

# Measuring of Non-reducing Terminal Glycosaminoglycan Fragments increases specificity and differentiates Mucopolysaccharidosis Type I (MPS I) from Mucopolysaccharidosis Type II (MPS II)



Taraka Donti<sup>1</sup>, Sara E. Smith<sup>1</sup>, Sasha Chernenkoff<sup>1</sup>, Madeline Ellgass<sup>1</sup>, Madhuri Hedge<sup>1</sup>  
1. Revvity Omics, Pittsburgh, PA 15275

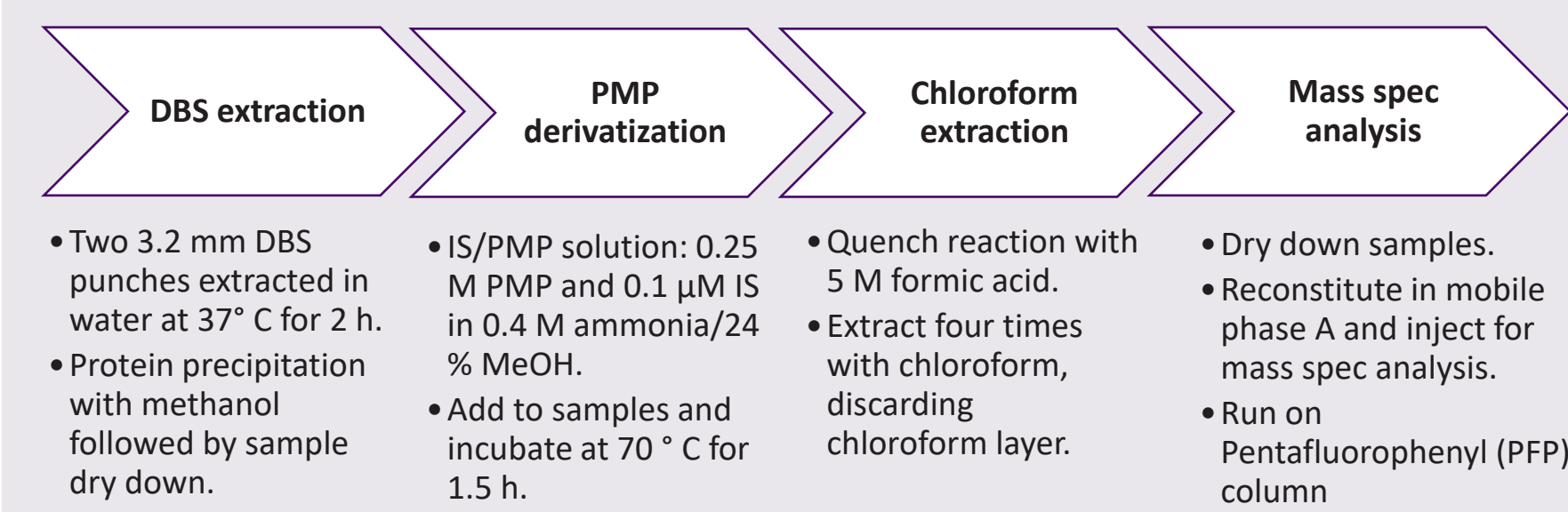
## INTRODUCTION

- Mucopolysaccharidosis (MPS) disorders are a category of lysosomal storage diseases that result in a disruption of the catabolism of glycosaminoglycans (GAGs), macromolecules consisting of long polysaccharide chains. There are a total of seven MPS disorders, with MPS I and II on the (Recommended Uniform Screening Panel) RUSP.
- Primary screening for MPS diseases is accomplished through enzyme activity testing<sup>1</sup>, however, additional methods are required to identify false positives.
- Second tier tests have traditionally relied on elevations of broad classes of GAGs such as dermatan sulfate and keratan sulfate.
  - This approach cannot differentiate between MPS I and MPS II disorders.
- Recent discovery of terminal non-reducing fragments cleaved from GAGs within affected patients<sup>2</sup> presented us with an opportunity to investigate the utility of measuring these markers in specific MPS disease subtype.

