

# Accelerating PROTAC and Mol Glue Drug Discovery

## Unleashing the Potential of TUBE Embedded Microtiter Plate Technology

**LifeSensors Inc.**

271 Great Valley Parkway

Malvern PA 19355

Phone: 610-644-8845 x 310

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

[www.lifesensors.com](http://www.lifesensors.com)

# LifeSensors Leadership in UPS Drug Discovery



Leading Biotech in PROTAC and Mol Glue Drug Discovery



500 Products supporting UPS research



E3 ligase and DUB compounds profiling



Discovery services for new E3 ligands, PROTACs and Mol Glues

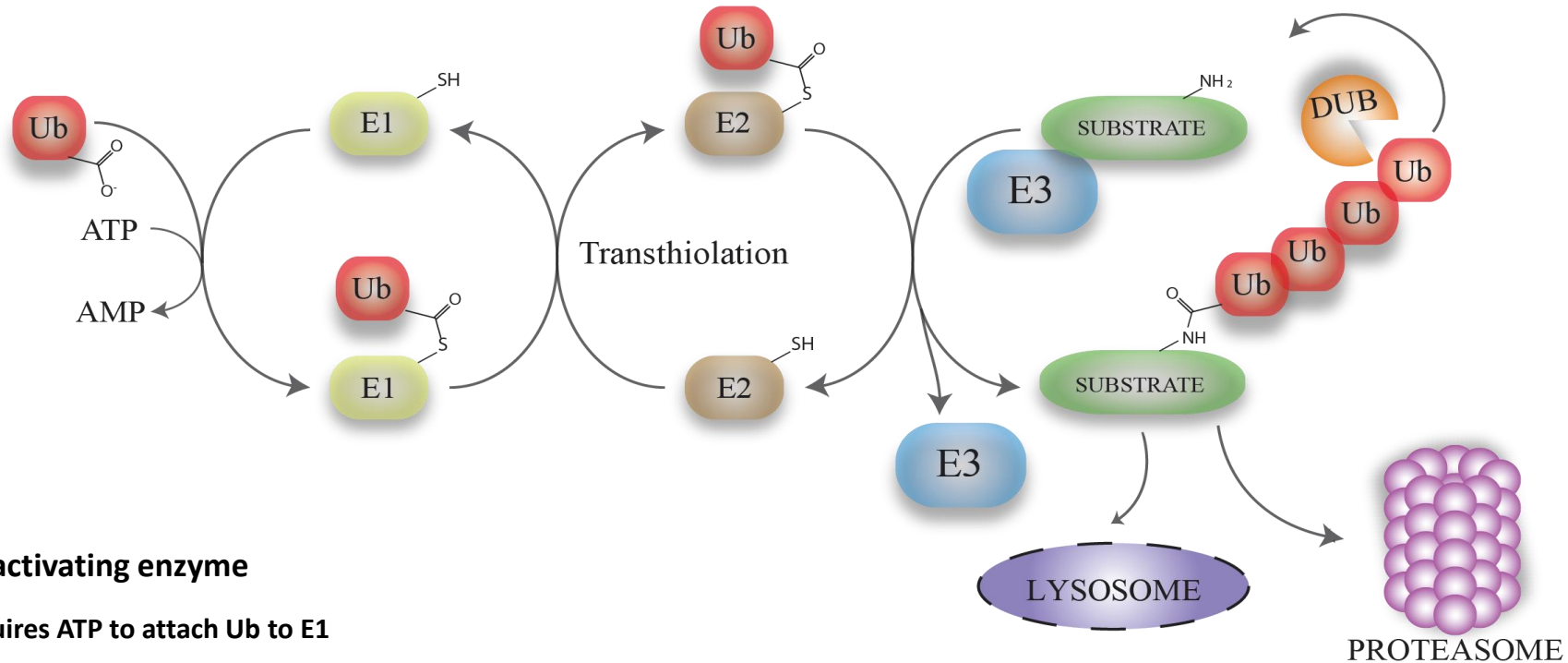


Collaborative and Independent Research

Malvern, Pennsylvania, USA



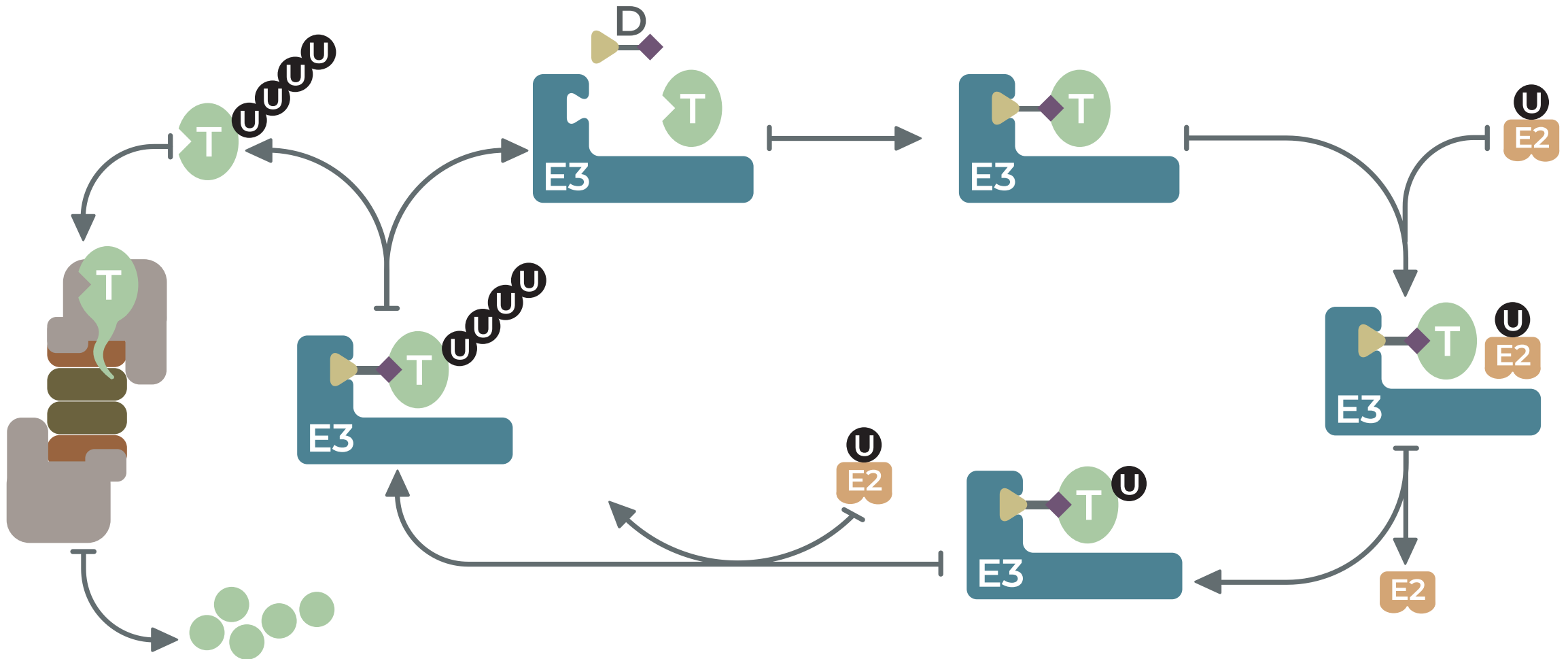
# Understanding Ubiquitin Proteasome System in Drug Discovery



- 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**

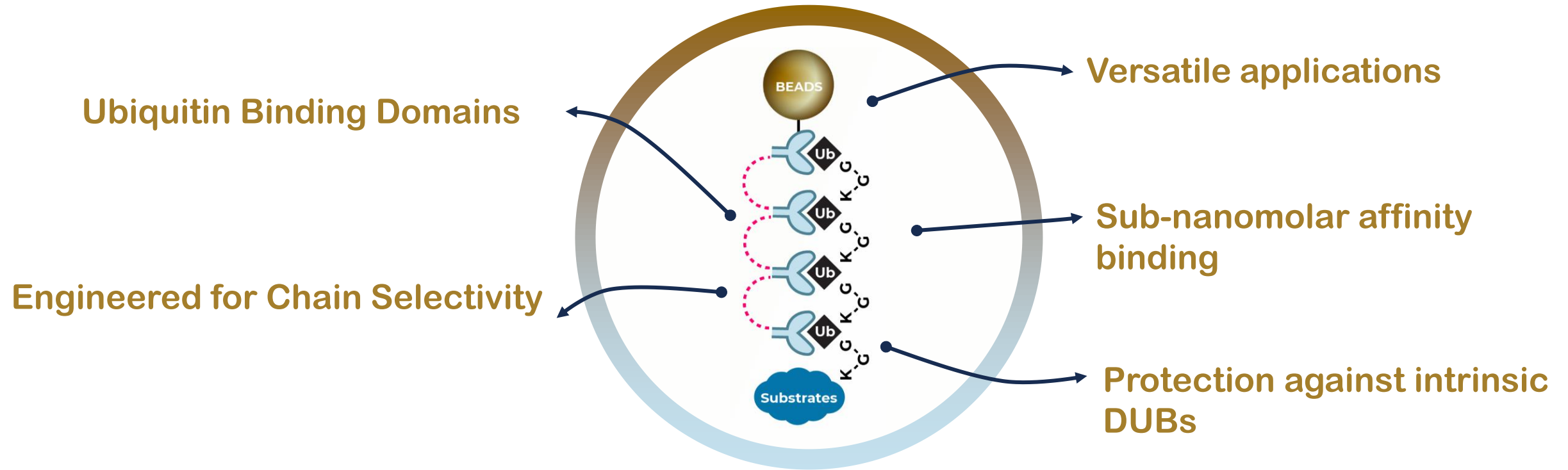
# PROTAC<sup>®</sup>s - Essential Catalytic Activators



