

PETH IMMUNOASSAY: A SEMI-QUANTITATIVE ELISA FOR ASSESSMENT OF ALCOHOL CONSUMPTION

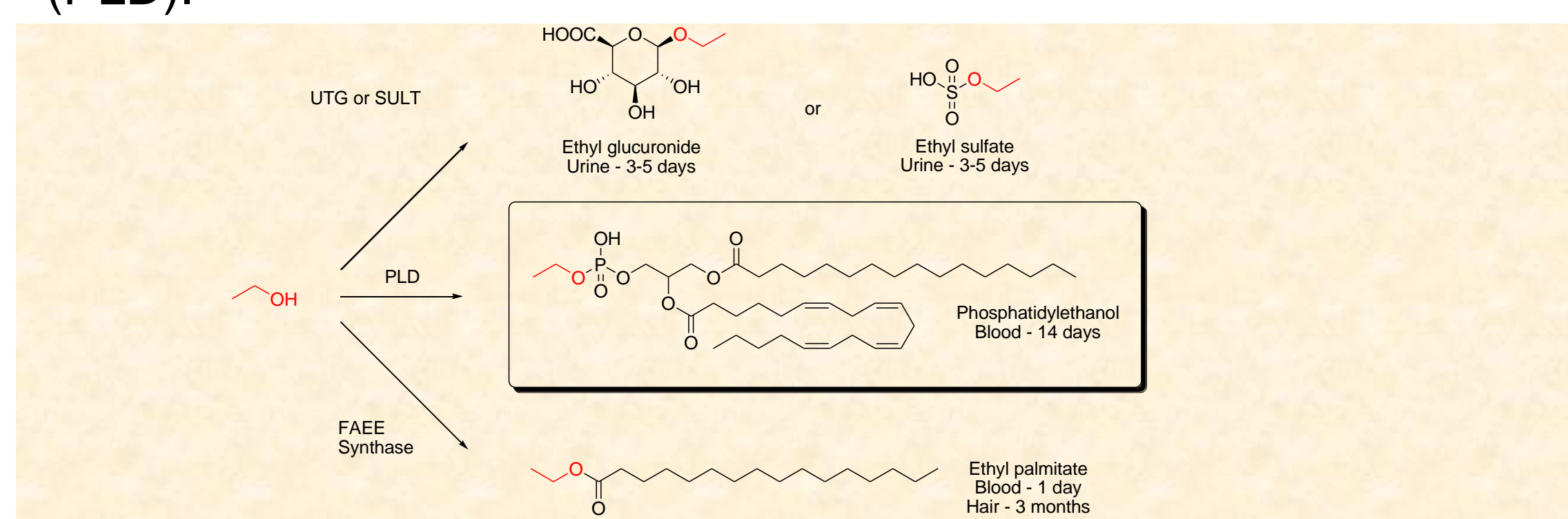
L. Jeffrey Johnson¹, Fok Yun Chan¹, Winnie R. Muyindike², Cameron Day¹, Judith A. Hahn³, Paul O. Neilsen¹
 Echelon Biosciences, Salt Lake City, UT¹; Mbarara University of Science and Technology²;
 University of California, San Francisco³

Overview

- Phosphatidylethanol (PEth) is accepted as a long-term direct biomarker for the detection and assessment of alcohol consumption.
- Utilizing a Semi-Quantitative PEth ELISA we analyzed a set of 120 clinical samples, compared four alcohol consumption categories, and identified a significant correlation between ELISA values and reported mass spectrometry (MS) values.
- Three blood collection devices were evaluated and confirmed to be compatible with the Semi-Quantitative PEth ELISA.

