

Deciphering the Ubiquitin Code with Poly-Ubiquitin Chain Selective Affinity Matrices

Rajesh K Singh¹, Cari Sagum², Jianji Chen², Apurva Chaturvedi¹, Peter Foote¹, Mark T Bedford²

¹Research & Development, LifeSensors Inc, Malvern, PA, ²Department of Epigenetics & Molecular Carcinogenesis, M.D. Anderson Cancer Center, University of Texas, Houston, TX

LifeSensors

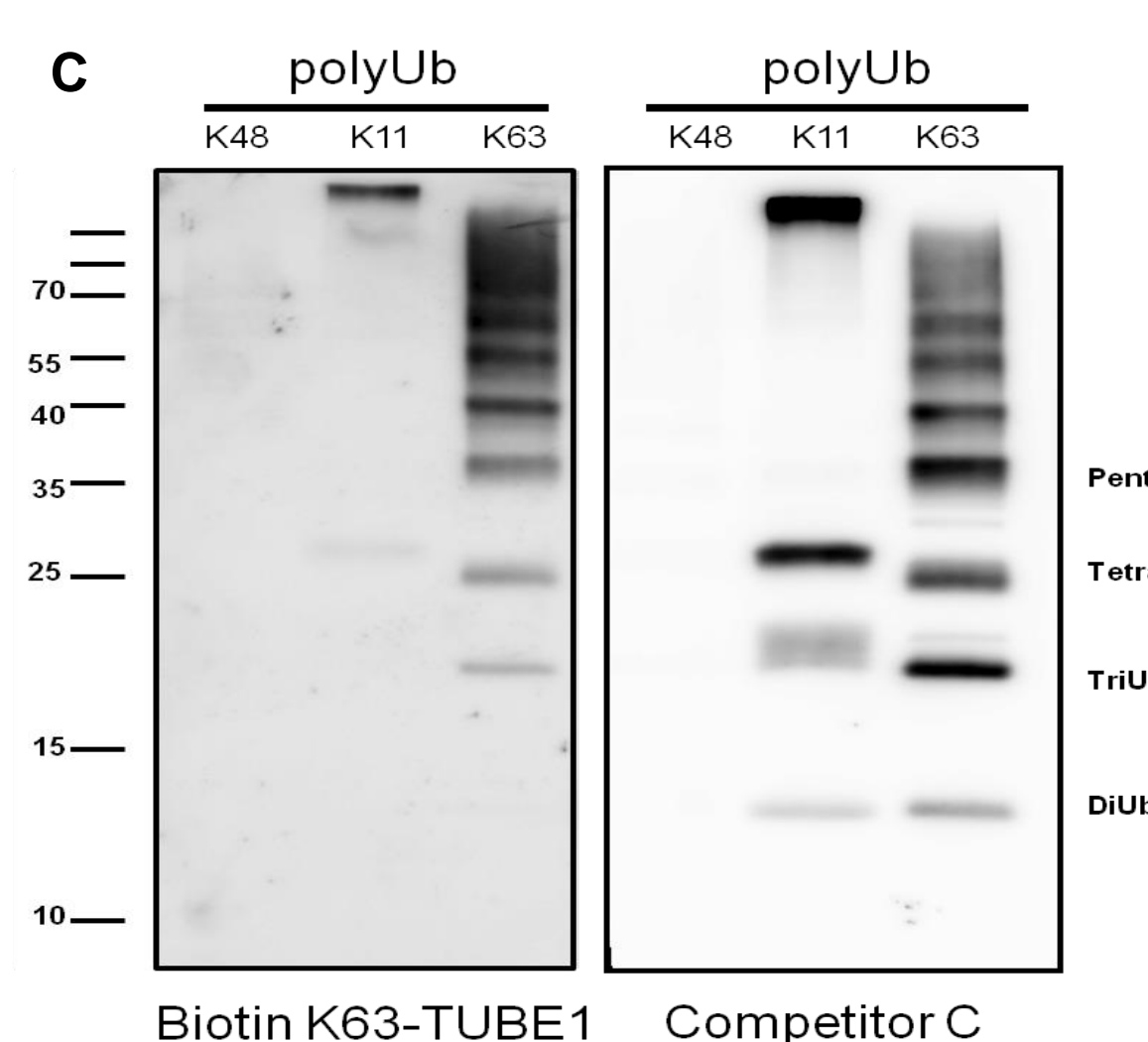
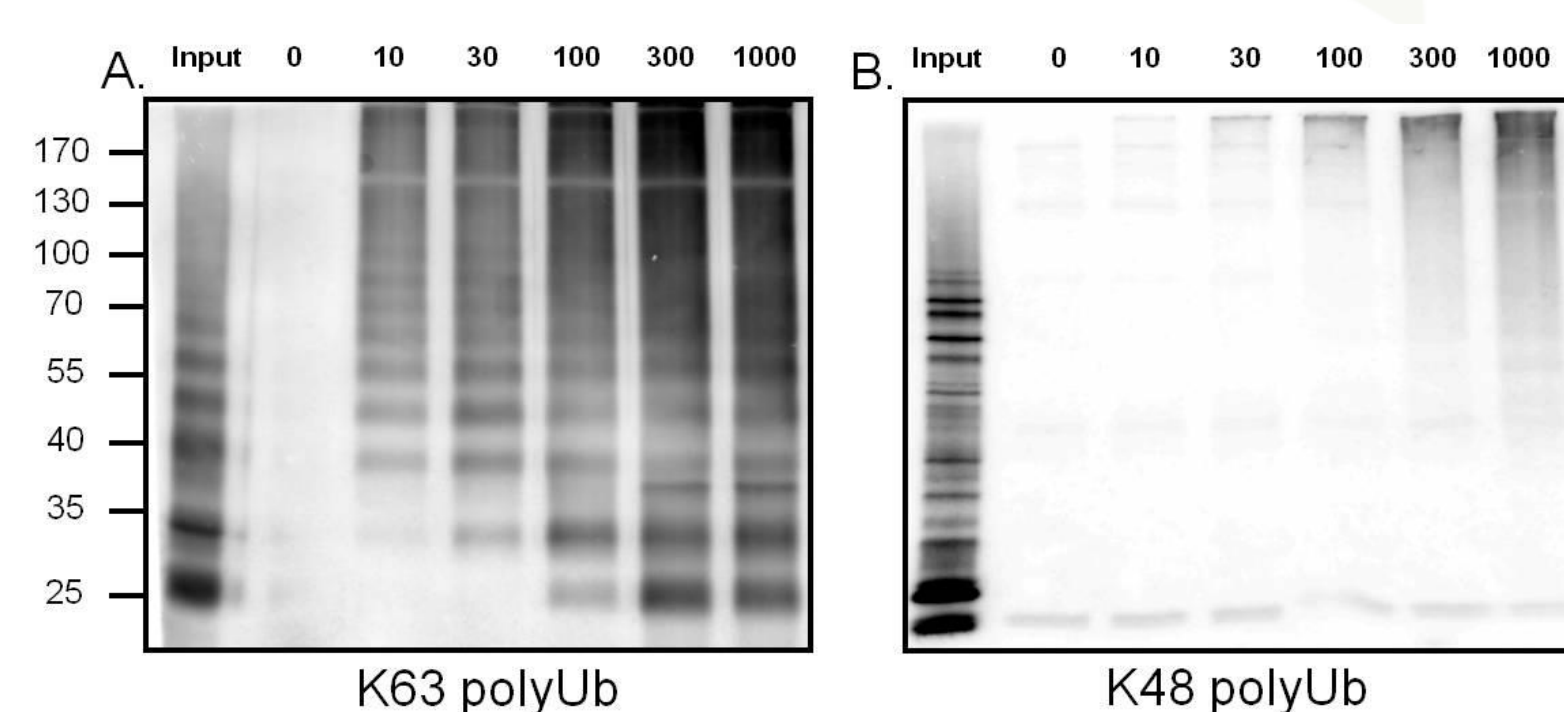
from genomics to proteomics

Abstract:

Purpose of study: The work presented here focuses on developing easy to use tools to distinguish between specific ubiquitin chain linkages, which are involved in a multitude of cellular functions.

