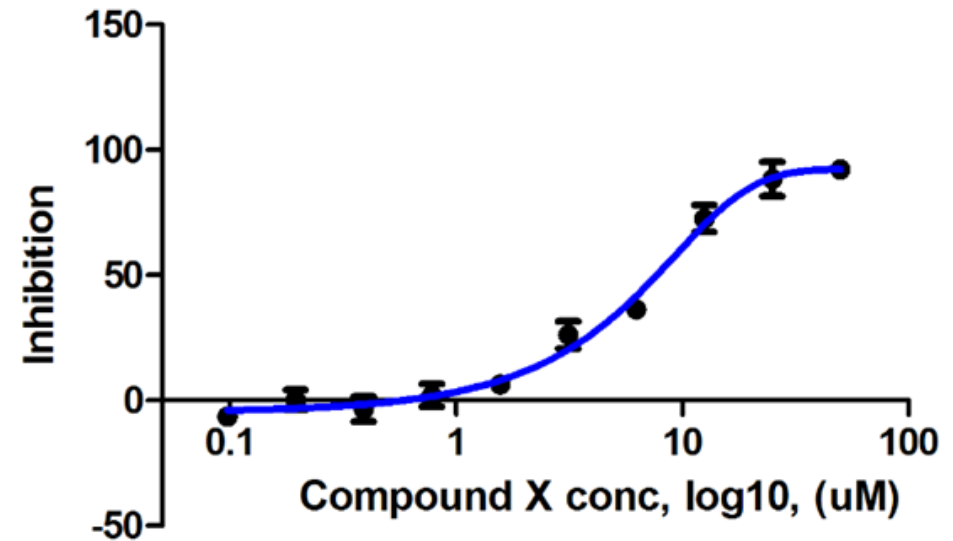
The **TR-FRET E3 Assay** involves Terbium-labeled TUBEs that bind to Fluorescein labelled polyubiquitin chains synthesized by the target E3 ligase. Terbium and Fluorescein are a FRET pair, so polyubiquitin chains containing Fluorescein-labeled ubiquitin yield a FRET signal when bound by a terbium-TUBE. This signal can be monitored over time in a homogenous, high-throughput format, making it ideal for small-molecule screening.

Example of TR-FRET E3 Ligase Assay

E3 Titration Protein X Ubiquitination



Protein X Ubiquitination Inhibition



E3 TR-FRET assay and inhibitor dose response curve

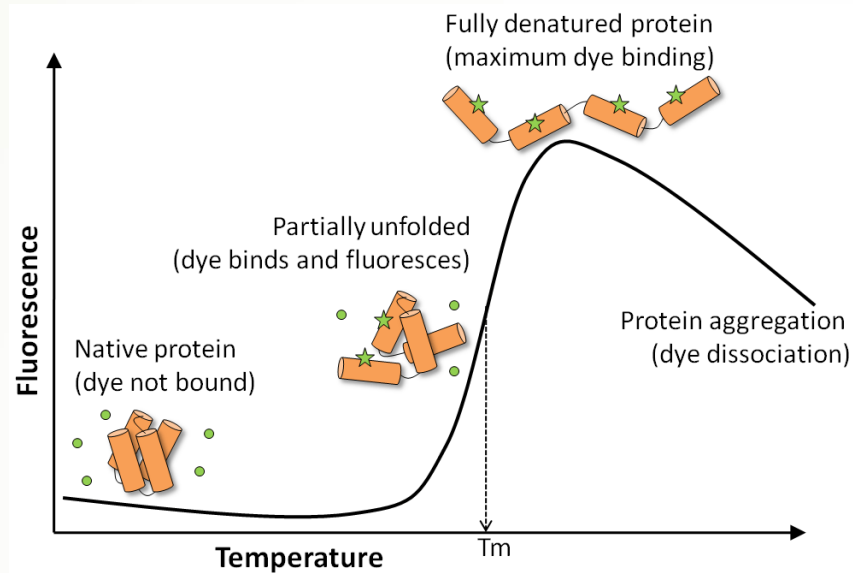
Protein X was used as a substrate for this E3 ligase. After initial TR-FRET high-throughput screening, selected compounds were used to determine IC₅₀ by titration assay.