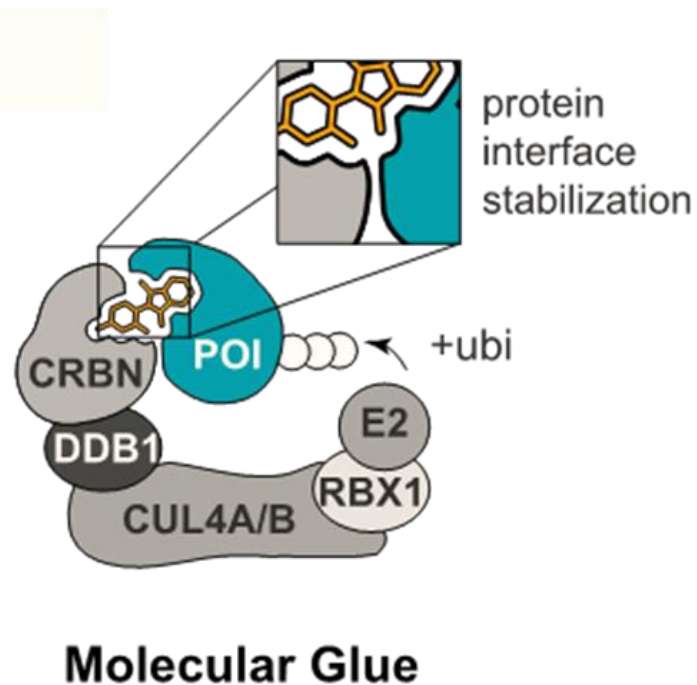
“ LifeSensors makes it easy to study Ubiquitination ”

# TANDEM UBIQUITIN BINDING ENTITIES

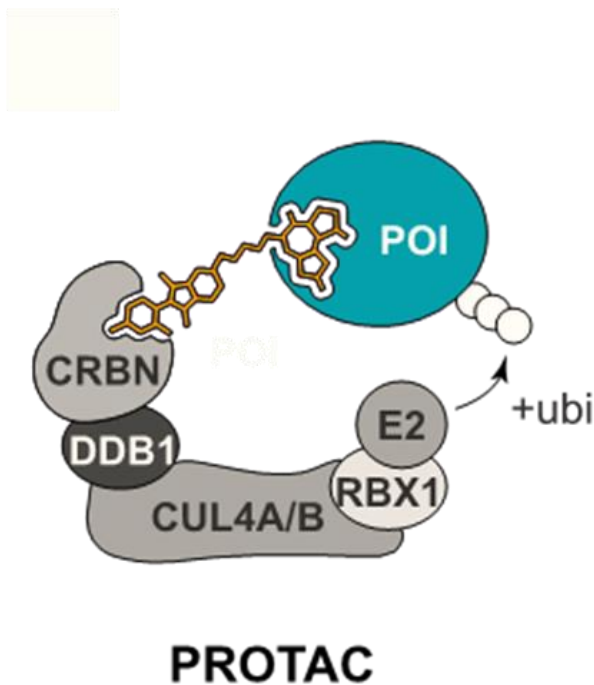
## TUBE<sub>s</sub>: A Versatile Tool in E3 Ligase Assays



# Targeted Protein Degradation



Immense promise for molecular glues challenged by **lack of reliable approaches** to rationally design them

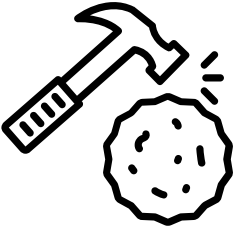


Expectations for PROTACs are very high, but recent clinical data suggest developmental challenges

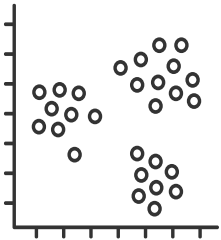
# TUBEs address critical TPD challenges



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



**Ternary complex rigidity**



**Ligand dependent ubiquitination, true functional HTS**

# LifeSensors Approaches Standout

## Traditional Approaches

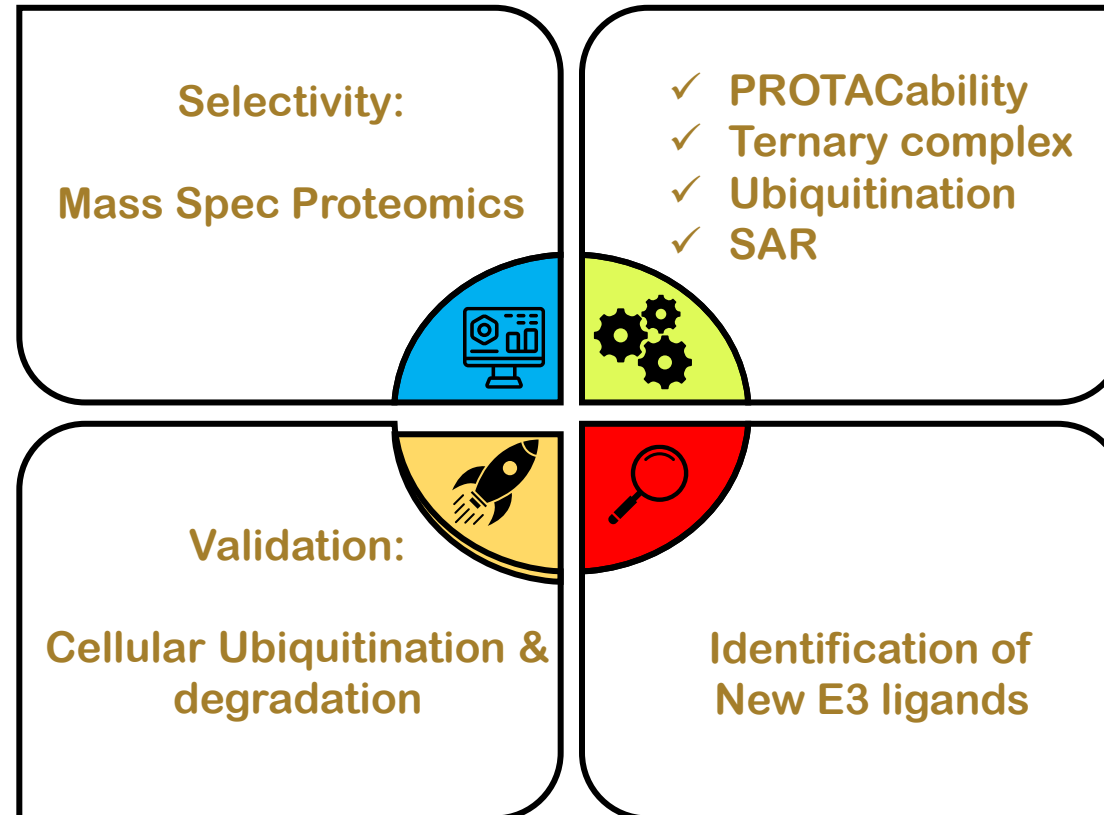
- **Reporter gene assays:**  
External tags (Off target ubiquitination)
- **Proximity Ligand Assays**  
protein's intrinsic factors - overlooked
- **Western Blotting**  
Low throughput and Irreproducible

## LifeSensors Approaches

- **Invitro Ubiquitination assays :**  
Ligand mediated ubiquitination, fast and functional approach
- **Cellular degradation assays:**  
No reporter tags required, applicable to clinical and preclinical studies "Biomarker"
- **HTS screening**  
Native lysines playing a role

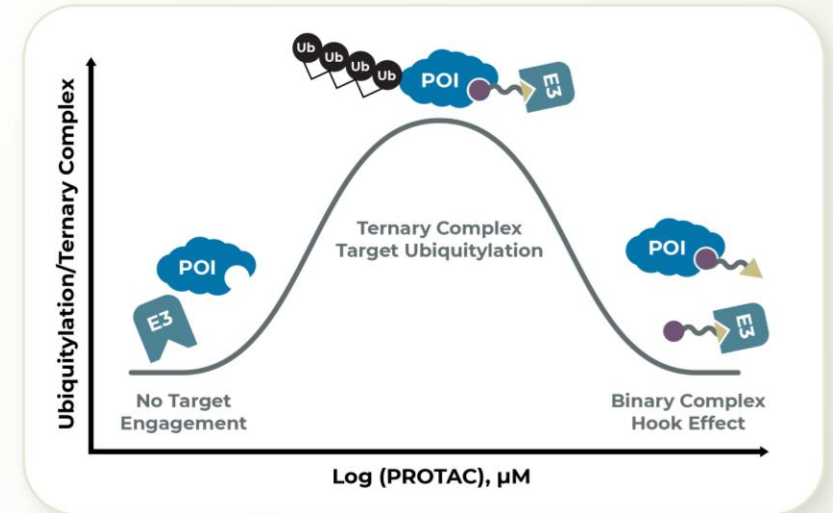
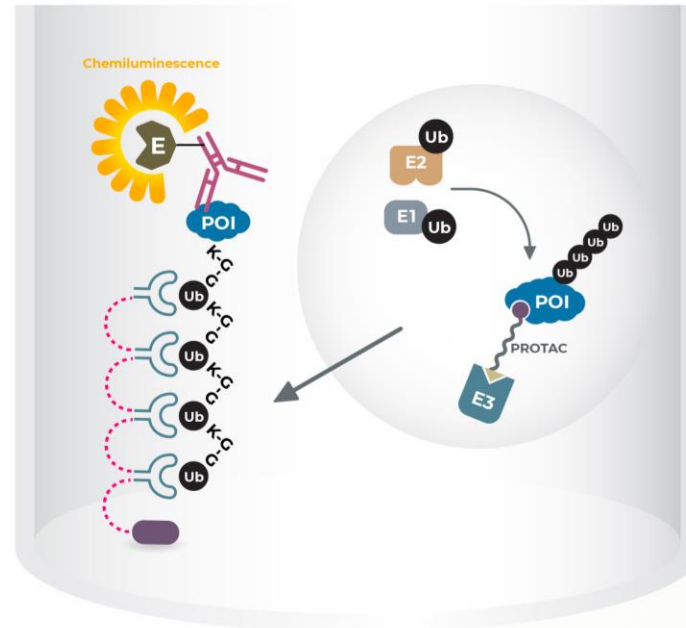
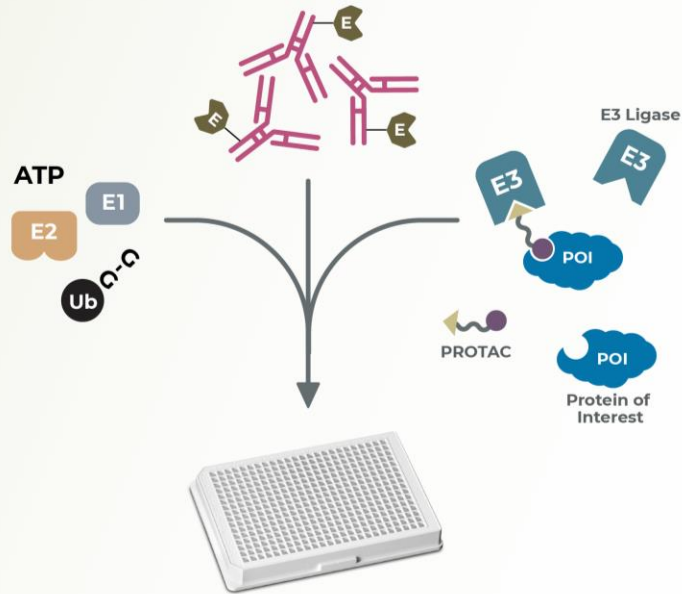