Introduction

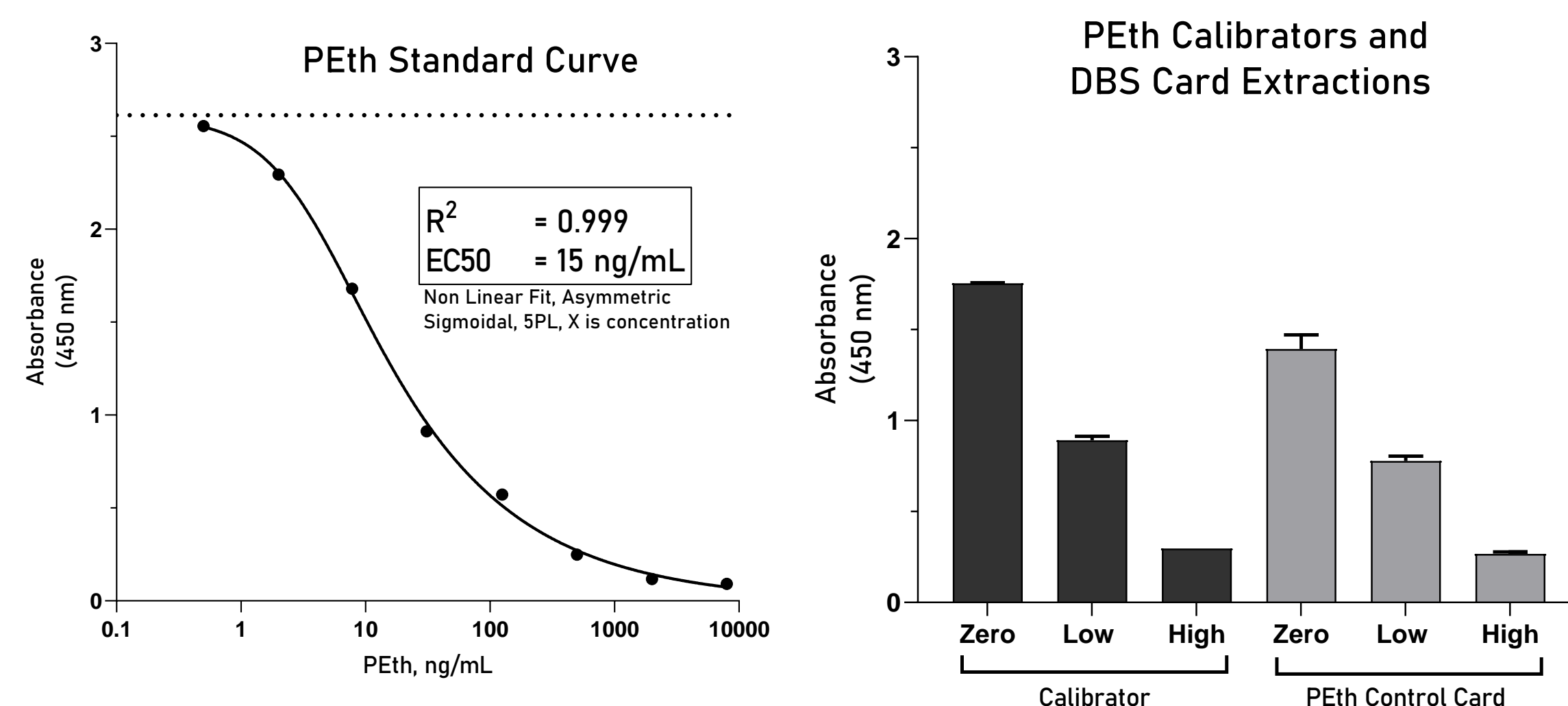
- Abuse and overuse of alcohol contributes to roughly 5% of the overall global health burden.
- Monitoring excessive binge and chronic drinking will benefit individual health outcomes as well as society at large.
- PEth is a phospholipid formed in the presence of EtOH by transacylation of Phosphatidylcholine (PC) by Phospholipase D (PLD).



- PEth is quite stable when stored, extracted, and measured from dried blood, allowing for flexibility in the collection process.
- We present the first immunoassay that can measure PEth levels in a semi-quantifiable manner to supplement LC/MS analysis.
- Our goal is to further simplify clinical laboratory processes, reduce time & cost, and expand the available options for clinical analysis of PEth.

