

# Automated online protein-ligand binding and its detection using native mass spectrometry

Weijing Liu<sup>1</sup>, Wilson Phung<sup>2</sup>, Wendy Sandoval<sup>2</sup>, Rosa Viner<sup>1</sup>

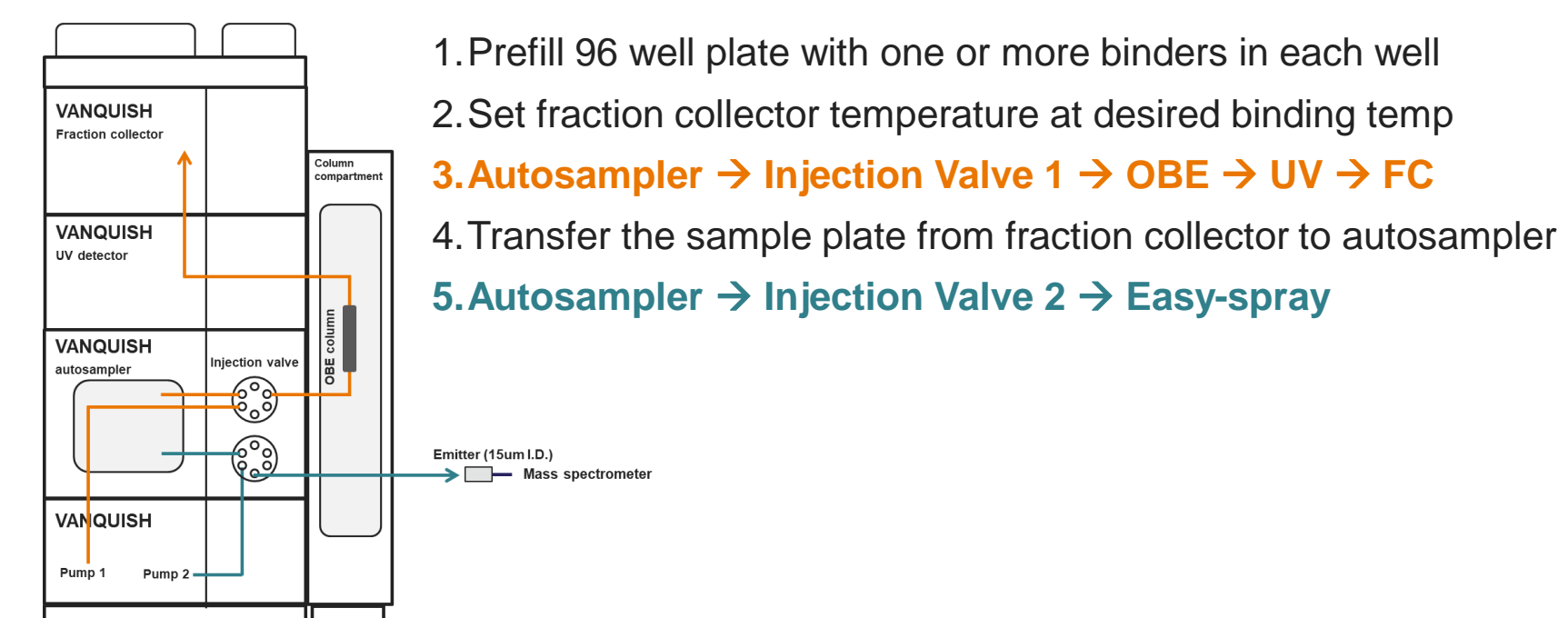
<sup>1</sup>Thermo Fisher Scientific, San Jose, CA; <sup>2</sup>Genentech, South San Francisco, CA

## Abstract

Many proteins involved in diseases are considered "undruggable" using traditional small-molecule inhibitors due to their lack of well-defined binding pockets. Molecular glue (MG), by strengthening weak intrinsic interaction between target and E3 ligase, enables the targeted protein degradation via the ubiquitin-proteasome system. Native mass spectrometry (nMS) preserving the noncovalent interactions allows the identification of E3-MG-target ternary complex. However, offline manual sample preparation and direct infusion limits the throughput. A newly designed integrated LC system coupled to native MS streamlines the rapid online buffer exchange, parallel compound binding, and direct nMS analysis of complexes. The novel workflow has been applied for MG screening to assess their efficacy by directly detecting the ternary complexes.