# LifeSensors Approaches for Targeted Protein Degradation



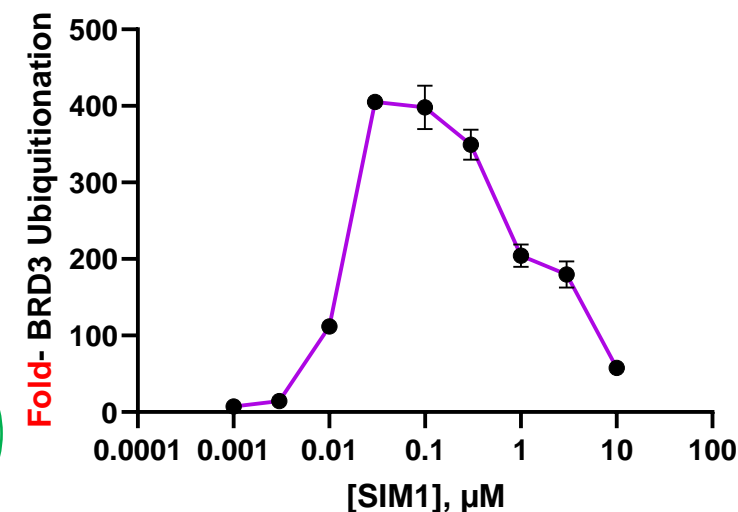
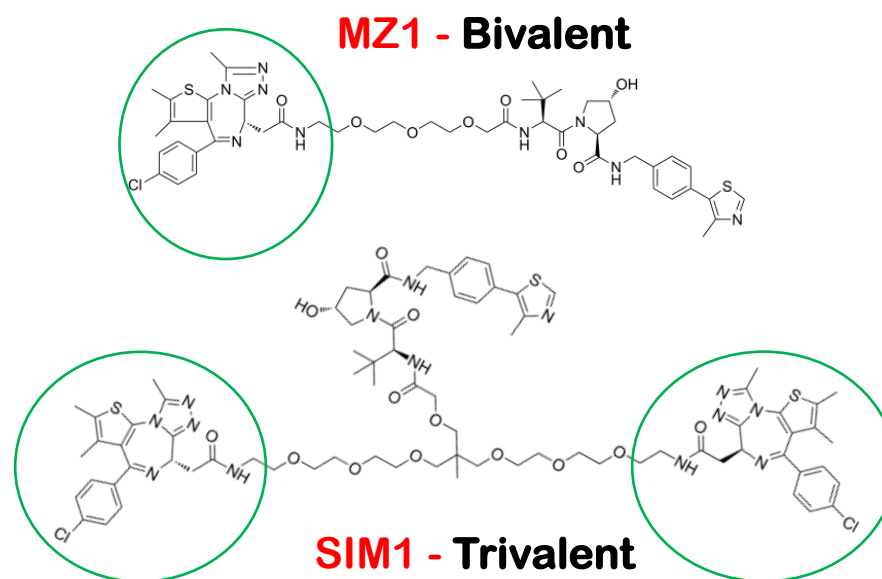
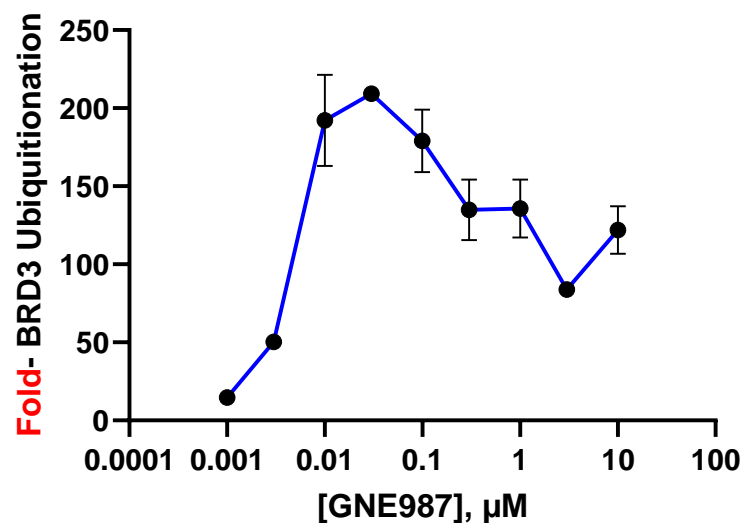
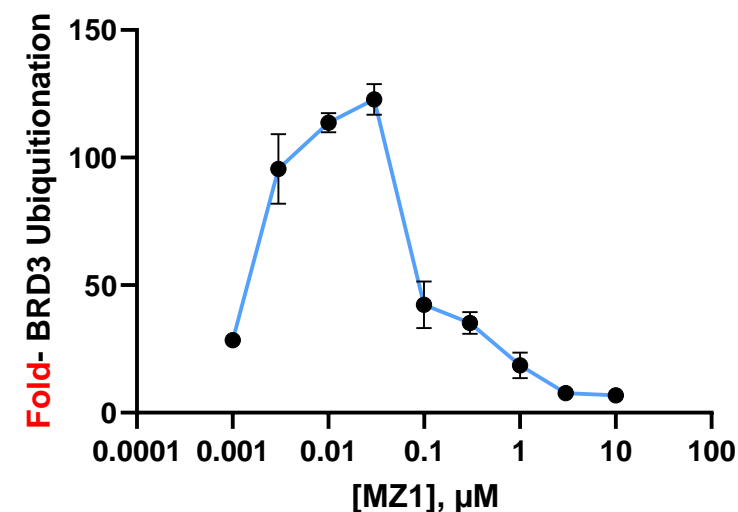
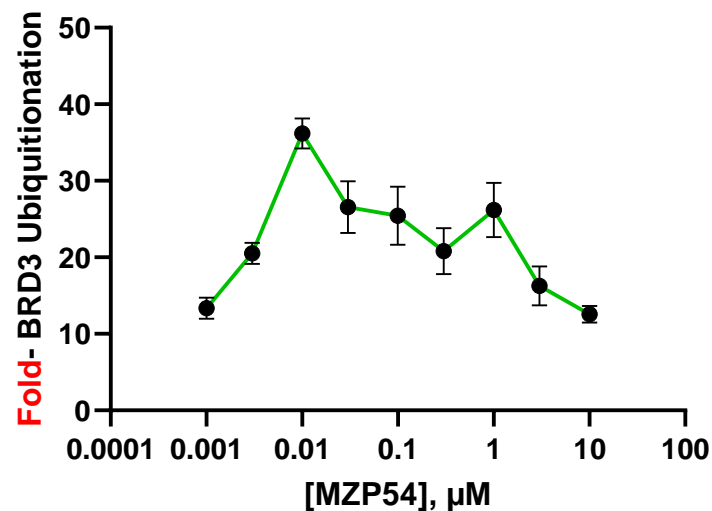
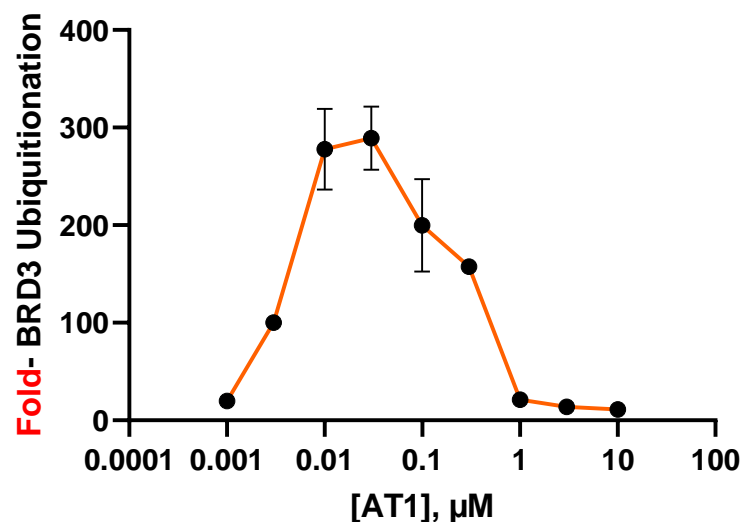
# HTS- In Vitro Biochemical Assay