## MATERIALS AND METHODS

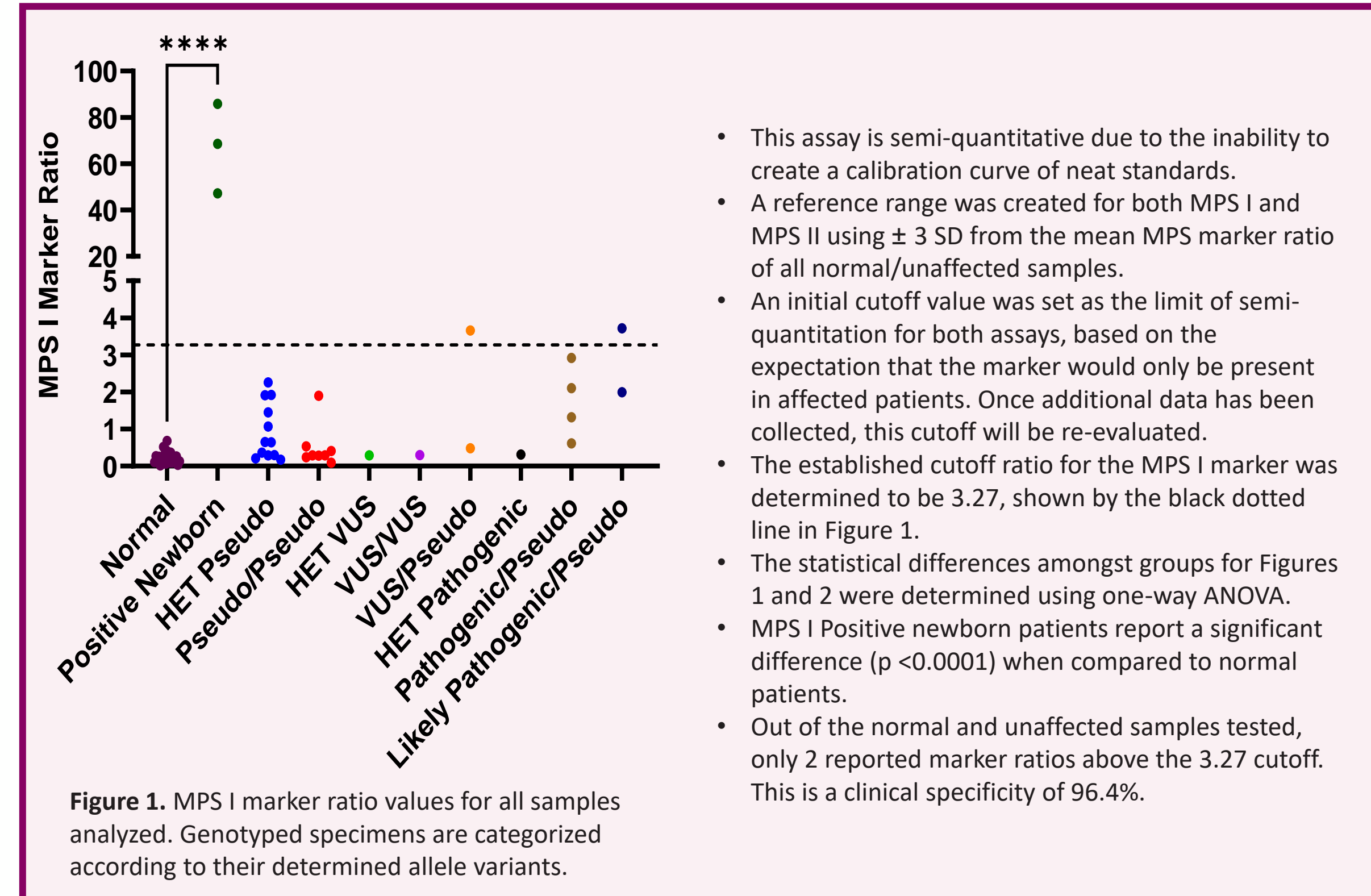
- MPS I
  - Whole blood DBS enriched with fibroblast generated MPS I biomarker served as the positive control (GelbChem)
  - Presumed normal whole blood DBS served as the negative control
- MPS II
  - MPS II positive patient urine served as the positive control during validation studies carried out and shown in Figure 1.
  - Whole blood DBS enriched with fibroblast generated MPS II biomarker was implemented as the positive control for clinical samples. (GelbChem)
  - Presumed normal whole blood DBS served as the negative control.
- Both biomarkers used Chondroitin disaccharide-d4 as the IS (Cayman Chemical) and 1-phenyl-3-methyl-5-pyrazolone (PMP) (Sigma-Aldrich) as the derivatizing agent.

