

# **Selection of Peptide Targets for Species-independent Quantitation of** Fish Allergens

### Introduction

Detection and quantitation of fish residues in foods for allergen control poses considerable challenges. The diversity of fish used in the food supply necessitates the use of analytical targets that are shared, and are quantitatively similar, in all potential contaminant species. Identifying such targets is challenging. Currently available methods, reliant on ELISA, targeting Parvalbumin, display unacceptable (often 200-fold) variation in performance across fish species rendering quantitation near-impossible. We designed a workflow employing untargeted high-resolution MS with label-free quantitation to identify peptide targets that are present in a broad range of consumed fish, that are similarly abundant, and that are not present in non-fish.

### Goals

- Determine the suitability of major allergen quantitation for broad-spectrum fish detection.
- Identify peptide targets that are present in a broad range of fish, are similarly abundant, and are not present in non-fish. Use these to develop a method to quantify the major allergens of fish (Parvalbumin, Enolase and Aldolase) across 9 fish species.

# Methods

Nine fish species, PCR speciated (cod, pollock, herring, salmon, tuna, skate, tilapia, grouper and halibut) were selected to represent the broad taxonomies of heavily consumed fish. Protein was extracted from 3 replicate samples (0.5 g) of the fillet using a reducing urea/thiourea extraction method (10 mL of 6 M Urea, 2 M Thiourea, 50 mM Tris-HCl, pH 8.6, 50 mM DTT).

Untargeted MS (3 replicate analyses of 3 replicate samples) was performed using a Thermo Scientific<sup>™</sup> Q Exactive<sup>™</sup> HF mass spectrometer coupled to a Thermo Scientific<sup>™</sup> UltiMate<sup>™</sup> 3000 nanoRS liquid chromatography (UPLC) system equipped with a Acquity UPLC M-Class Peptide CSH C18 Column 130A, 1.7µm, 75µm x 250mm (Waters Corp). Peptide, protein ID and label-free quantitation was performed using PeaksQ v8.5 with a combined protein database of all sequences available (UniProt) for analyzed fish species. Subsequent analysis of peptide quantitation data was performed using Microsoft Excel. BLAST P searches were performed against the whole NCBI database. Figure 1 shows an overview of the methodology and peptide selection strategy.



# Justin T. Marsh<sup>1</sup>, Charles Yang<sup>2</sup> and Philip Johnson<sup>1</sup>

<sup>1</sup>Department of Food Science and Technology, University of Nebraska-Lincoln, NE; <sup>2</sup>Thermo Fisher Scientific, San Jose, California \*Corresponding author: Philip.Johnson@unl.edu

# Label-free Quantitation of Allergens from Multiple Fish Species

### Parvalbumin

- 51 peptides from 14 Parvalbumin isoforms were detected.
- Label-free quantitation achieved by a single unique peptide to each Parvalbumin isoform family (See alignment in Figure 2a).
- Only 4 of the isoform families are present in all fish studied. Parvalbumin shows a variable abundance (6-fold difference between Tilapia and Tuna)
- across the fish species (See Figure 2b).





# Identification of "Universal" Fish Peptide Targets

- CoV all fish.

# allergens.

- peptides detected.

Our DDA-dependent strategy of peptide target selection appears to be valid work-flow, and could well be applied to other cross-species studies. However, one limitation to this strategy has been the limited proteome data from some fish. Thus, our quantitation may be biased to the fish species with comprehensive databases (e.g. Salmon and Tilapia). Theses considerations do not impact out target selection strategy though. Further work will include screening the peptide candidates against each fish, using a PRM methodology and finally in relevant fish spiked food matrices.

This research is part of a collaboration between FARRP/UNL and Thermo Fisher Scientific<sup>™</sup>. This work is supported by Improving Food Safety 1019035 from the USDA National Institute of Food and Agriculture.





• 200 peptides met our limitation of <150

 After screening these peptides with BLAST P against NCBI, only 25 % of the peptides were shown to be fish-specific. • The majority of the candidate peptides are, as expected, derived from muscle associated proteins (See Figure 5). Interestingly, only 1 peptide from the major allergens and 1 peptide from the minor allergens fit our filtration criteria.



## Discussion

• Quantitation of the major allergens of fish reveals high variation of abundance in the

• A complex Parvalbumin isoform distribution exists in fish.

• Concluding from this work, parvalbumin is likely not an ideal analytical target for fish residue quantitation. This is consistent with the known analytical performance of ELISA methods based on parvalbumin detection.

• Multiple "universal" fish peptide target candidate peptides, derived from non-

allergens, have been identified. This pool of candidate peptides comprises 0.5 % of the

• As might be expected, the need to quantify cartilaginous fish (here, skate) complicates the development of a 'universal' fish detection method.

## **Conclusions and Future Work**

### Acknowledgements