To study functional ternary complex and PROTACability

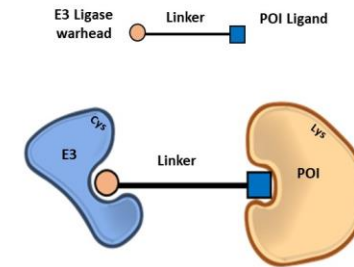
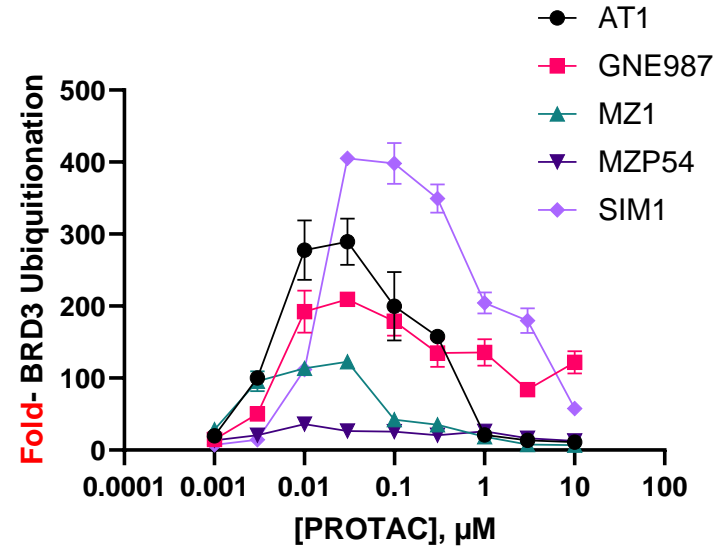
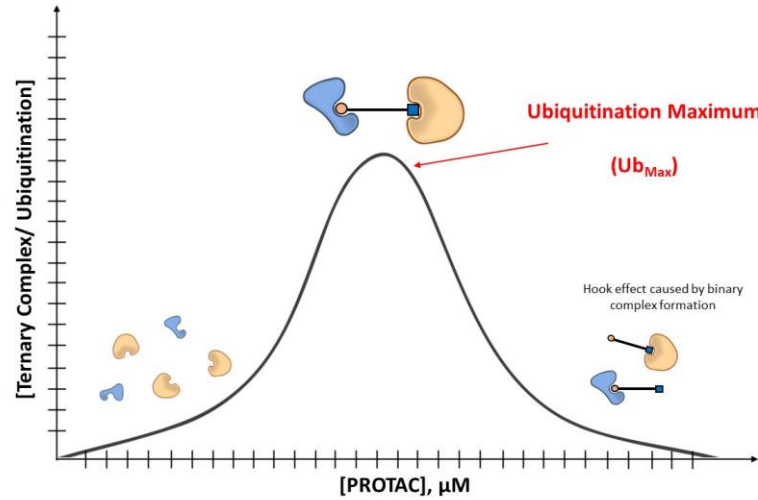


TUBE Capture & PROTAC Mediated Ubiquitination of POI Detection

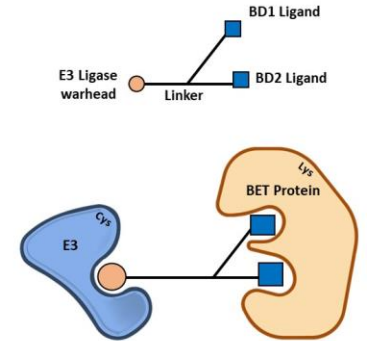
# VHL Bromodomain PROTACs



# VHL Bromodomain degraders



**MZ1 - Bivalent**

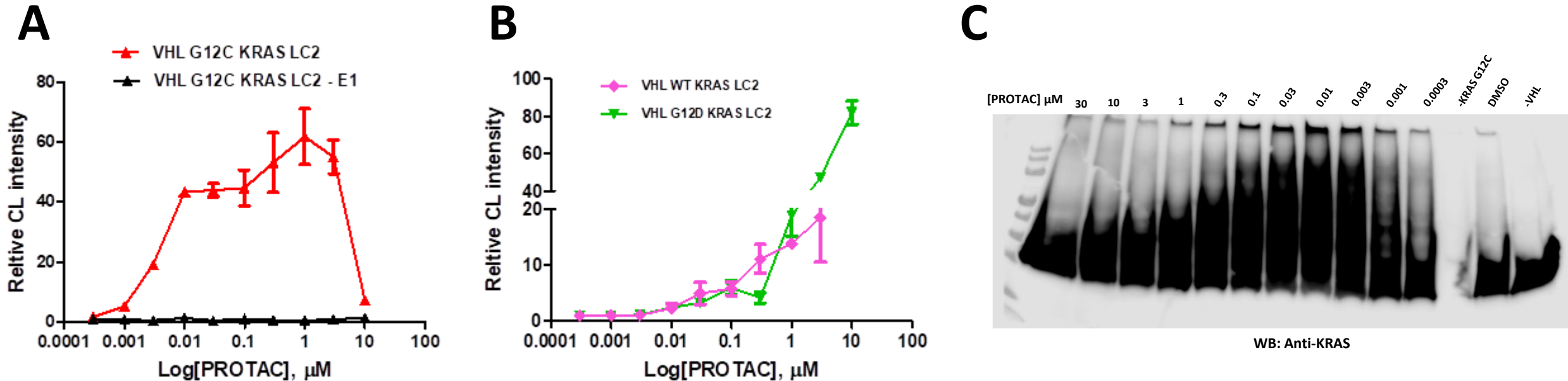


**SIM1 - Trivalent**

PROTAC	Cellular EC <sub>50</sub>	Area Under the curve
AT1	~100.0 nM	244.4
GNE987	9.9 nM	1084.0
MZ1	189.9 nM	112.4
MZP54	~10.0 nM	167.3
SIM1	1-10 nM	1519.0

# K-RAS PROTACs

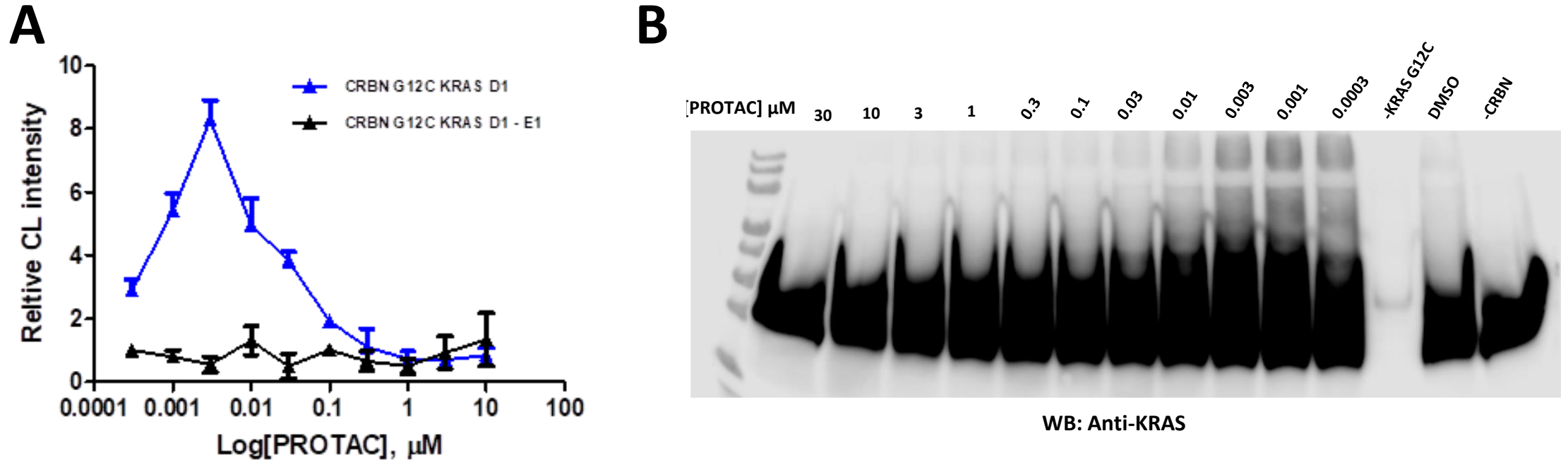
## VHL PROTAC mediated in vitro ubiquitination



*In vitro* ubiquitination assay with KRAS G12C VHL based degraders: (A & B) VHL-based PROTAC LC2 in a dose response study to monitor PROTAC mediated ubiquitination of KRAS G12C, G12D and wildtype. CL intensities plotted in response to  $\frac{1}{2}$  log dose response demonstrates PROTAC mediated ubiquitination. (C) Western blot analysis to confirm PROTAC mediated ubiquitination via characteristic poly-ubiquitination smears with anti-KRAS antibody.

# K-RAS PROTACs

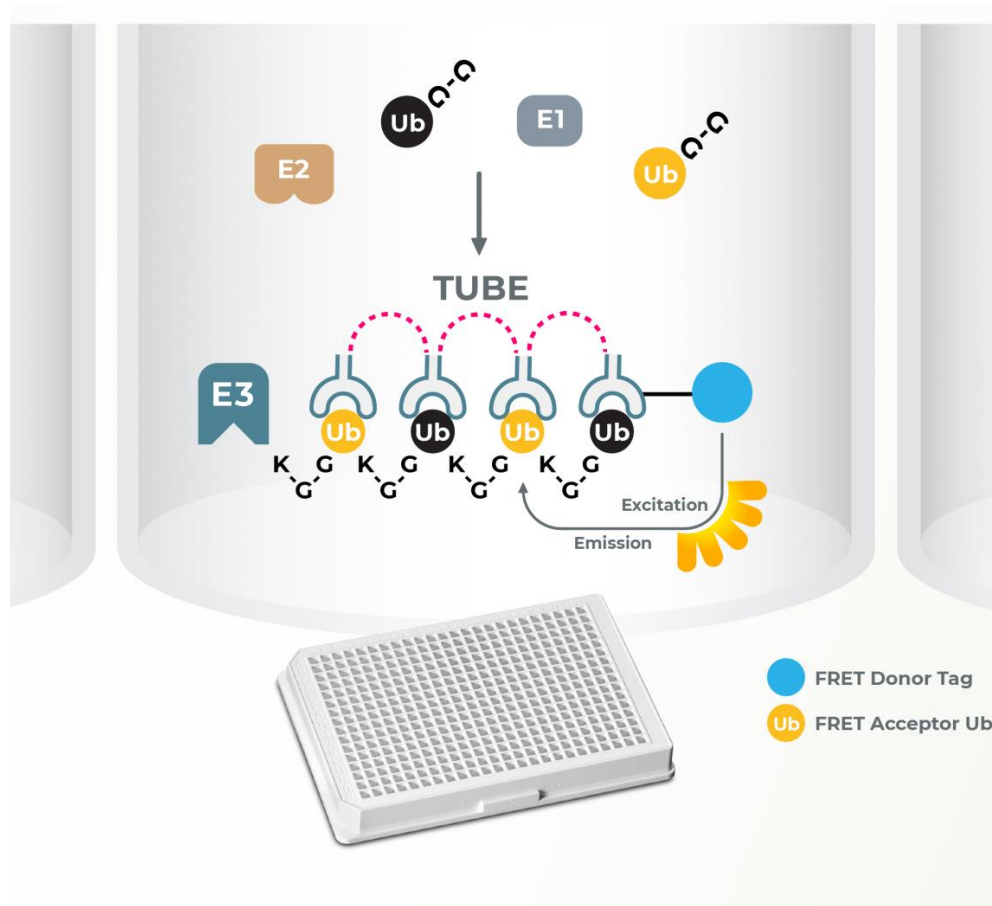
## CRBN PROTAC mediated in vitro ubiquitination