Step Two: Hit-to-lead optimization

Validation Assays

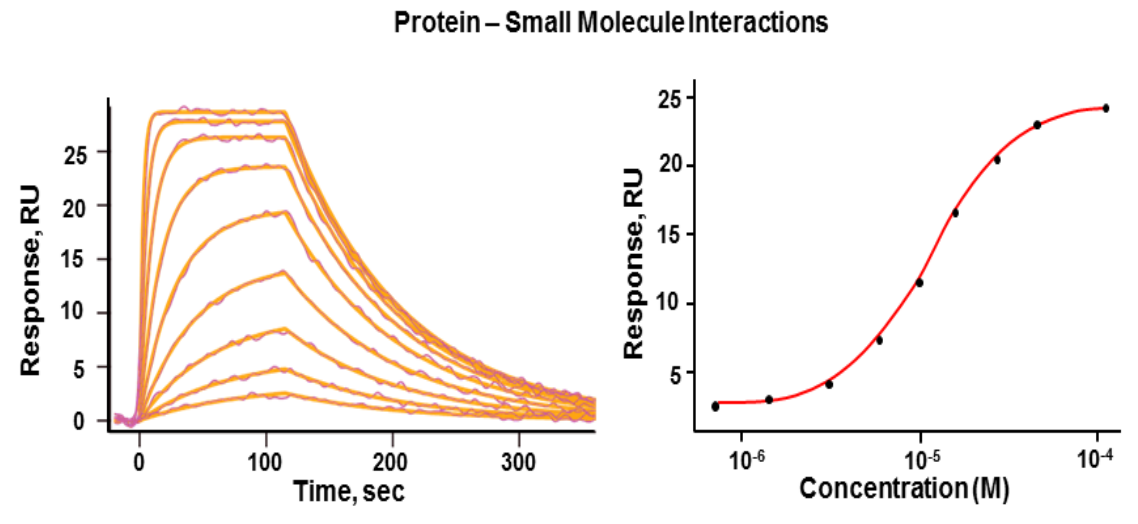
Thermal Shift Assay

HTS assay to detect compound binding to a target



Surface Plasma Resonance

Determination of a small molecule affinity to a target



Step Two: Hit-to-lead optimization

Selectivity Assays

E3 Ligase Panel	Representative E3s
Panel I (5 E3 ligases)	CRBN, CARP2, gp78, CHIP, Nedd4L
Panel II (10 E3 ligases)	CRBN, VHL, HDM2, cIAP2, CARP2, gp78, CHIP, Nedd4L, Praja1, Cbl-b
Panel III (29 E3 ligases, includes E3 from panel I as well)	CRBN, VHL, Hdm2, RNF4, CARP2, TRIM32, TRIM47, Cbl-b, c-Cbl, cIAP2, IDOL, SIAH, MURF1, MURF2, MURF3, Praja1, TRAF6, Parkin, E6-AP, Itch, Nedd4L, WWP1, WWP2, MARCH5, Hrd1, gp78, CHIP, RNF114, Nedd4

Each ligase assay has been validated in TR-FRET assays regarding E2 pairing.

LifeSensors profiles inhibitory or activation properties of every compound in Panel I followed by Panel II.