## Materials and methods

### Methods



### Instruments

Thermo Scientific™ Vanquish™ LC: UV detector, fraction collector, dual injection autosampler, and dual Flex pumps

Thermo Scientific™  
Q-Exactive™ UHMR



Thermo Scientific™  
Orbitrap™ Ascend™  
Structural Biology edition



- m/z 350-80,000
- Quad up to m/z 25,000
- Native MS
- Native top-down
- Direct Mass Technology

- m/z 50-16,000
- Quad up to m/z 8,000
- Native MS
- Native top-down
- PTCR/HCD/CID/ETD/UVPD

### Data Analysis

Data analysis was performed with Thermo Fisher™ BioPharma Finder™ 5.0

## Results

### 1. Carbonic anhydrase - Ligand binding

➤ We initially utilized Carbonic Anhydrase (P) as the target to evaluate ligand binding efficacy of the proposed workflow. Initial tests employed 4 ligands with varying  $K_d$  values, ranging from <100nM to >10uM.

Table 1. Carbonic Anhydrase ligands

Compound	MW	Ligand#	$K_d/\mu\text{M}$
Sulfanilamide	172.20	L1	13.2
Benzenesulfonamide	157.19	L2	1.44
1,3-Benzenedisulfonamide	236.27	L3	~1.1
4-Sulfamoylbenzoic acid	201.20	L4	0.27