Post-translational modification of proteins by ubiquitin (Ub) is a versatile process, highly dynamic, and involved in nearly all aspects of biological functions in eukaryotes. Dysregulation of ubiquitylation has been implicated in plethora of pathological conditions such as developmental abnormalities, neurodegenerative diseases, and cancers. Ub is attached via isopeptide bonds to lysine residues in the target protein or to another Ub to form poly-Ub chains. The reversibility, heterogeneity, and diversity of these modifications combined with the lack of suitable tools have made it difficult to properly isolate and characterize poly-ubiquitylated cellular proteins. However, in 2009, a novel technology called TUBEs (Tandem-repeated Ub-Binding Entities) was developed. TUBEs have revolutionized the Ub field by allowing poly-ubiquitylated proteins to be enriched/purified from cellular extracts. Development of TUBEs that selectively bind K48-, K63-, and M1-linked poly-ubiquitylated proteins has helped understand the role of modified proteins in cell physiology. However, the roles of other Ub-linkages remain obscured mainly due to the lack of tools that specifically recognize them. To overcome this, we recently developed a novel microarray platform that facilitates the identification of unique Ub binding domains (UBDs) for rare Ub-linkages. This array contains GST-tagged versions of 140 known and predicted UBDs from over 8 different Ub interacting families. Our initial screen with mono-Ub as well as K48-linked diUb revealed several unique binders. For example, Ub-binding motif (UBM) of REV1L bound uniquely to mono-Ub. On the other hand, Ub-associated (UBA)-like domain of NSFL1C bound preferably to K48-linked diUb. Moreover, UBA domains of RAD23A and RAD23B interacted uniquely with K48-linked diUb, which is concordant with the published data. We also identified several proteins that non-selectively bind mono-Ub and K48-linked diUb, such as the UBA domain of NBR1.

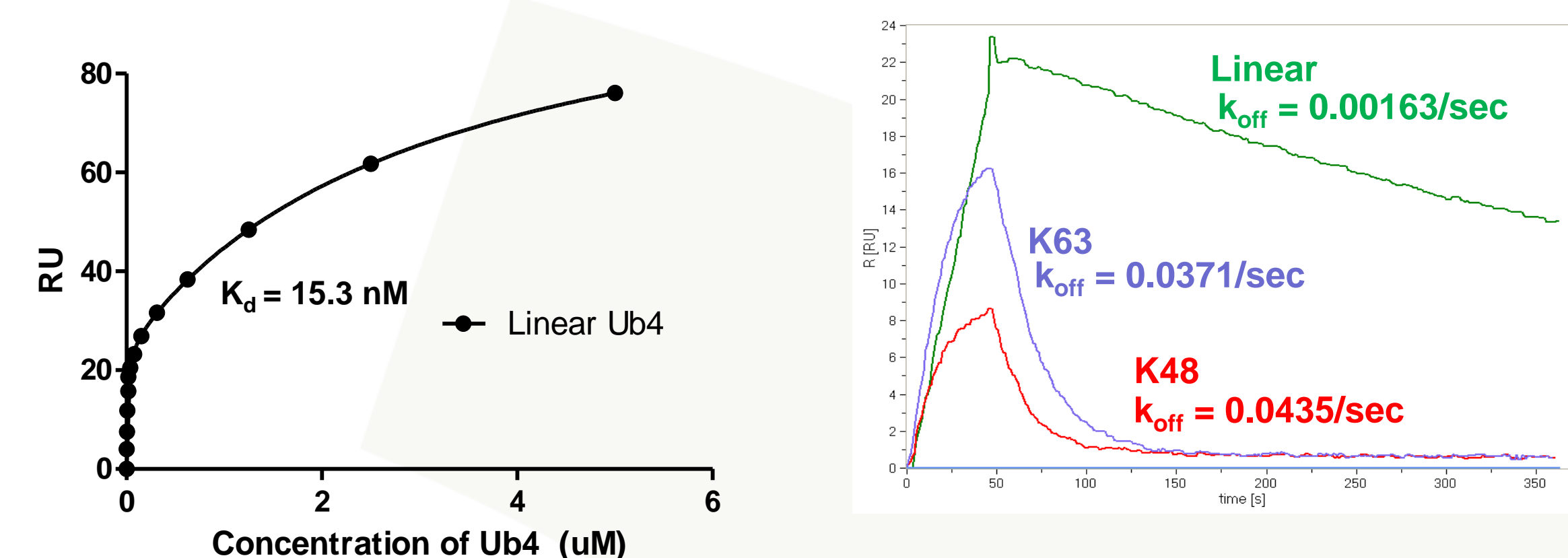
Enrichment of Cellular PolyUb with K63 TUBE