*In vitro* ubiquitination assay with KRAS G12C CRBN based degraders: ((A) CRBN-based PROTAC degrader 1 (compound 518) in a dose response study to monitor PROTAC mediated ubiquitination of KRAS G12C. CL intensities plotted in response to  $\frac{1}{2}$  log dose response demonstrates PROTAC mediated ubiquitination. (B) Western blot analysis to confirm PROTAC mediated ubiquitination via characteristic poly-ubiquitination smears with anti-KRAS antibody.



# Novel E3 Ligase for KRAS

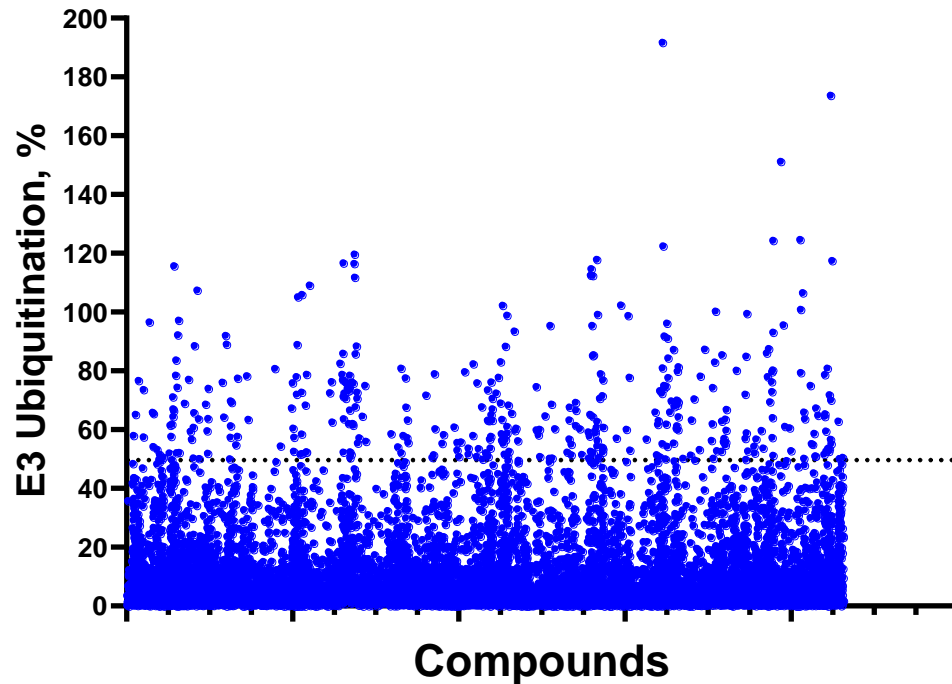


- ✓ **High throughput screening for E3 ligase activators**
- ✓ **Homogenous assay for library screening**
- ✓ **Identification of novel E3 ligands**
- ✓ **SPR / TSA based confirmation and PROTACability**