Figure 1. Comparison of pre & post-column binding

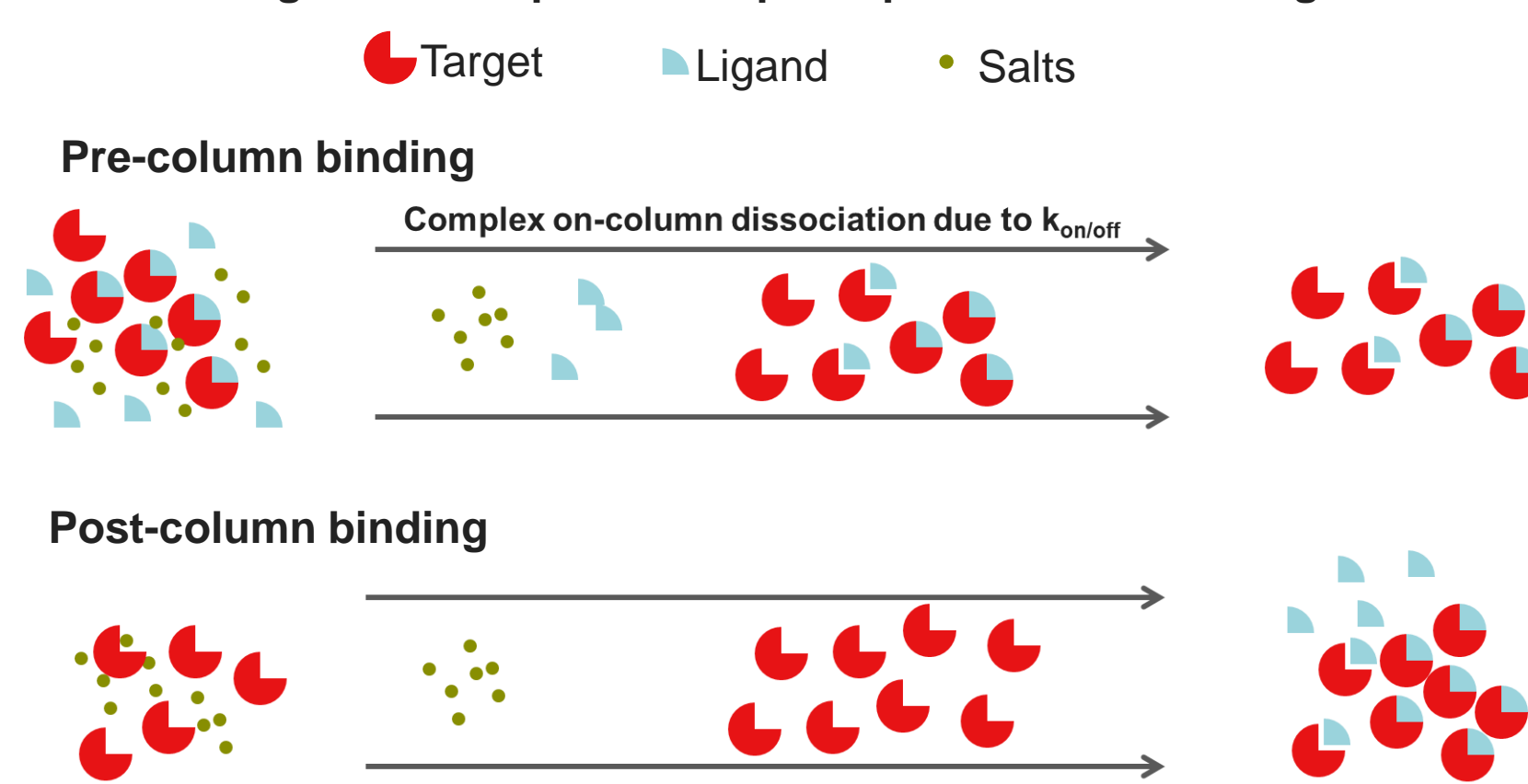
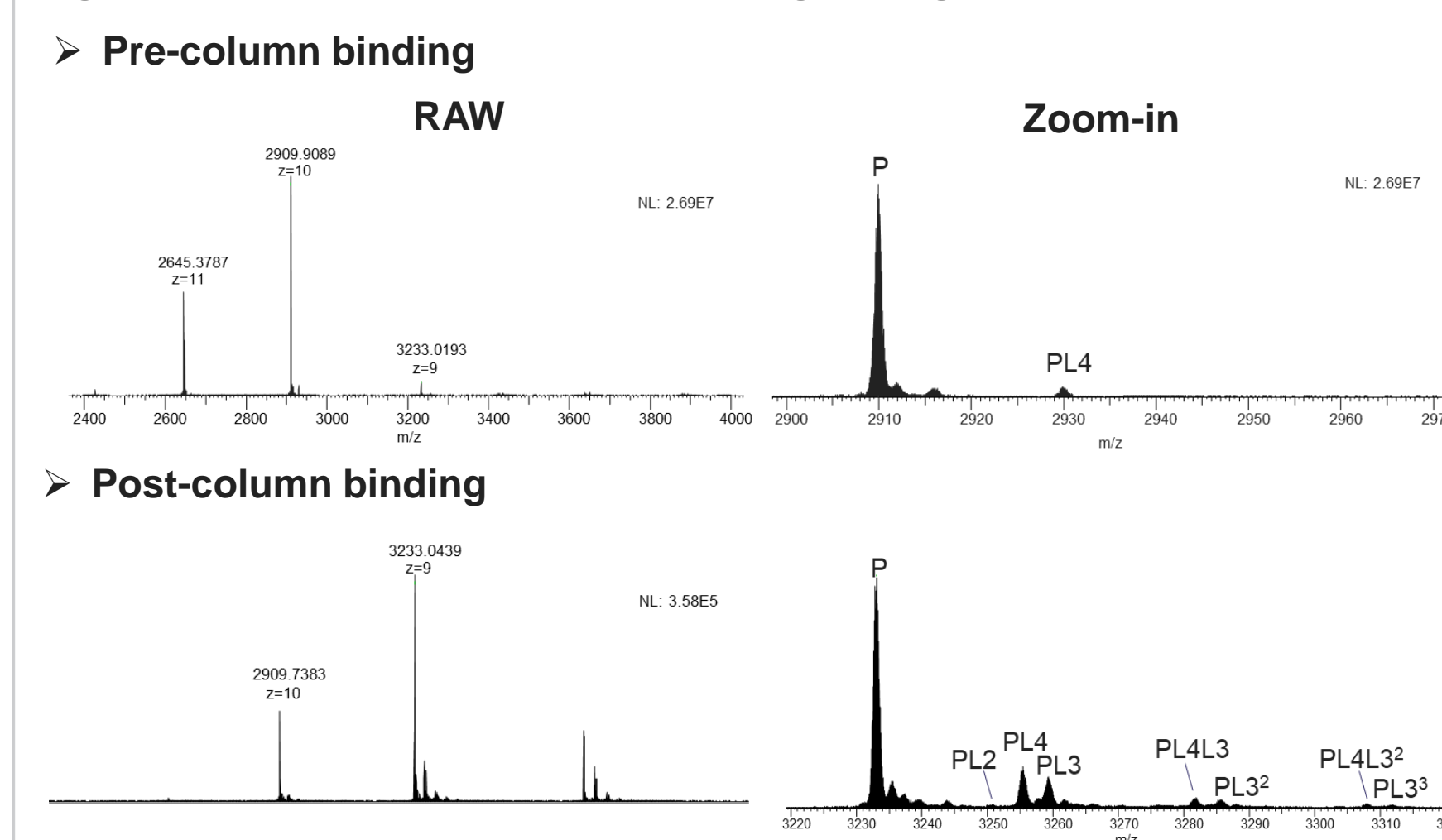