A-B. Pull-down Experiments using FLAG K63 TUBE. K63 linked PolyUb chains (10 µg) were generated by yMMS2/Ubc13 (Panel A) and K48 chains were generated by E2-25K (Panel B). The ubiquitin chains were pulled-down with indicated amounts of FLAG® K63-TUBE (nM) for 1.5 hours (4°C) and M2 FLAG Affinity Resin (10 µl, Sigma-Aldrich).

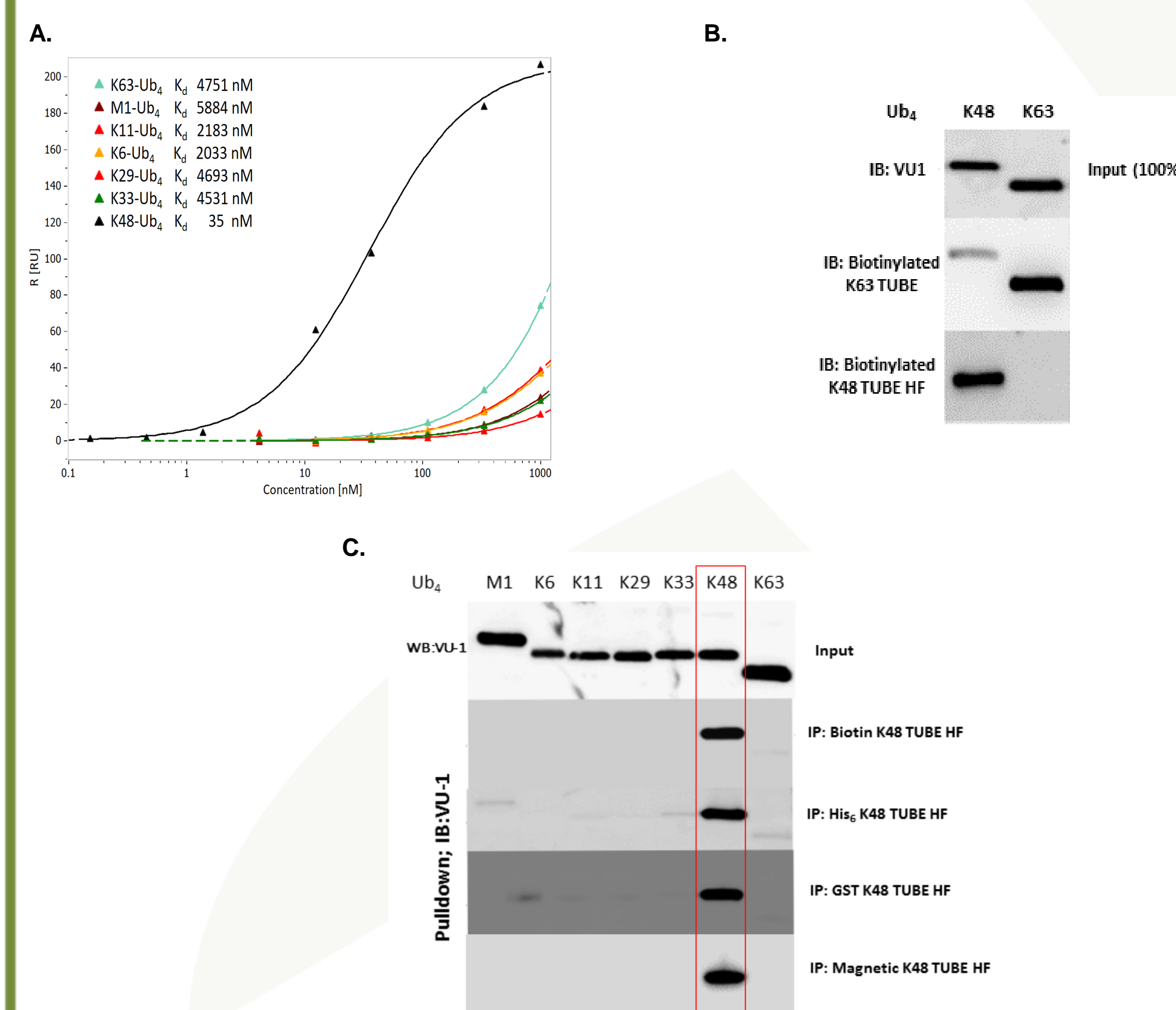
C. Comparison of K63 TUBE Specificity to Monoclonal Antibody. PolyUb chains generated with either E2-25K (K48), yMMS2/Ubc13 (K63), or UBE2S (K11) were subjected to duplicate loading, SDS-PAGE analysis and Western blot. The blot was sectioned and probed as described above with either Biotin K63 TUBE or a commercial monoclonal antibody specific for K63 polyUb chains to further investigate sensitivity and specificity for these two reagents.