Semi-Quantitative PEth ELISA

- Competitive Semi-Quantitative Immunoassay utilizing a synthetic 16:0/18:1 PEth standard curve, allowing the user to assign samples a PEth concentration (ng/mL).
- Free PEth in solution, either extracted from samples, controls, or provided as a standard, competes with the PEth ab against PEth bound to the wells of a 96 well plate.
- Decreased PEth ab binding to the plate indicates the presence of PEth. A lower signal indicates higher levels of PEth.

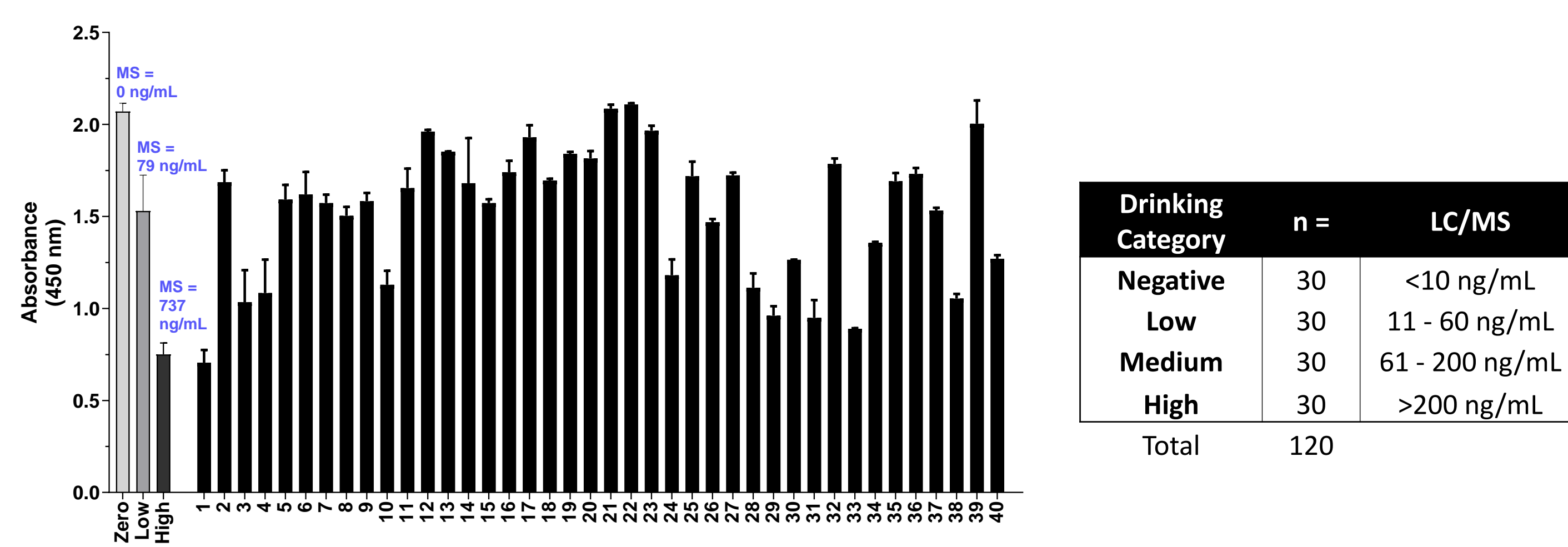


Sample Type	Reported LC/MS Value	Semi-Quantitative PEth ELISA (Interpolated Value)	
		Calibrator	DBS Card Extraction
Zero	< 8 ng/mL	7 ng/mL	13 ng/mL
Low	79 ng/mL	37 ng/mL	50 ng/mL
High	737 ng/mL	414 ng/mL	521 ng/mL