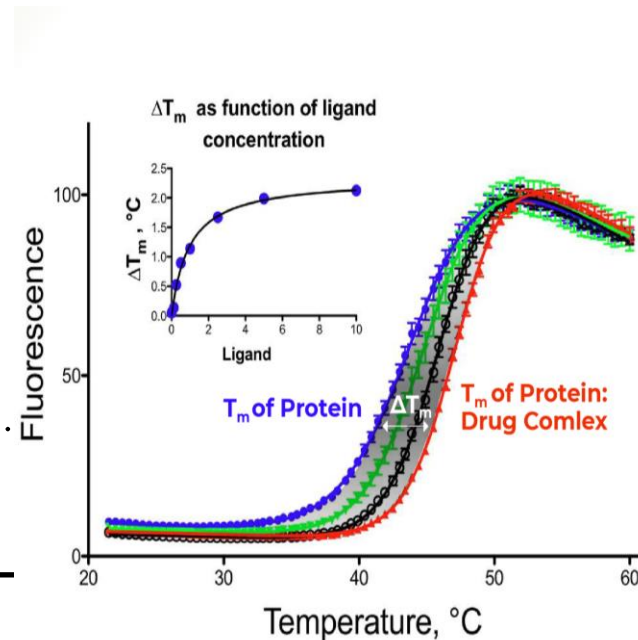
# Discovery of Novel E3 Ligase for KRAS

## Superior Properties as compared to VHL and CBRN

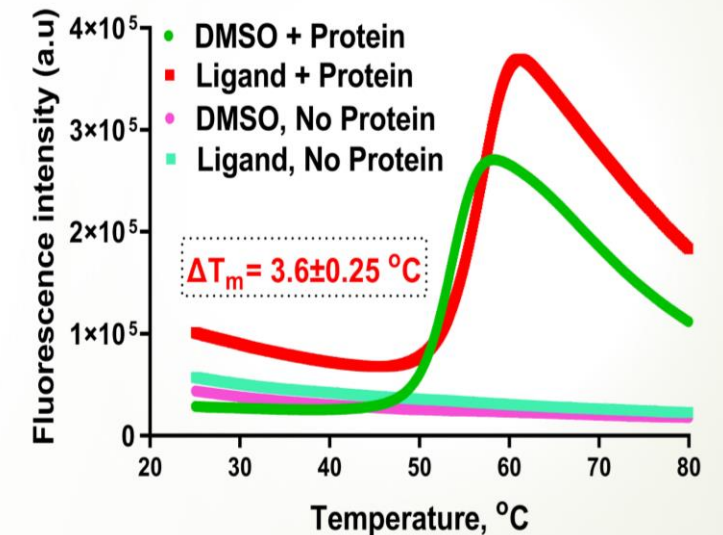
50K compound library screening



Thermal Shift Assay Validation

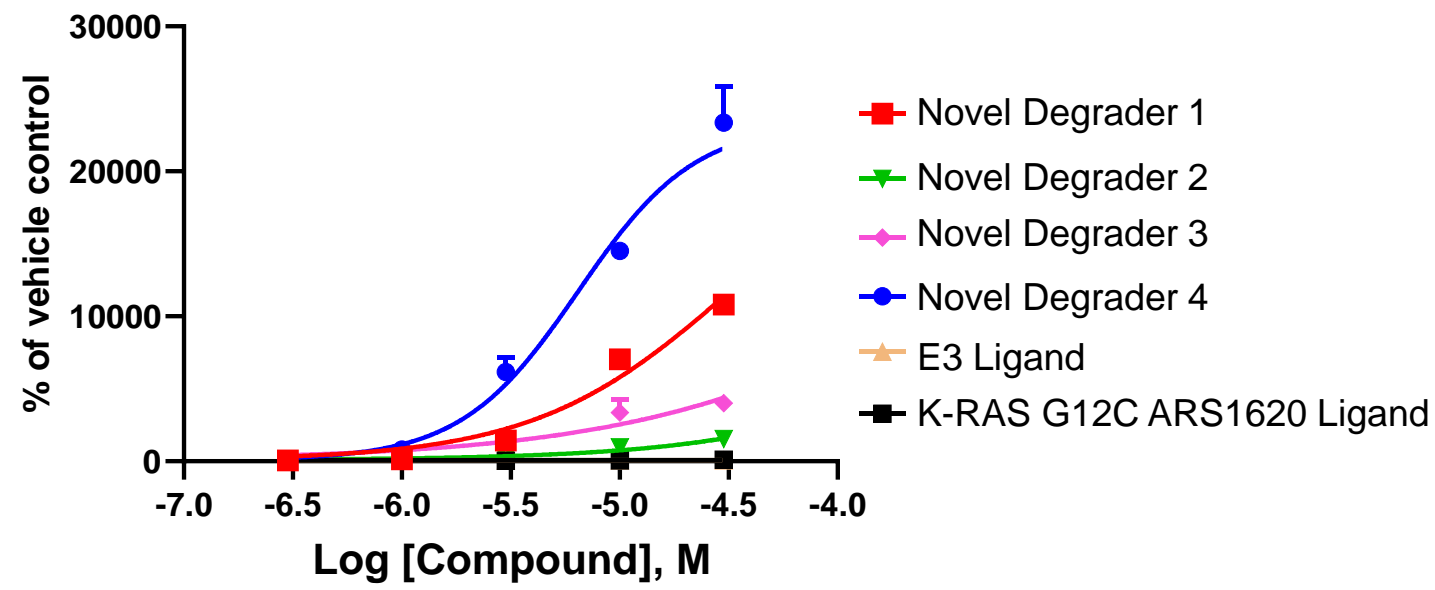


Thermal Shift Assay Validation: Ligand Binding

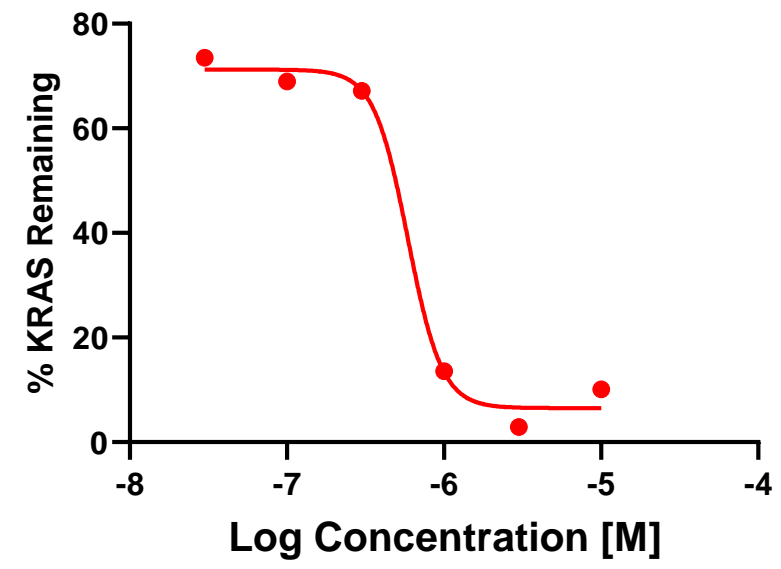
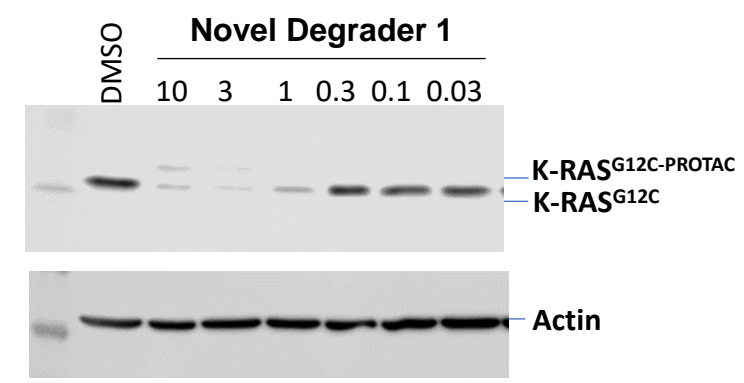




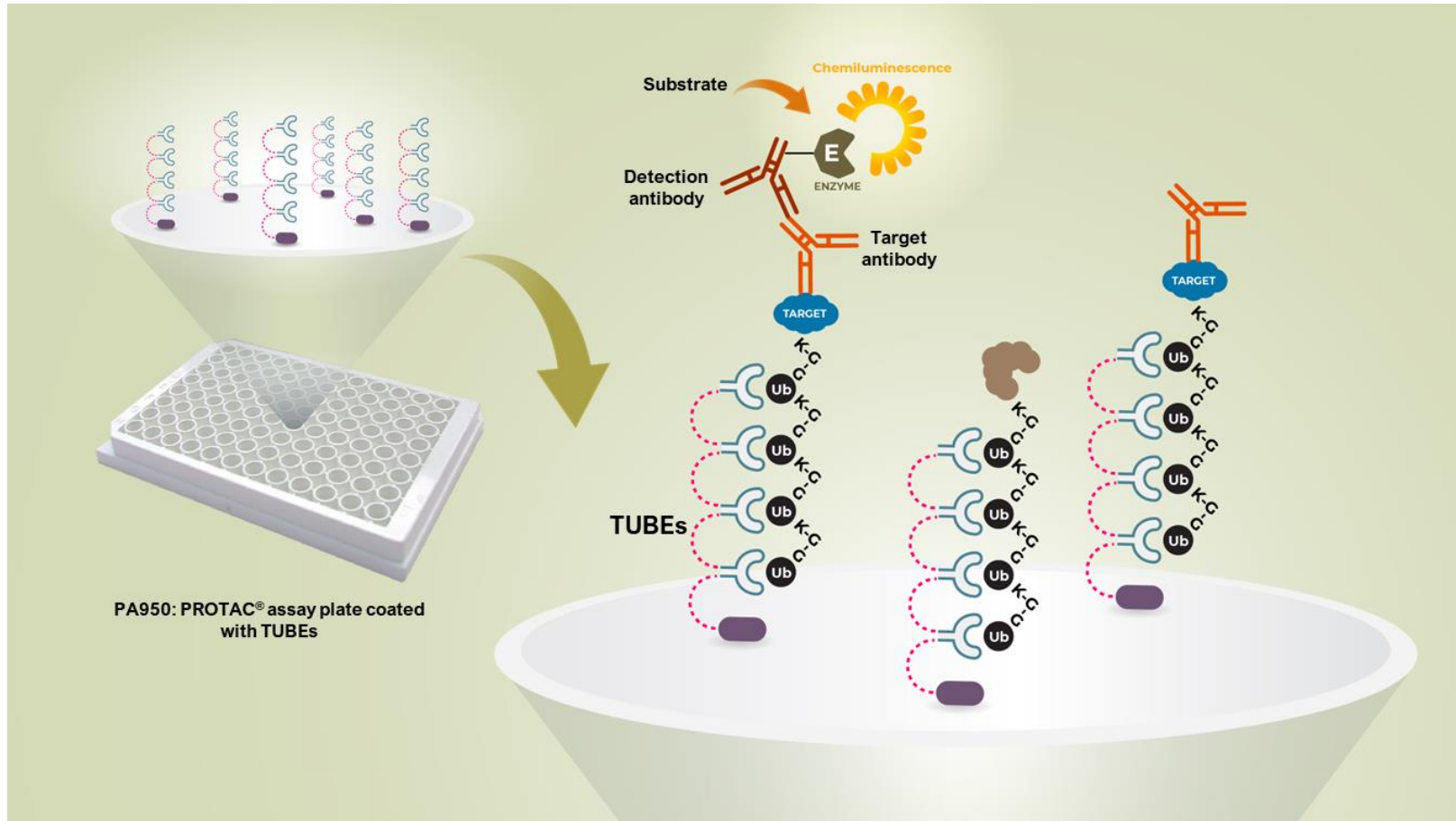
# Novel E3 Ligand with ARS-1620 for KRAS Degradation



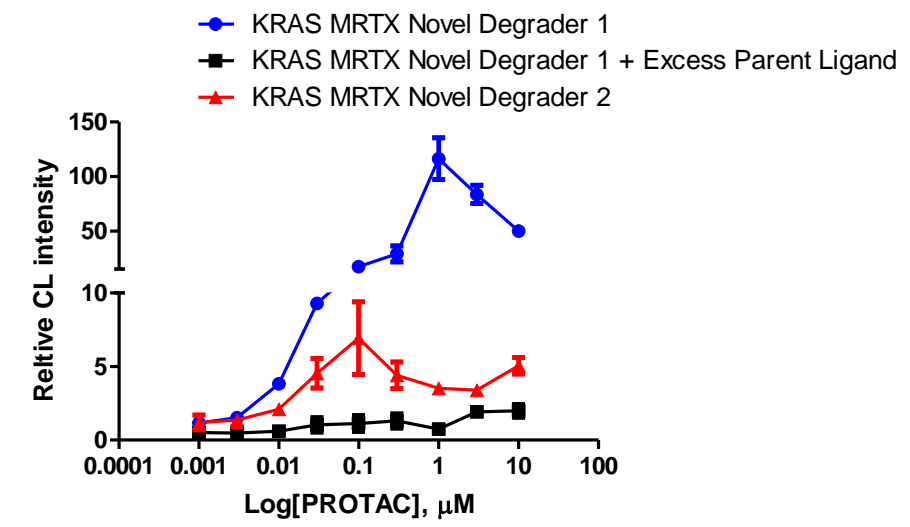
(A) *In vitro* ubiquitination assay with ARS 1620 KRAS G12C Novel E3 based degraders : dose response study to monitor PROTAC mediated ubiquitination of KRAS G12C. CL intensities plotted in response to 1/2 log dose response demonstrates PROTAC mediated ubiquitination. (B) Cell based assay to monitor degradation of KRAS G12C in H358 cells - PROTAC mediated degradation in a dose response study. (C) Representation of D<sub>max</sub> as dose dependent degradation.



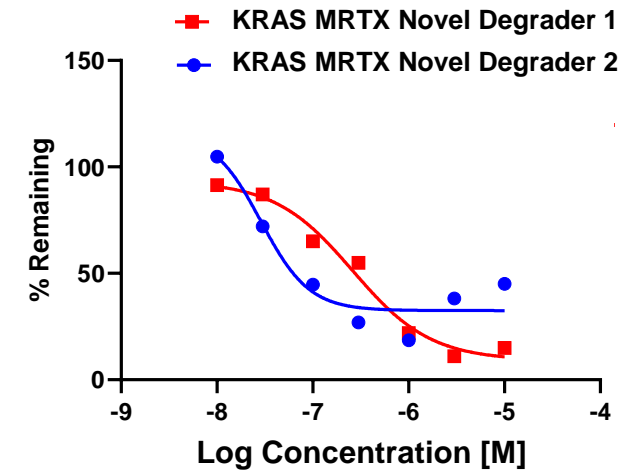
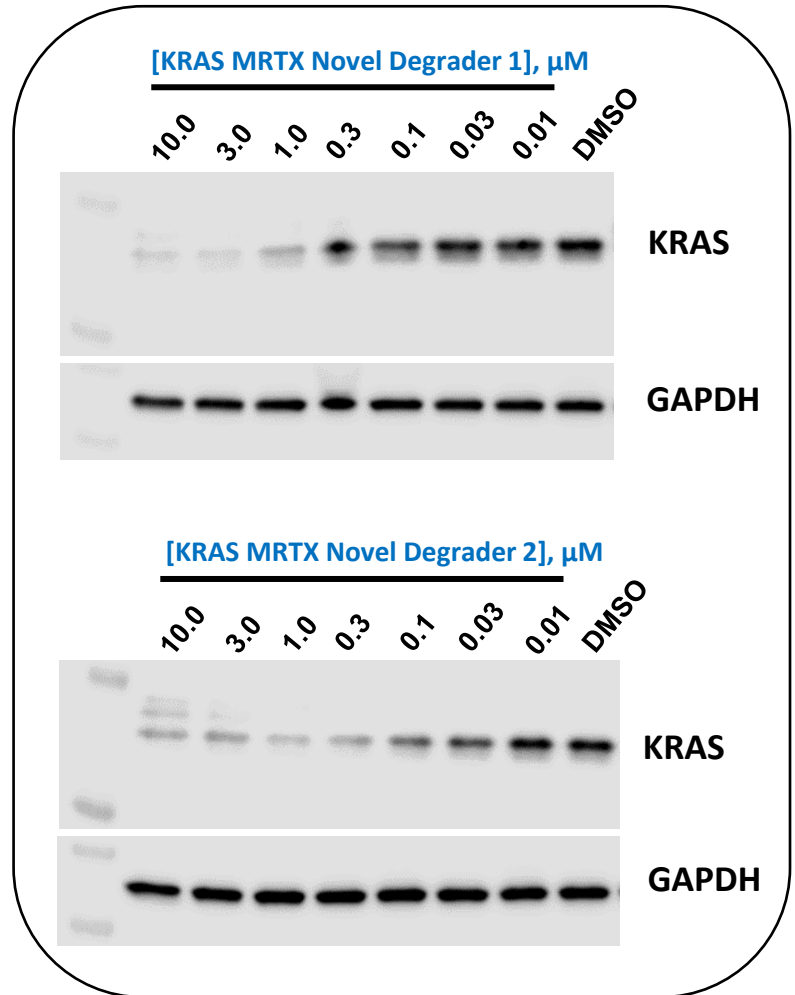
# K-RAS Cellular Ubiquitination and Degradation