Step Two: Hit-to-lead optimization

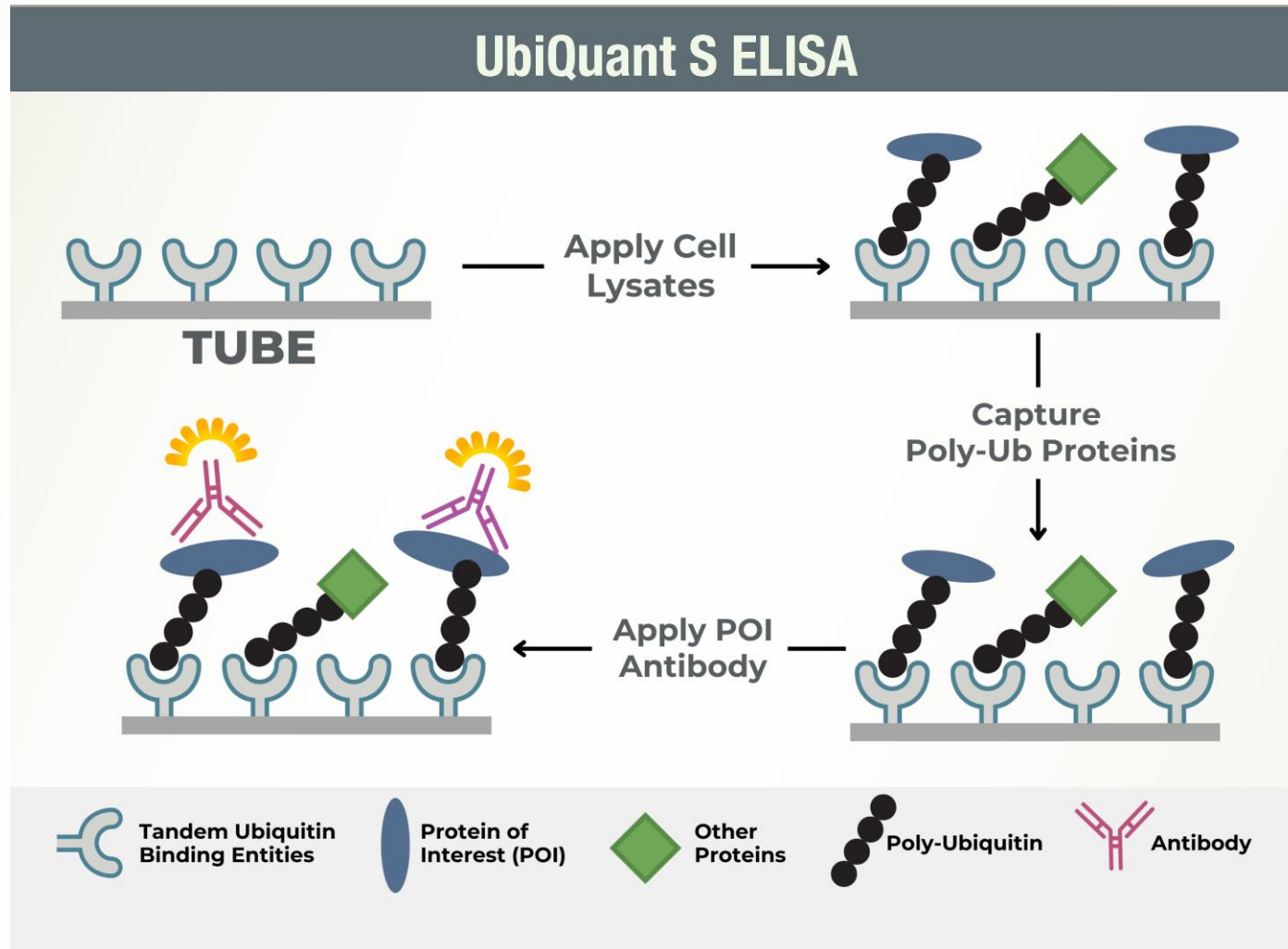
Mechanistic Validation Assays

**Secondary screens to deconvolute hits from E3 screening
(eliminating compounds that affect E1-E2 conjugation)**

- E3 Lite - Measures E3 activity
- E1 Lite - Measures E1 activity
- E1/E2 transfer - Measures transfer between E1 to E2
- E2 Profiling and Selection - Finds the best E2 for your E3

Step Three: Validate hits in cellular ubiquitination assays

Enables accurate determination of **cellular substrate (POI) ubiquitination** for monitoring the effects of various treatments



Step Three: Validate hits in cellular assays

TUBE-based Mass Spec Ubiquitin Proteomics

- **TUBE-based proteomics to identify ubiquitylation patterns specific to drug treatment**
 - **Optimized for cell and tissue lysates**
 - **Customer provides cell pellets, we do the rest**
- **Superior to Di-Gly proteomics method**
- **Assess specificity of E3 ligands, inhibitors, PROTACs and Molecular Glues**
- **Identify polyubiquitylation site(s) (number & position) on the protein sequence**
- **Fee for service model, defined milestone-based agreement**

Identification E3 ligase modulators for Clients:

Example #1

➤ E3 ligase X: Assay development, validation and HTS

50K small molecule library screen using TR-FRET E3 Assay

1600: number of primary hits, $Z' > 0.5$

64: number of confirmed hits with selectivity

10: number of compounds with IC50s sub micromolar to nM

➤ Step Two: Hit-to-lead optimization

Hit expansion (with medicinal chemistry team)

Extended selectivity panel, compound profiling

➤ Step Three: Confirm hits in cellular assays

10 hits transferred to client for cellular validation

Identification E3 ligase modulators for Clients:

Example #2

- **E3 ligase X: Assay development, validation and HTS**

 - Client's compounds screened using TR-FRET E3 Assay

 - 10: number of confirmed compounds with selectivity

 - 10: number of compounds with IC50s sub micromolar to nM

- **Step Two: Hit-to-lead optimization**

 - Hit expansion (with medicinal chemistry team)

 - Extended selectivity panel, compound profiling

- **Step Three: Confirm hits in cellular assays**

 - Confirmation through [Ubiquitin Mass Spec Proteomics](#)

Identification of E3 ligase target for Molecular Glue degraders:

Example #3

- **E3 ligase X degrading target Y: Assay development, validation**

 - Validation of target degradation and ubiquitination in cells (kinetics)

 - Rescue of degradation using Proteasome/Lysosome inhibitors

 - Determine the optimal dose and time needed to robustly ubiquitinate target

- **Step Two: Mass Spec Proteomics for E3 identification**

 - Pull down target protein ubiquitination complex from cells treated with degrader

 - Perform proteomics to identify interacting E3 ligases

- **Step Three: Validate hits in in vitro and cellular assays**

 - Use recombinant E3s to confirm molecular glue mediated ubiquitination of target in vitro

 - Validate the role of E3 in cells using CRISPR/Cas knock-out system

E3 Ligase Screening & Profiling Services

- **We help customer discover E3 ligase ligands, inhibitors and activators**
 - **Express & purify biologically active E3 Ligases and substrates**
 - **Develop and optimize HTS assay for E3 ligase**
 - **Screen in house libraries or customer libraries at LifeSensors**
 - **Confirmation and counter screening to eliminate off-target compounds**
 - **Biophysical and biochemical assay development for target engagement**
- **Cell-based assays to determine target engagement by compound**
- **All IP and data belong to the customer**
- **Work performed under CDA and Master Service Agreement**
- **Fee for service model, defined milestone-based agreement**

Contact Us!

We are your partner for E3 Ligase drug discovery

Contact Information

Research & Product Inquiries

R&D

info@lifesensors.com

610-644-8845 (ext 339)

Custom Service & Assays BD

bd@lifesensors.com

610-644-8845 (ext 354)