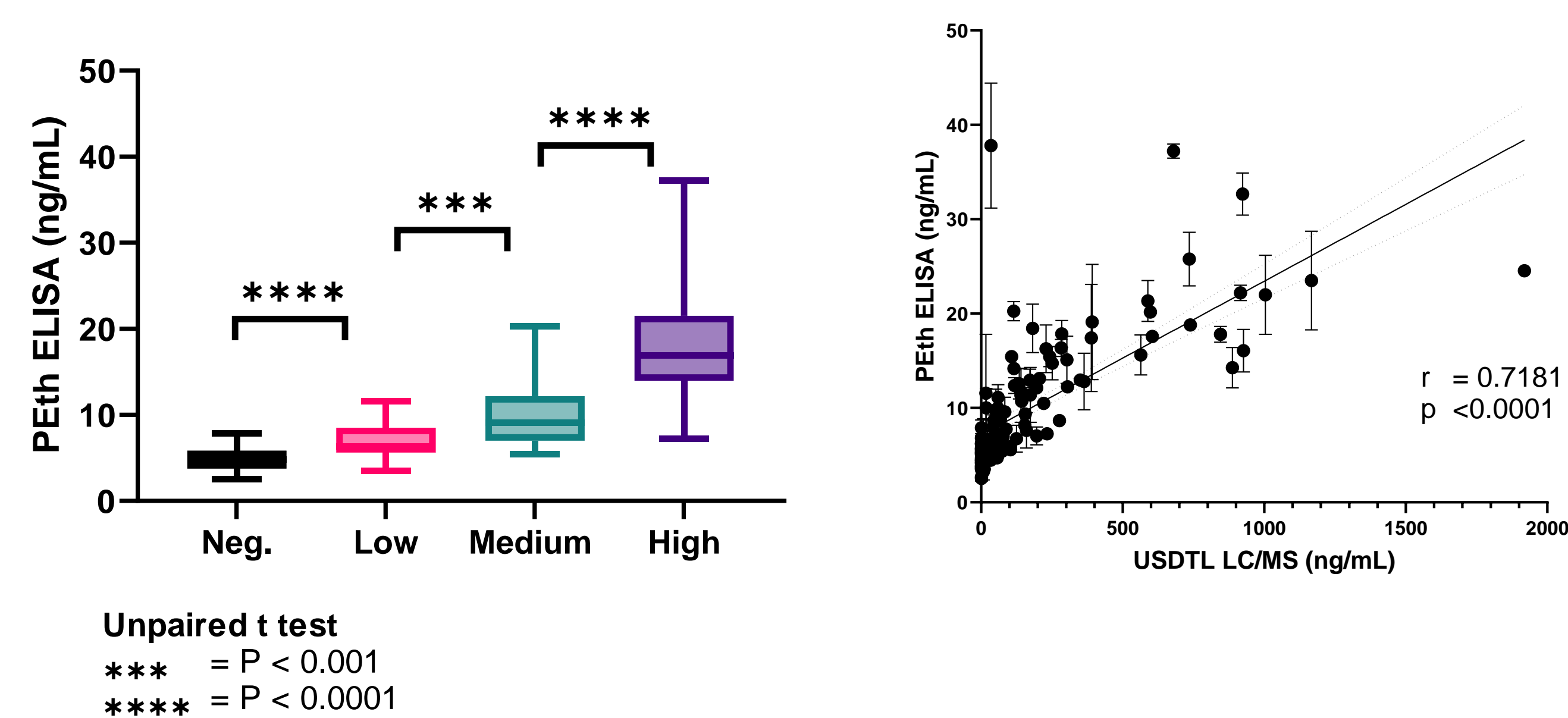
*The observed PEth ELISA concentrations (ng/mL) are correlative but are not equivalent to their LC/MS counterparts. This is likely due to differential matrix effects.