# Novel E3 Ligand with MRTX-849 for KRAS Degradation



*In vitro* Ubiquitination

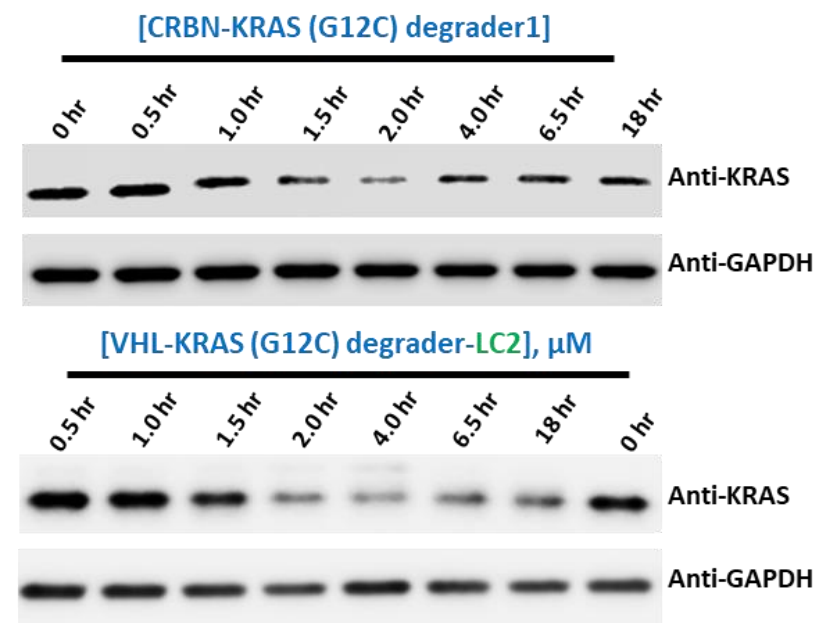
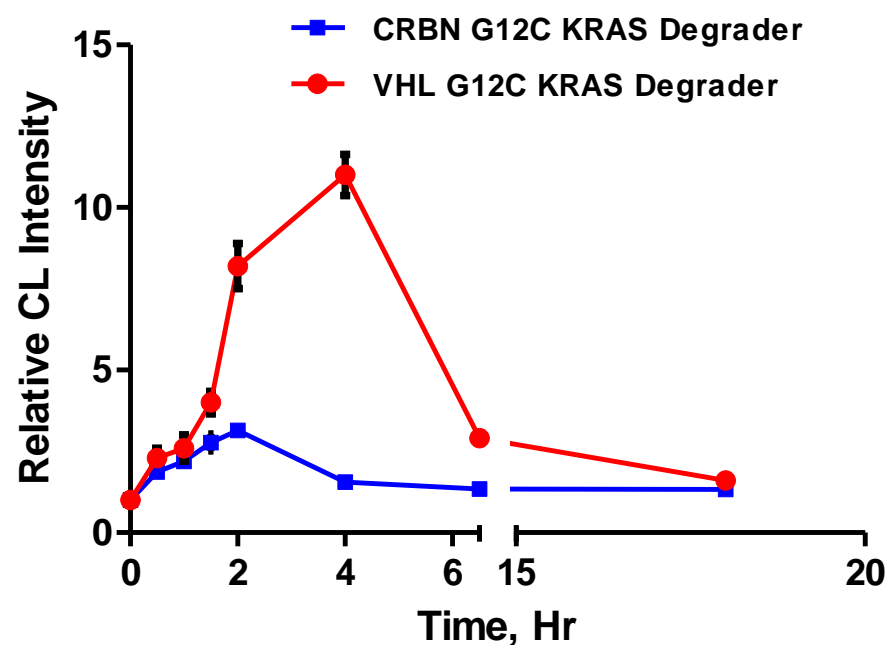


Cellular Degradation

# CRBN & VHL KRAS Degraders

Monitor PROTAC mediated Cellular Ubiquitylation and Degradation

## Time Course Study - HTS

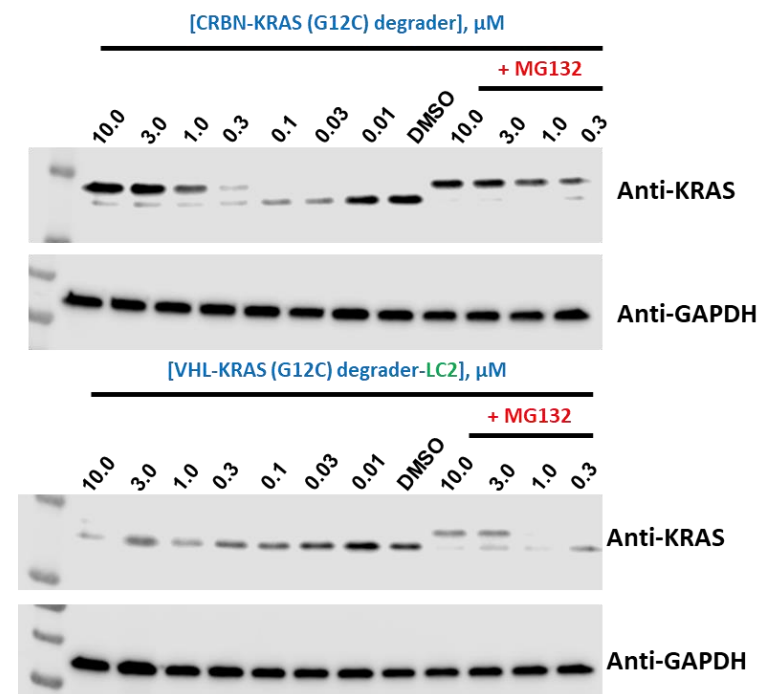
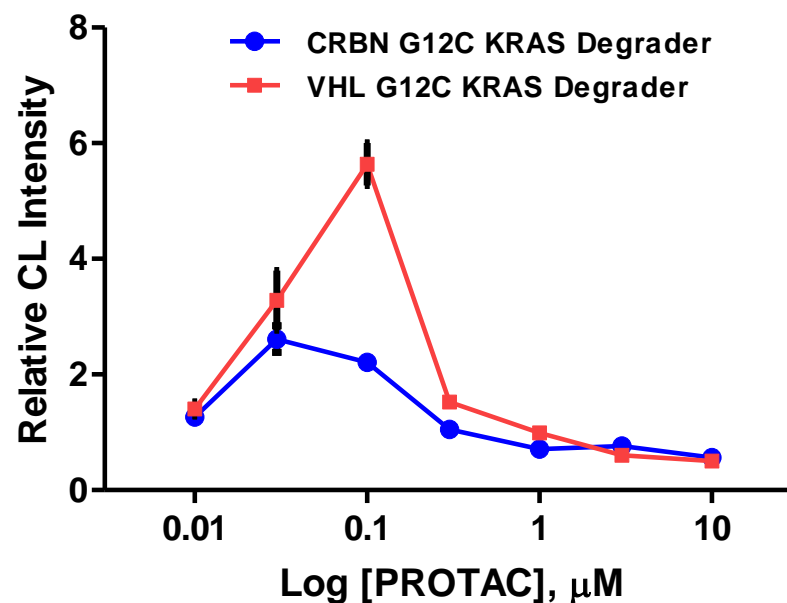


Monitoring Ubiquitination Kinetics – changes in PROTAC mediated ubiquitination profiles of endogenous KRAS and subsequent degradation in H358 cells. VHL and CRBN PROTACs designed with covalent ligands to engage KRAS G12C and successfully ubiquitinate and degrade within 2-4hrs of treatment.