MPS I Patient set	MPS II Patient set
<ul style="list-style-type: none"> <li>25 apparently normal</li> <li>27 MPS I pseudo-deficient or carrier</li> <li>4 MPS I VUS</li> <li>3 known MPS I positive</li> </ul>	<ul style="list-style-type: none"> <li>25 apparently normal</li> <li>4 known MPS II positive in treatment</li> <li>3 known MPS II positive not in treatment</li> </ul>



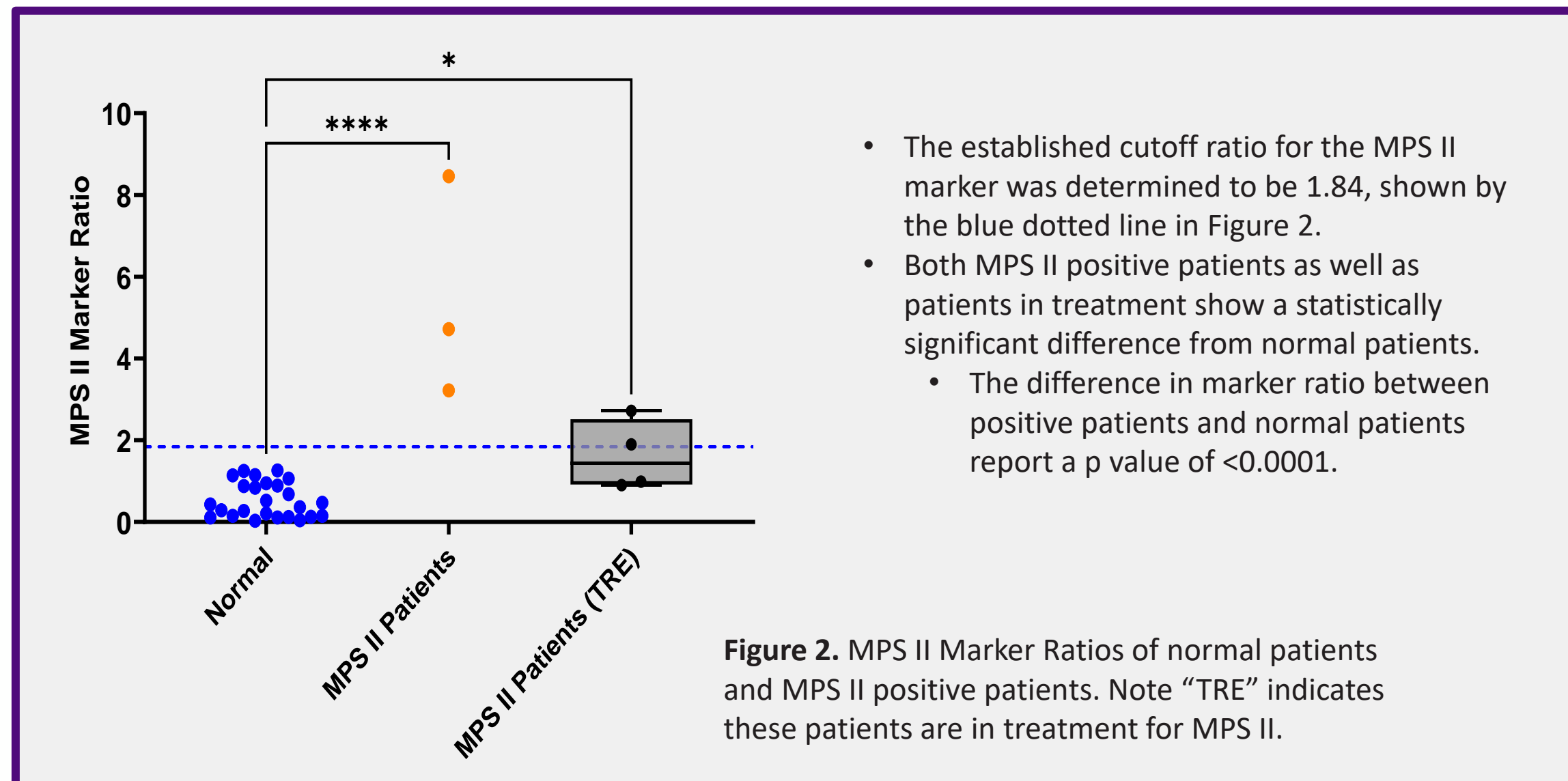
Sample preparation methods and instrument parameters were designed from methods previously reported by Herbst *et al.*<sup>3,4</sup>