Figure 2. MS spectra of Carbonic binding to 4-ligand mix pre & post-column



- ✓ Pre-column binding shows on-column dissociation
- ✓ Apparent  $K_d$ : L1>L2>L3>L4, aligning with published  $K_d$
- ✓ Strongest ligand L4 shows most abundant primary binding

➤ Fraction collection of online desalted Carbonic Anhydrase prior to ligand binding

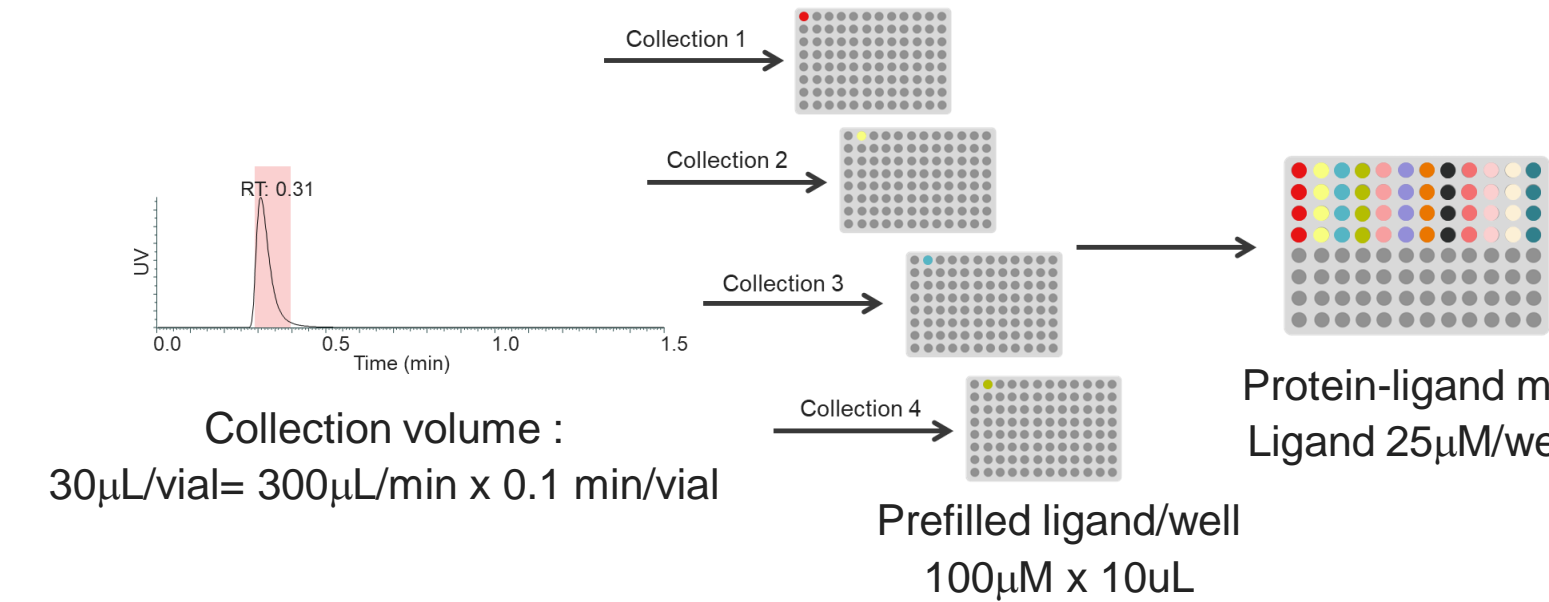
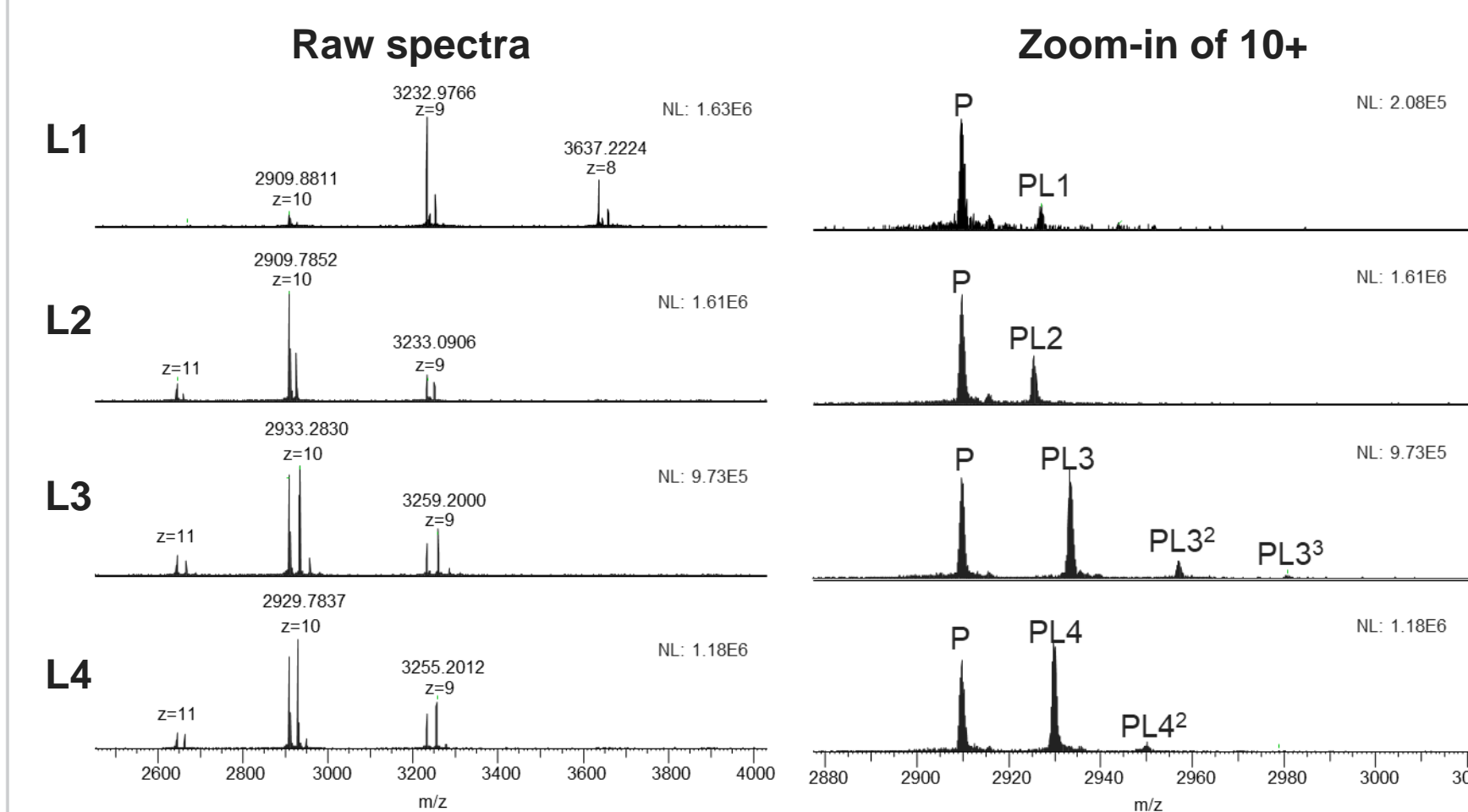


Figure 3. Carbonic Anhydrase binding to 4 ligands respectively



- ✓ PL fractional abundance%: L1<L2<L3<L4
- ✓ Apparent  $K_d$ : L1>L2>L3>L4
- ✓ Rank of  $K_d$  aligns with published results

### 2. Molecular glues screening

➤ Molecular glues stabilize interactions between E3 ligases and therapeutically relevant substrates to initiate substrate degradation via the ubiquitin-proteasome system.