# CRBN & VHL KRAS Degraders

Monitor PROTAC mediated Cellular Ubiquitylation and Degradation

## Dose Response Study - HTS

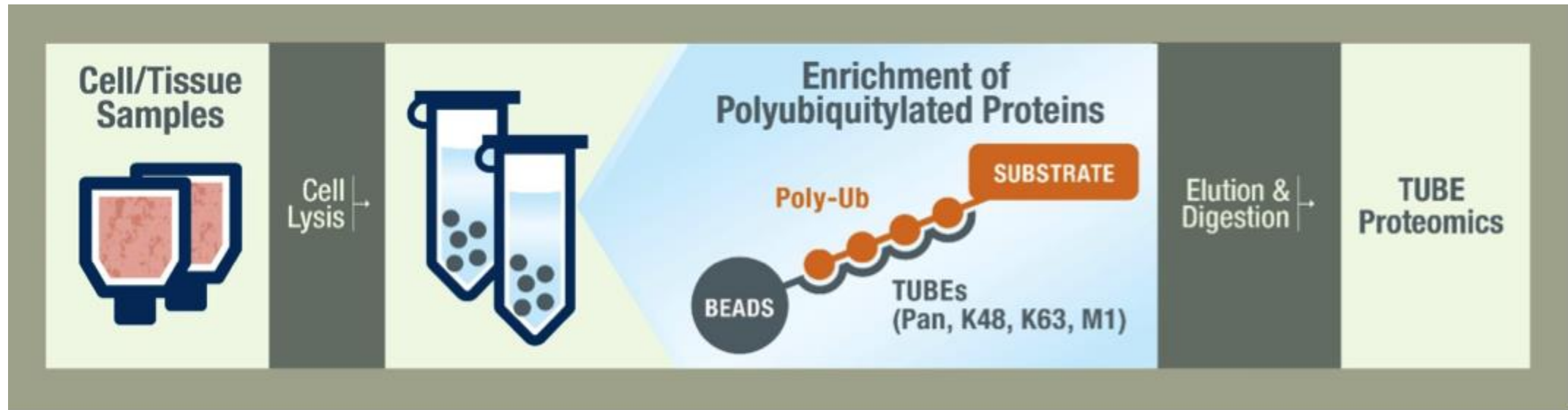


**Monitoring Ubiquitination – Dose Response** : changes in ubiquitination profiles of endogenous KRAS and subsequent degradation in H358 cells with changes in dose of both VHL and CRBN KRAS degraders. VHL and CRBN PROTACs designed with covalent ligands to engage KRAS G12C and successfully ubiquitinate and degrade with 3hrs of treatment between 30-100 nm.

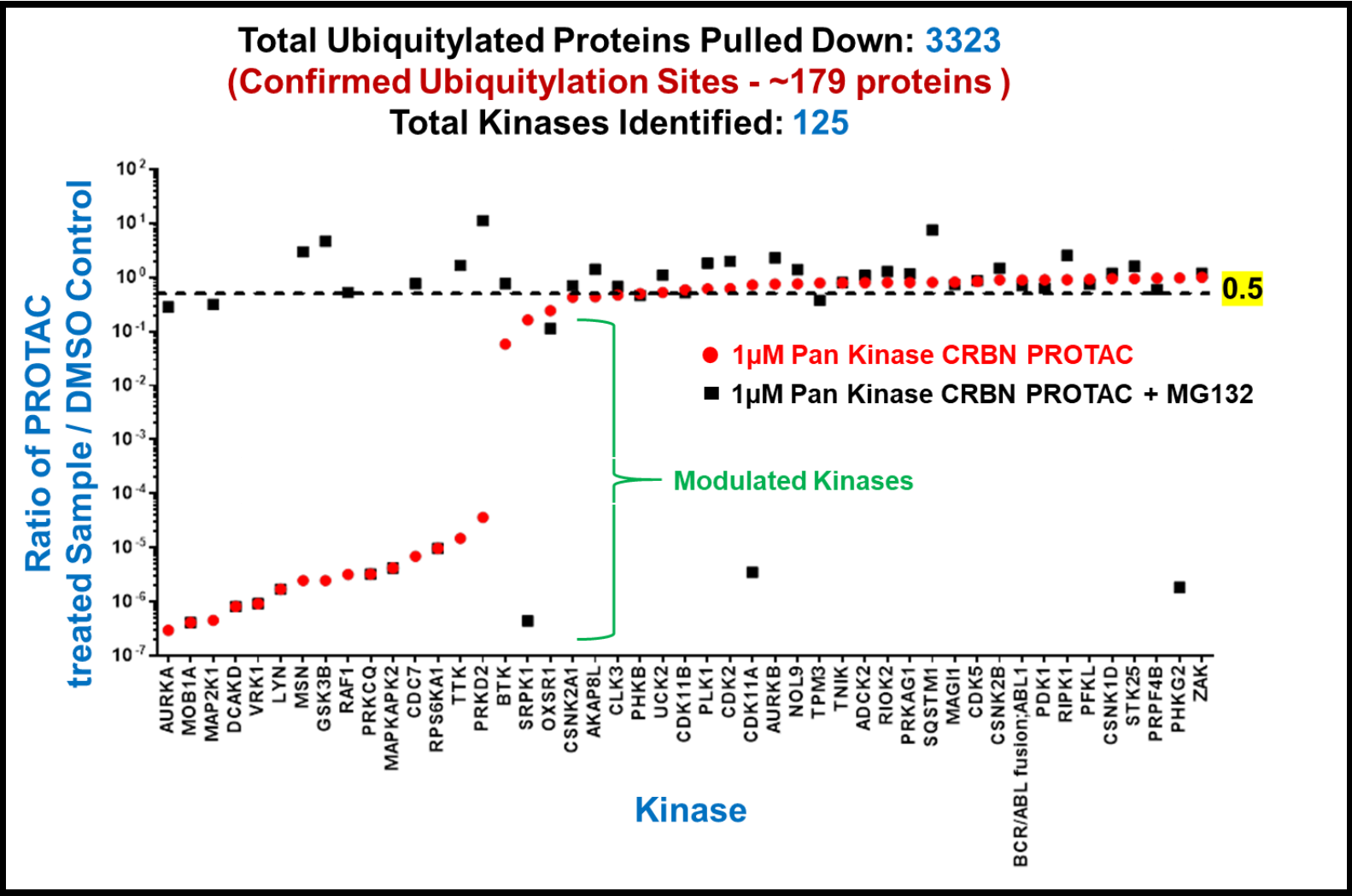


# TUBE Based Ubiquitin Mass spec Proteomics

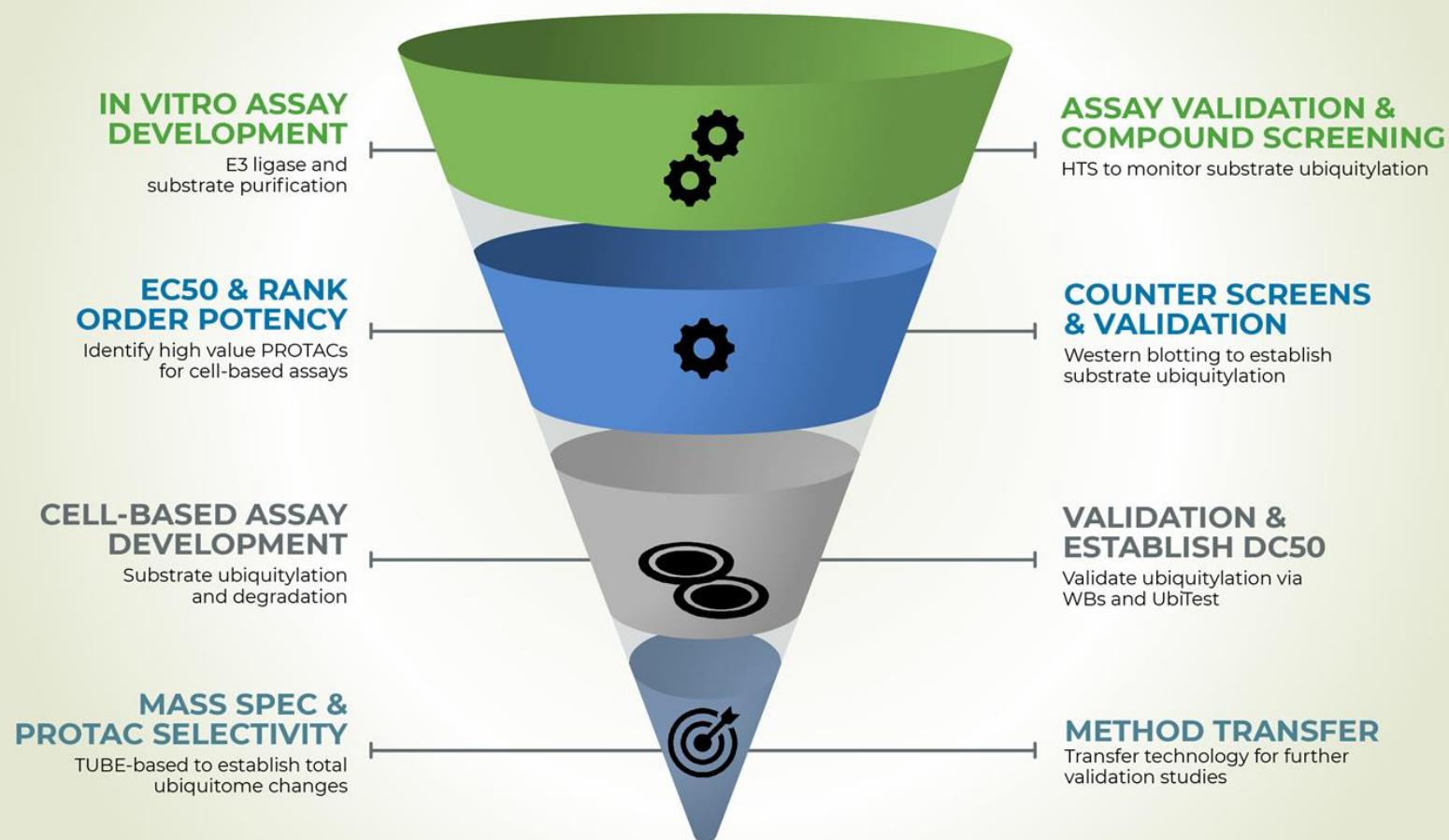
## To Validate PROTAC Mediated Ubiquitination and Degradation



# Mass Spectrometry Studies with Pan Kinase Degradator



# Pathway to PROTAC Drug Discovery



- ✓ **TUBEs based PROTAC® Assays** provides a link between ubiquitination and degradation
- ✓ **Robust SAR with true function of ubiquitination** for development of potent degraders



# Contact Us!

**We are your partner for PROTAC drug discovery**

## Contact Information

**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 354)**