Analysis of 120 Clinical Samples

PEth ELISA Quantifies PEth levels in clinical samples



Strong correlation between ELISA and MS PEth concentrations

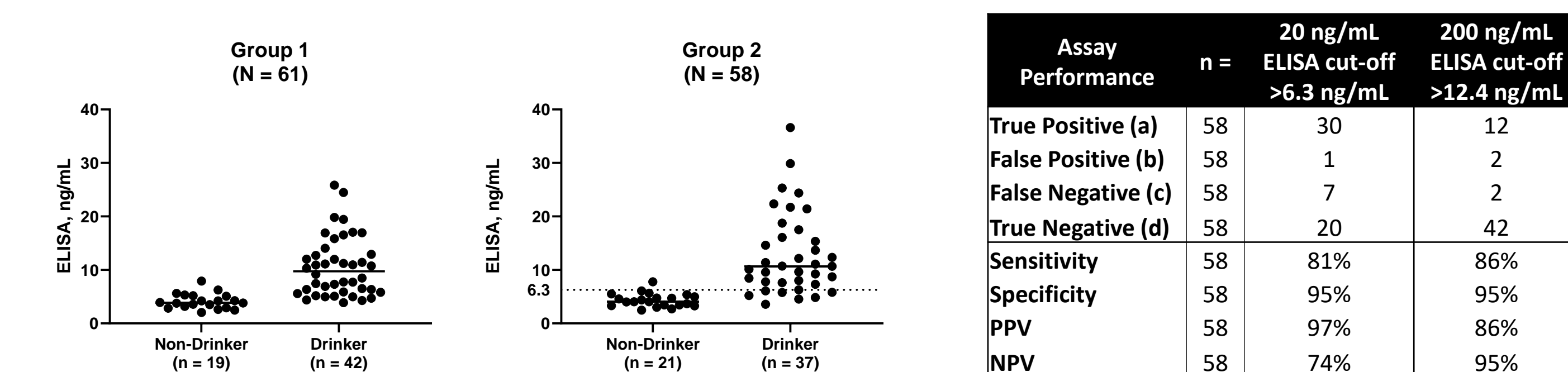


Sensitivity, Specificity, and Determination of Cut-Off Values

ELISA Cut-Off Values are contingent on specificity and sensitivity. Our goal was to define the ELISA limits for testing both Abstinence and Excessive alcohol use.

Approach:

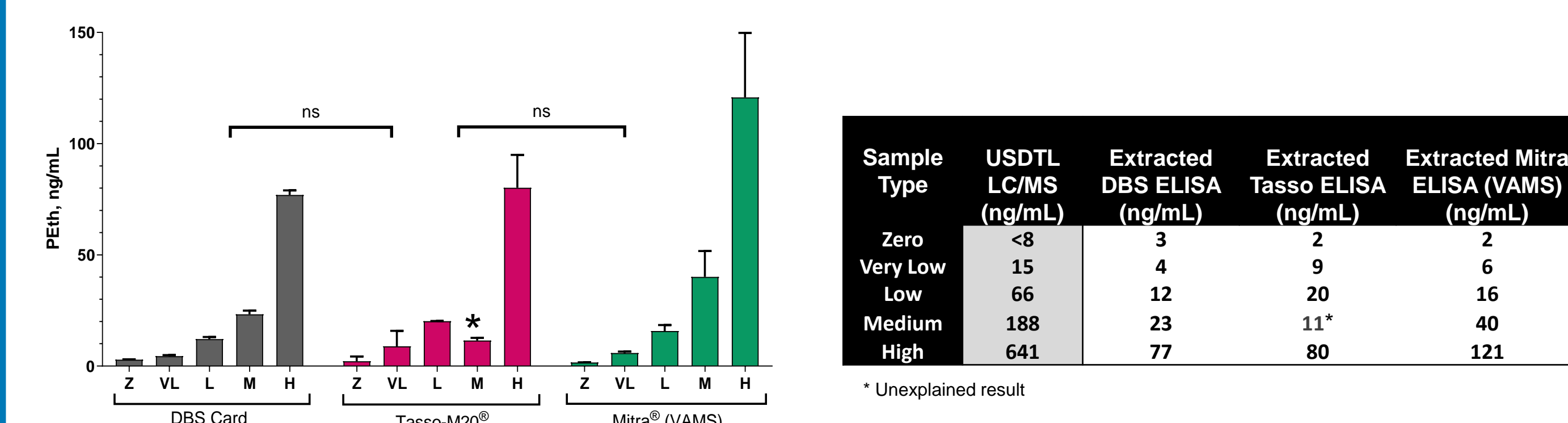
- We randomized half of the clinical samples (n=61) and created two groups, non-drinkers (n=27) & drinkers (n=34), defined by MS values less than or greater than 20 ng/mL for the first analysis. For the second we used 200 ng/mL.
- Using ROC (receiver operating characteristics) and a maximized likelihood ratio to determine sensitivity and specificity, we established ELISA cut-off values of >6.3 ng/mL for the 20 ng/mL, and >12.4 ng/mL for the 200 ng/mL groups.
- We analyzed the ELISAs performance using the second half of the samples (n=58) with the above determined cut-off values to define values for assay sensitivity, specificity, positive, and negative predictive values (PPV & NPV).