## MPS I MARKER



- This assay is semi-quantitative due to the inability to create a calibration curve of neat standards.
- A reference range was created for both MPS I and MPS II using  $\pm 3$  SD from the mean MPS marker ratio of all normal/unaffected samples.
- An initial cutoff value was set as the limit of semi-quantitation for both assays, based on the expectation that the marker would only be present in affected patients. Once additional data has been collected, this cutoff will be re-evaluated.
- The established cutoff ratio for the MPS I marker was determined to be 3.27, shown by the black dotted line in Figure 1.
- The statistical differences amongst groups for Figures 1 and 2 were determined using one-way ANOVA.
- MPS I Positive newborn patients report a significant difference ( $p < 0.0001$ ) when compared to normal patients.
- Out of the normal and unaffected samples tested, only 2 reported marker ratios above the 3.27 cutoff. This is a clinical specificity of 96.4%.

## MPS II MARKER



- The established cutoff ratio for the MPS II marker was determined to be 1.84, shown by the blue dotted line in Figure 2.
- Both MPS II positive patients as well as patients in treatment show a statistically significant difference from normal patients.
  - The difference in marker ratio between positive patients and normal patients report a p value of  $< 0.0001$ .

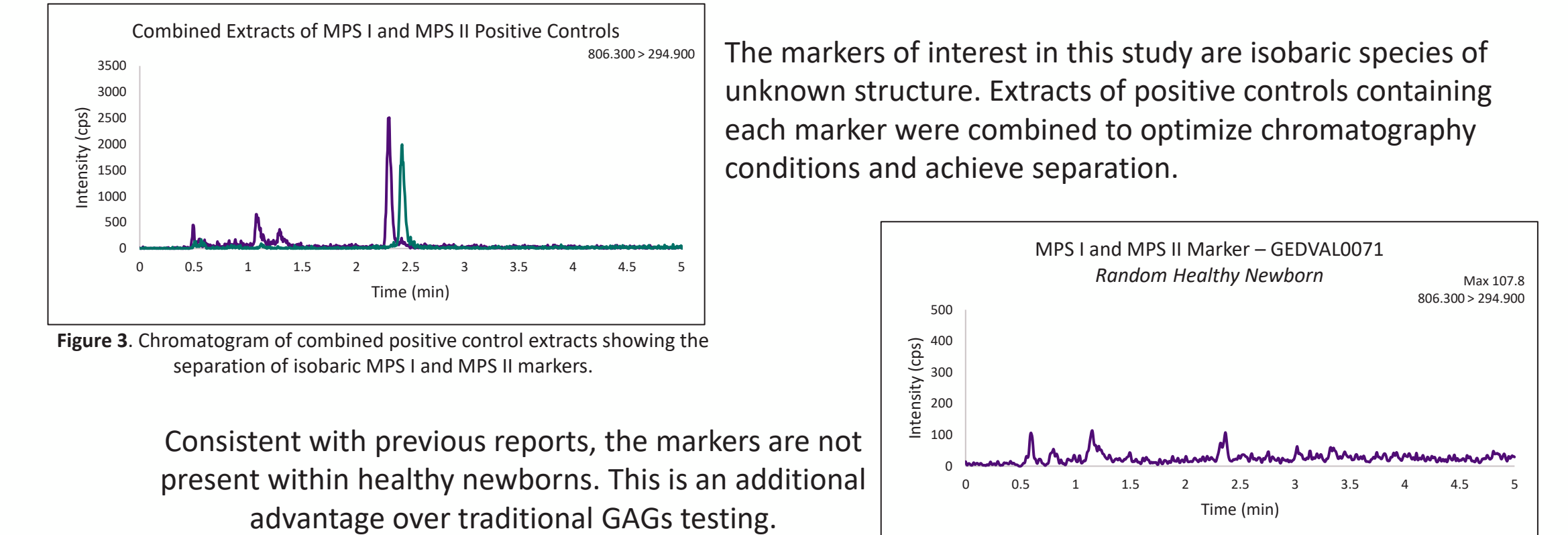
## ACKNOWLEDGEMENTS

We would like to thank GelbChem for the supplied DBS control and the support of Michael Gelb, Zackary Herbst, and Hamid Khaledi throughout this project.