Binding and Kinetics of M1 (Linear) TUBE



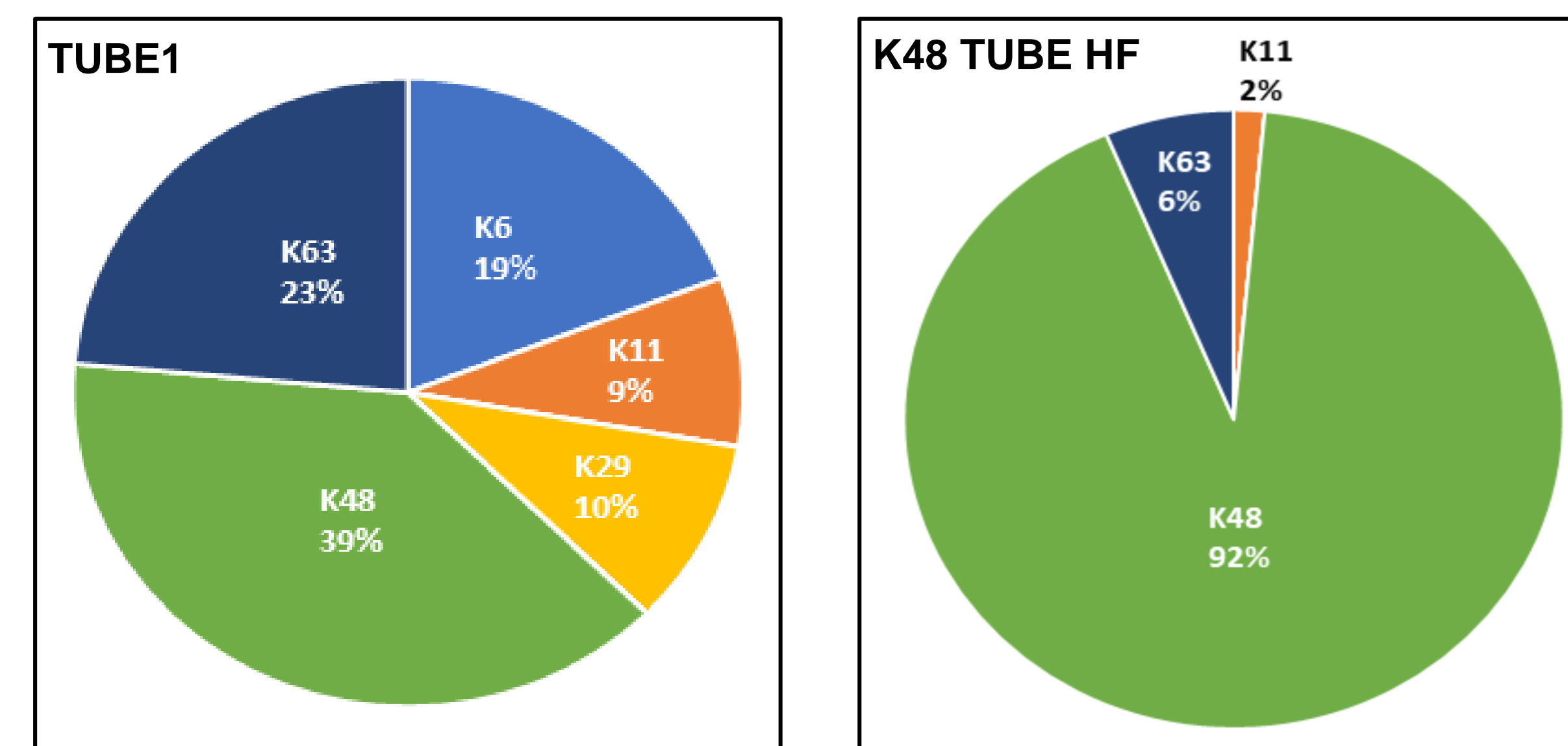
Determining Binding Affinity of M1 (Linear) TUBE. His₆-tagged TUBEs were immobilized on a His Capture chip. His₆-SUMO was loaded onto the control surface. Injections of increasing amounts of tetra-ubiquitin (Ub₄) were analyzed. The binding curve and K_d was generated in PRISM. (Right) Injections of 9.5 nM tetra-ubiquitin were performed on the same surface of Linear (M1) TUBE. Kinetic analysis was performed using Sierra Sensors SPR software.