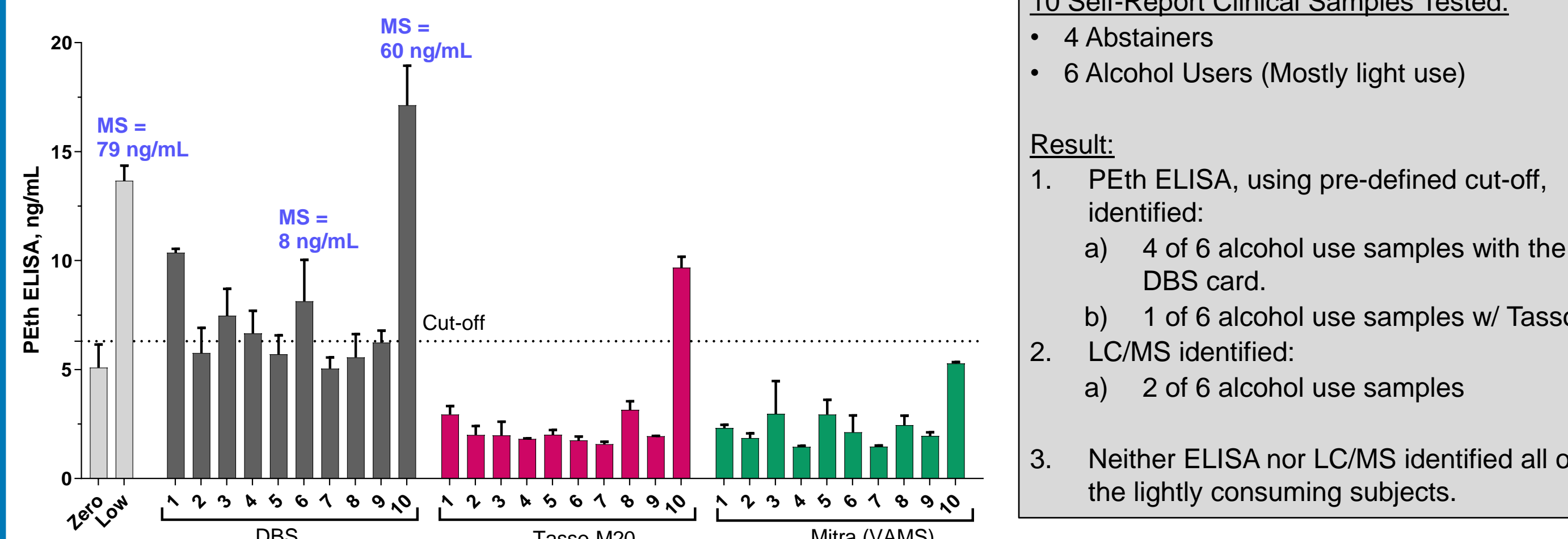
Comparison of Blood Collection Devices

Blood Collection Devices:	Whole blood equivalent	Sample Type
1. Whatman® 903 DBS Card	(15 µL, 5 x 3.2 mm punches)	Capillary and venipuncture
2. Tasso-M20® Device	(17.5 µL eq.)	Capillary and venipuncture
3. Mitra® Device (VAMS®)	(20 µL eq.)	Capillary and venipuncture