- Peck *et al.*, *Int. J. Neonatal Screen*, **2020**, 6 (1), 10.
- Saville *et al.*, *Genetics in Medicine*, **2019**, 21 (3), 753-757.
- Herbst *et al.*, *Int. J. Neonatal Screen*, **2020**, 26 (6), 69.
- Herbst *et al.*, *Int. J. Neonatal Screen*, **2022**, 21 (8), 9

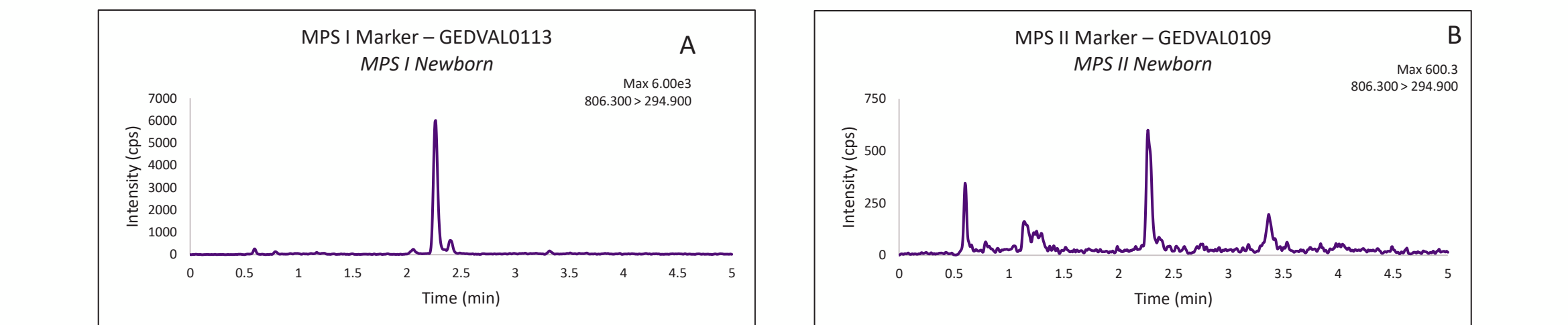
Pursuant to applicable federal and/or state laboratory requirements, Revvity Omics establishes and verifies the accuracy and precision of their testing services.

## SPECIFICITY OF MPS I & II MARKERS



Consistent with previous reports, the markers are not present within healthy newborns. This is an additional advantage over traditional GAGs testing.

- MPS I positive newborn samples display a large response for the MPS I marker. We have consistently observed a small signal at the MPS II marker retention time within MPS I positive patients, however this peak is resolved from the true MPS I marker peak and is well below the threshold for the MPS II marker.
- We have not observed the presence of any signal at the MPS I retention time within MPS II affected patients.



**Figure 5.** Chromatograms from newborns affected by (A) MPS I and (B) MPS II.

## CONTROL STABILITY

MPS Control	Timepoint	Marker Ratio	% Difference from t = 0	Pass/Fail
MPS I Positive Control -80 °C	t = 0	18.84		
	t = 3 weeks	19.95	5.87	Pass
	t = 1 month	21.08	11.89	Pass
	t = 2 months	17.16	8.94	Pass
	t = 3 months	16.41	12.92	Pass
MPS II Positive Control -80 °C	t = 0	31.90		
	t = 3 weeks	31.90	0.00	Pass
	t = 1 month	28.20	11.60	Pass
	t = 2 months	24.60	22.90	Pass
	t = 3 months	25.20	21.00	Pass
	t = 6 months	30.10	5.6	Pass

**Table 1.** Shows MPS I and MPS II control long-term stability data collected thus far. Samples were considered passing if the calculated marker ratio was  $< 25\%$  difference from time = 0.

## CONCLUSION AND NEXT STEPS

- Our validated methods allow for increased specificity between MPS I and MPS II disease subtypes from measuring a characteristic fragment of accumulated GAGs compound.
- The MPS I marker test was performed retroactively on a set of samples that tested low for IDUA during primary newborn screening yet were confirmed unaffected by gene sequencing. The MPS I marker was found to be within normal limits for all 28 samples tested in this set, further supporting the benefit of performing this second-tier test as part of a newborn screening algorithm.
- Continuation of the MPS I allele study to further elucidate any possible correlation with genotype and MPS I marker ratio values.
- An allele and biomarker correlation study for MPS II.
- Validation of MPS I and MPS II biomarkers in urine samples.