Ubiquitin Binding Properties of the Newly Developed Highly Specific K48 TUBE HF



Determination of Binding Affinity and Specificity of K48 TUBE HF. (A) SPR experiments using increasing amounts of Ub₄ were performed using His₆-K48 TUBE and His₆-M1-Ub₂ on the control surface. The binding curves and K_ds were generated using Sierra Sensors analyzer software. (B) The detection of K48- and K63-Ub₄ by ligand blotting (far-Western) using biotinylated K63-TUBE and K48-TUBE HF. (C) 1 µg of the indicated Ub₄ were pull-down with 1.05 nmol of either biotinylated-, His₆-, GST- or Magnetic- K48 TUBE HF for 2 hours at 4°C. 20 % of the pulled-down samples were loaded onto SDS-PAGE followed by Western blot. The blots were probed with anti-Ub (VU-1) antibody.

Ubiquitin Proteomics using TUBEs



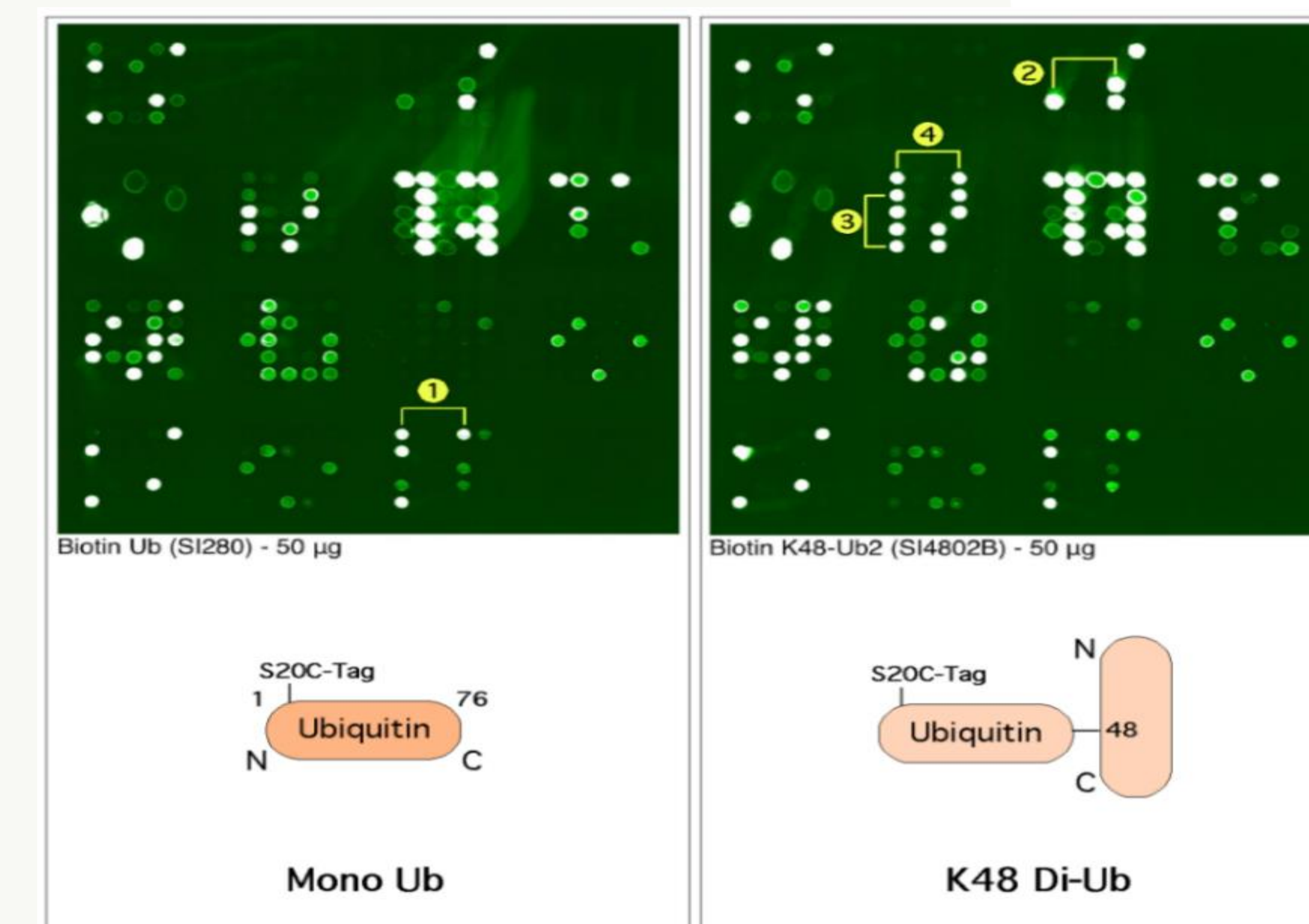
Ubiquitin proteomics utilizing TUBEs. Treated cells were enriched using magnetic TUBEs (TUBE1 and K48 TUBE HF). The eluted polyubiquitylated proteins were separated by SDS-PAGE and pixelated into 4 sections followed by digestion with trypsin. Individual ubiquitylated proteins were identified by the presence of at least 2 peptides in at least 2 of 3 replicates. Ub linkage types were identified by K-ε-GG peptides corresponding to covalent Ub-Ub modification at a particular lysine.