Figure 4. Molecular glue (MG) strategy premise

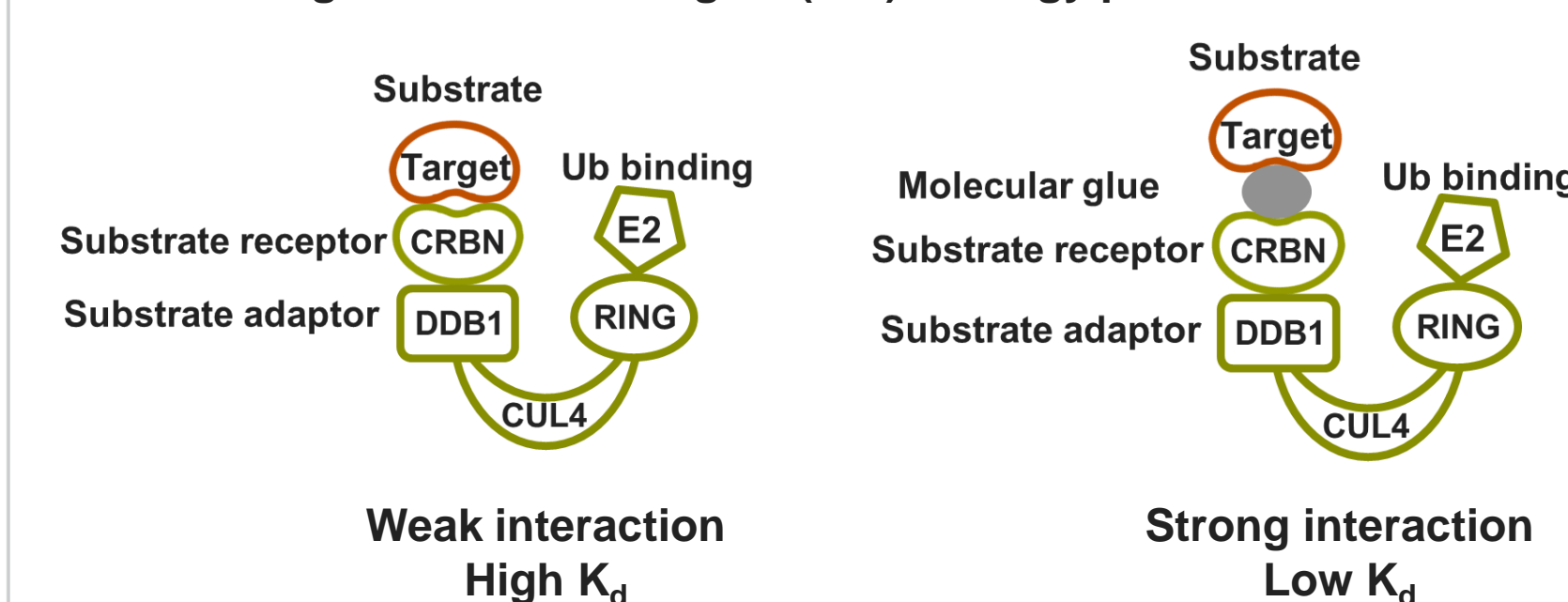


Table 2. Molecular glues of CRBN-DDB1 and target

Glue	MW	Ternary complex $K_d$ ( $\mu\text{M}$ )	Comments
A1	483.46	0.0941	Control, strong $K_d$
B1	584.02	0.003	Top Glue with low $K_d$
C1	577.00	0.035	Top Glue with low $K_d$
D1	456.51	0.913	High $K_d$
E1 (-Ctrl)	441.86	>1	Negative control binds CRBN