ELISA differentiates between five ex vivo generated PEth controls and displays no significant difference with blood collection devices

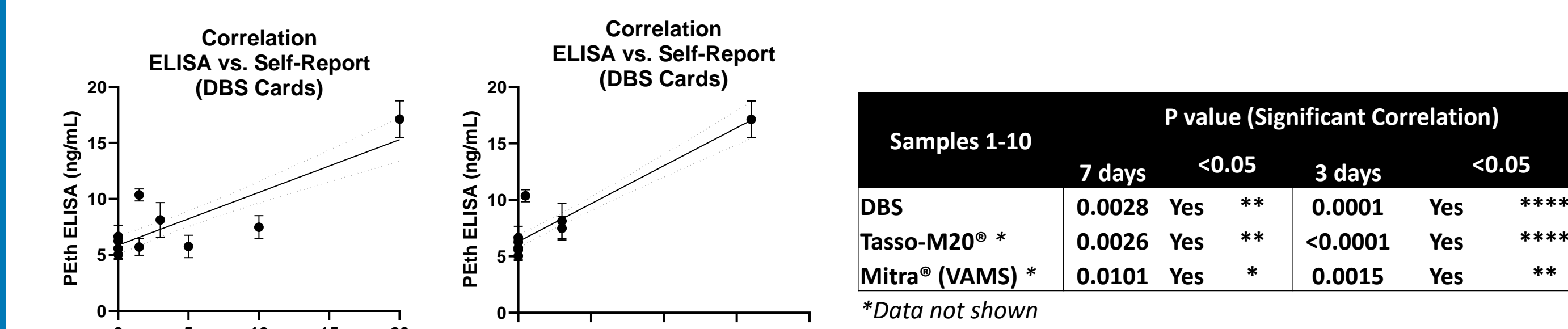


ELISA and MS analysis of Self-Reported Alcohol Use Samples are Comparable



Clinical Samples	n	Positive		Negative		Positive		Negative	
		w/ DBS	w/ DBS	w/ DBS	w/ DBS	w/ Tasso	w/ Tasso	w/ Mitra	w/ Mitra
Non-Drinking	4	0	4	1	3	0	4	0	4
Drinking	6	2	4	4	2	1	5	0	6
Total	10	2	8	5	5	1	9	0	10

ELISA Alone Correlates with Self-Report of Alcohol Use



- The Tasso-M20® and Mitra®(VAMS®) further simplify the blood collection process and are compatible with the PEth ELISA.
- The PEth ELISA identified drinking in multiple samples with DBS Cards and in 1 of 6 drinkers in all three collection devices.
- As expected, ELISA results demonstrate stronger correlation to self reported drinking when drinking events are closer to the analysis date.

Conclusions, Future, & Acknowledgements

Conclusions

- The Semi-Quantitative PEth ELISA:
 - Differentiates between four alcohol consumption categories in a quantitative fashion.
 - Demonstrates significant correlation with reported LC/MS values.
 - Is compatible with the Tasso-M20® and Mitra®(VAMS®) collection devices.
- This 96-well immunoassay expands the availability of PEth analysis and is poised to become a valuable clinical tool in screening large numbers of samples in high-throughput workflows.

Future Directions:

- Explore additional extraction methods and assay formulations to refine assay sensitivity and specificity.
- Evaluate the stability of samples collected and stored in different conditions & increase assay ruggedness and robustness.

The authors wish to thank:

- Funding: NIH R01AA024990 (funding to support Hahn and Muyindike to collect the 120 clinical samples)
- The Echelon Team for support, suggestions, and analysis
- Colin Ferguson and Bo Ye for synthesizing phospholipid analogs
- Study volunteers and blood donors