Ubiquitin Binding Domains Array

Ubiquitin-Binding Domains Array			
CUE A1) AMFR A2) ASCC2 A3) AUP1 A4) CUEDC1 A5) CUEDC2 A6) SMARCA4(1-2) A7) TAB2 A8) TAB3 A9) TOLLIP	JAB/MPN B1) MPN B2) MYSM1 B3) PRPF8 B4) PSM14 B5) PSM17 B6) STAMBPL1	UBA-like C1) C6orf106 C2) Cezanne/OTUD7B C3) DCUN1D1 C4) DCUN1D2 C5) FAM100A/UBALD1 C6) FAM100B/UBALD2 C7) NSFL1C C8) OTU07A C9) TTRAP/TDP2 C10) USP25	UBA D1) LATS1 D2) LATS2 D3) MARK1 D4) blank D5) MARK3 D6) MARK4 D7) NACA D8) NACA2 D9) NBR1 D10) NICE4/UBAP2L D11) NYREN18/NUB1(1-3) D12) p62/SQSTM1
UBA E1) RAD23A(1-2) E2) RAD23B(1-2) E3) RHBD3 E4) STS2/UBASH3A E5) TRKCC E6) UBAC1(1-2) E7) UBAC2 E8) UBE2K E9) UBL7 E10) UBL7 E11) UBL7 E12) UBL7	UBA F1) UBAP2 F2) UBQLN1 F3) UBQLN2 F4) UBQLN3 F5) UBQLN4 F6) UBXN1 F7) UBXN7 F8) USP13(1-2) F9) USP24 F10) USP5(1-2) F11) VPS13D F12) VPS13D F13) USP25	UBA G1) CBL G2) CBLB G3) EFTS/TFM G4) ETEA/FAF2 G5) FAF1 G6) TRD3 G7) STS-1/UBASH3B G8) TDRD3	UBA-Tudor H1) TDRD3
UIM H1) ANKIB1 H2) ANKRD13A(1-4) H3) ANKRD13D(1-4) H4) ATXN1(1-2) H5) ATXN3(1-3) H6) DNAJB2(1-2) H7) EPS15(1-2) H8) Epsin-1/EPN1(1-3) H9) Epsin-2/EPN2(1-2) H10) Epsin-3/EPN3(1-2) H11) PSM4(1-2) H12) RAB8/UMC2(1-2)	UIM I1) RNF166 I2) UBXN7 I3) USP25(1-2) I4) USP37(1-3) I5) ZFAND2B(1-2) I6) Hrs/HGS I7) blank I8) STAM2 I9) STAM1/STAM I10) STAM2 I11) UIM+UBX I12) UIM+UBX I13) UBXN7	GAT J1) GGA1 J2) GGA2 J3) GGA3 J4) Srcasm/TOM1L1 J5) TOM1 J6) TOM1L2 J7) GGA1 J8) GGA2 J9) GGA3 J10) Srcasm/TOM1L1 J11) TOM1 J12) TOM1L2	VHS K1) GGA2 K2) GGA3 K3) Hrs/HGS K4) Srcasm/TOM1L1 K5) STAM1/STAM K6) STAM2 K7) TOM1 K8) TOM1L2 K9) STAM1/STAM K10) STAM2
UIM+MIU L1) RNF168(UIM+MIU1) L2) RNF168(2) L3) RNF169(2) L4) PLAA L5) ETEA/FAF2 L6) FAF1 L7) NSFL1C L8) UBXN1 L9) UBXN7 L10) UEVLD	Znf M1) TAX1BP1(1-2) M2) USP13 M3) USP20 M4) USP3 M5) USP33 M6) USP39 M7) USP44 M8) USP49 M9) BRAP2 M10) USP5	UBM N1) hPol iota (1-2) N2) REV1L N3) hPol eta N4) hPol kappa (1-2) N5) PAD18 N6) SPARTAN Other N7) hRPN13/ADRM1 FL	gGST 1:1000