Figure 4. MS spectra of CRBN-DDB1 and target binding to 5 MGs

CRBN-DDB1:Target:MG = 2.3 $\mu\text{M}$ :2.3 $\mu\text{M}$ :23 $\mu\text{M}$

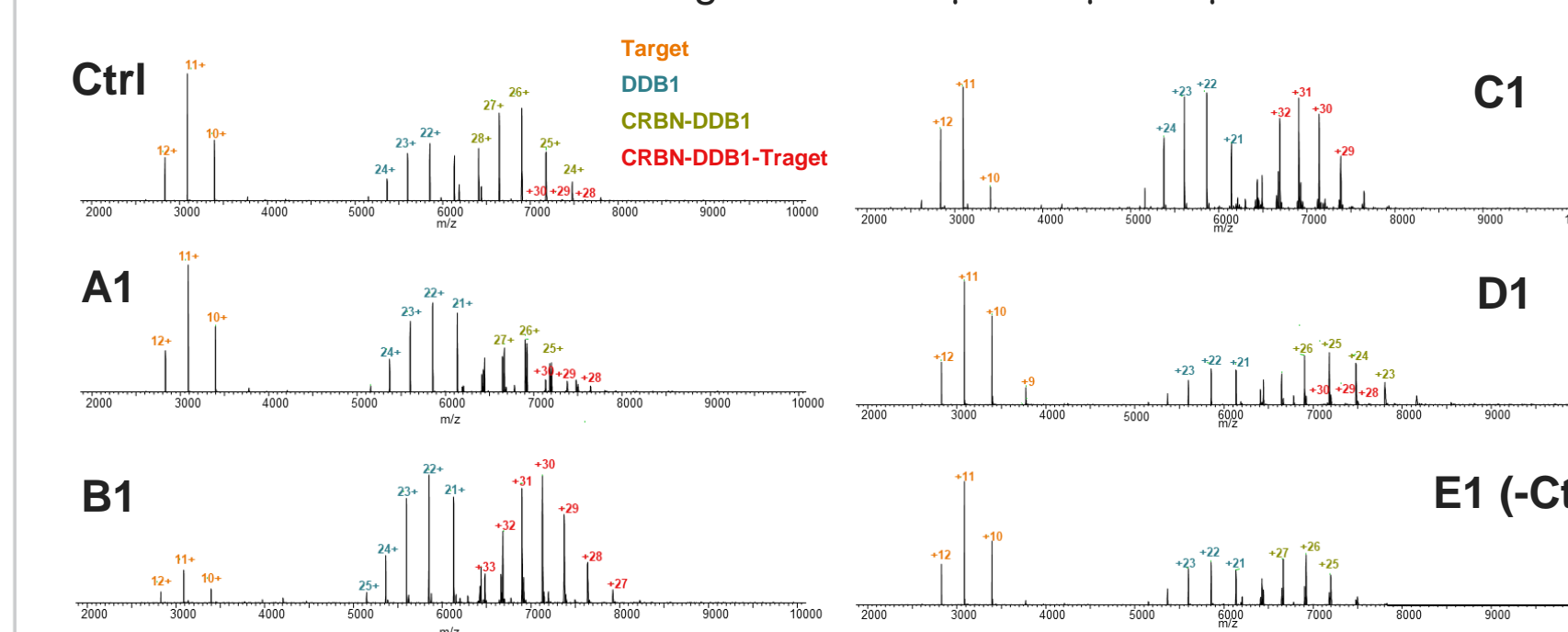
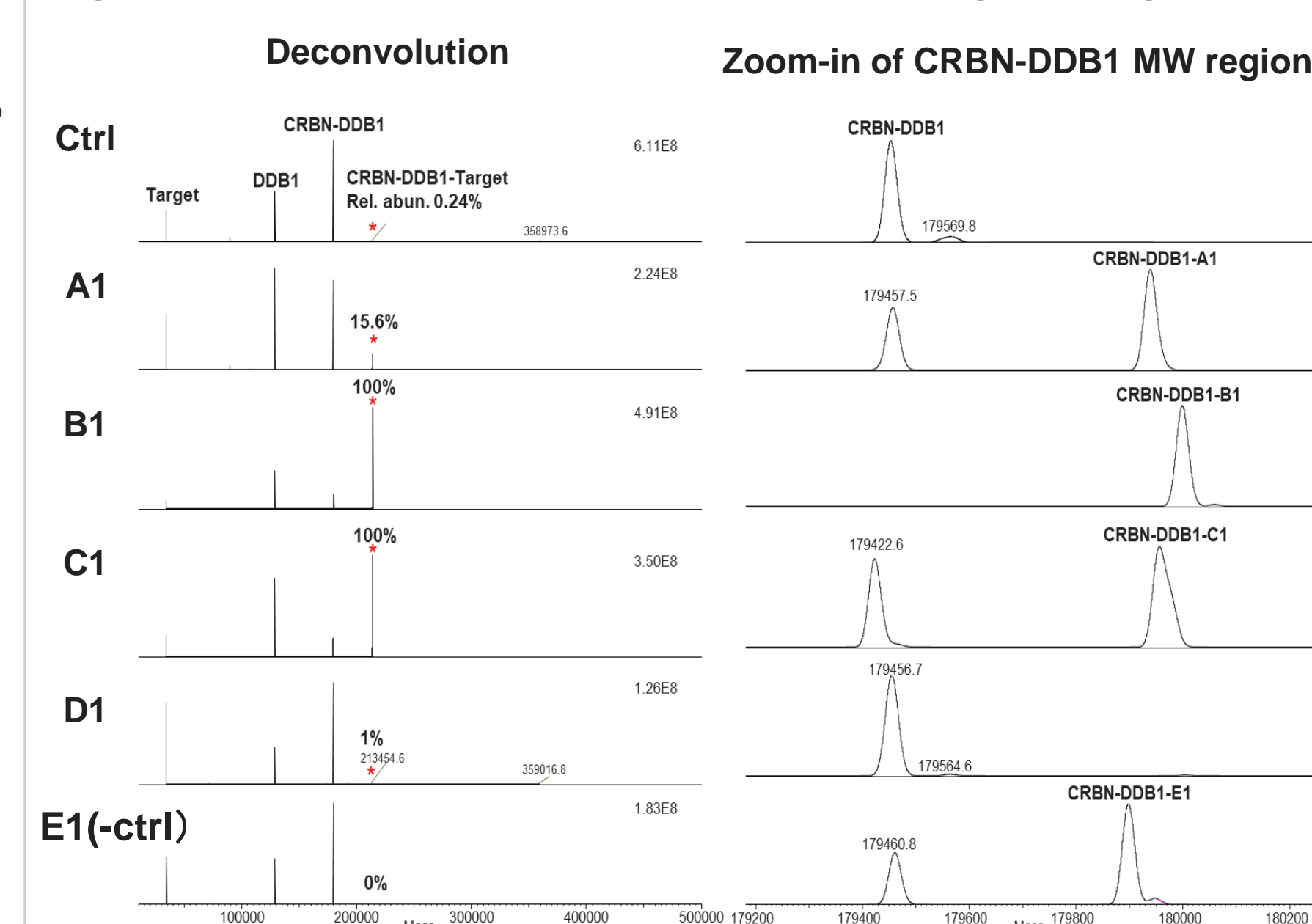