Microarray layout of ubiquitin readers. GST-tagged ubiquitin binding domains are spotted in duplicate. The arrangement is such that no two duplicates are immediately next to each other. Bottom right: anti-GST was used to ensure equal spotting of each domain. Note that different grids have different number of proteins as described in the family of proteins in the table. These domains are selected after thorough searches and analysis of families of proteins that contain UBDs.

Screening of UBD Microarray with Mono-Ub and K48-linked diUb



Selective binding of mono-ubiquitin and K48-linked di-ubiquitin to UBD Protein Microarray. Monoubiquitin (left) and K48-linked di-ubiquitin (right) were incubated with the microarray of UBDs and tested for proof-of-concept experiments. The microarray grid is the same as shown in the middle panel. Numbered spots (yellow dots) indicate ubiquitin binding domains that show preliminary specificity (each domain is spotted in duplicate and shown by connected line).

Conclusions:

A collection of linkage specific TUBEs will dramatically improve current methods for studying ubiquitin chain linkages and their regulation/prevalence in disease states. The array allowed us to identify unique binding partners for mono-Ub and K48-linked diUb, and we strongly believe that our future screens with other Ub-linkages will identify several selective binding partners for rare Ub-linkages. Domains that display selective binding specificity will be utilized to construct linkage-specific TUBEs. The expansion of the TUBE toolkit to include recombinant proteins that can discriminate between each of the Ub linkages would dramatically accelerate the pace of discovery in this important area of biology.