Figure 5. Deconvoluted spectra of CRBN-DDB1 and target binding to 5 MGs



- ✓ Apparent  $K_d$ : C1<B1<A1<D1<Ctrl<E1 (-Ctrl) aligns with results from biophysical assays
- ✓ Strong MGs show more binding on CRBN-DDB1

Table 4. Stability of CRBN-DDB1-Target complex

Glue	$K_d$	Frac. Abun. @ RT 10min	Frac. Abun. @4°C overnight
B1	Low	100%	97%
C1	Low	100%	22%
D1	High	1%	1.5%
E1 (-Ctrl)	Negative control	0%	0%

- ✓ Relative abundance of CRBN-DDB1 and CRBN-DDB1-Target dropped overnight
- ✓ Samples are unstable in 200mM AmAc, reflecting the importance of rapid screening after buffer exchange

## Conclusions

1. On-column dilution and dissociation prevent the detection of weak binders.
2. Post-column binding facilitates the observation of weak binding event  $K_d$ >10uM.
3. Apparent  $K_d$  aligns well with published results or measurements from other assays.
4. Fraction collection coupled to direct infusion enables high-throughput ligand screening.
5. This workflow has been applied for MG screening to assess their efficacy by directly detecting the ternary complexes.

## Acknowledgements

- Jennie Lill at Genentech
- Dennis Koehler at Thermo Fisher Scientific
- Mark Tracy at Thermo Fisher Scientific

## Trademarks/licensing

© 2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others. PO154-2024-EN