

Analytical Validation of an Ultrasensitive Neuroinjury Multiplex Assay Panel

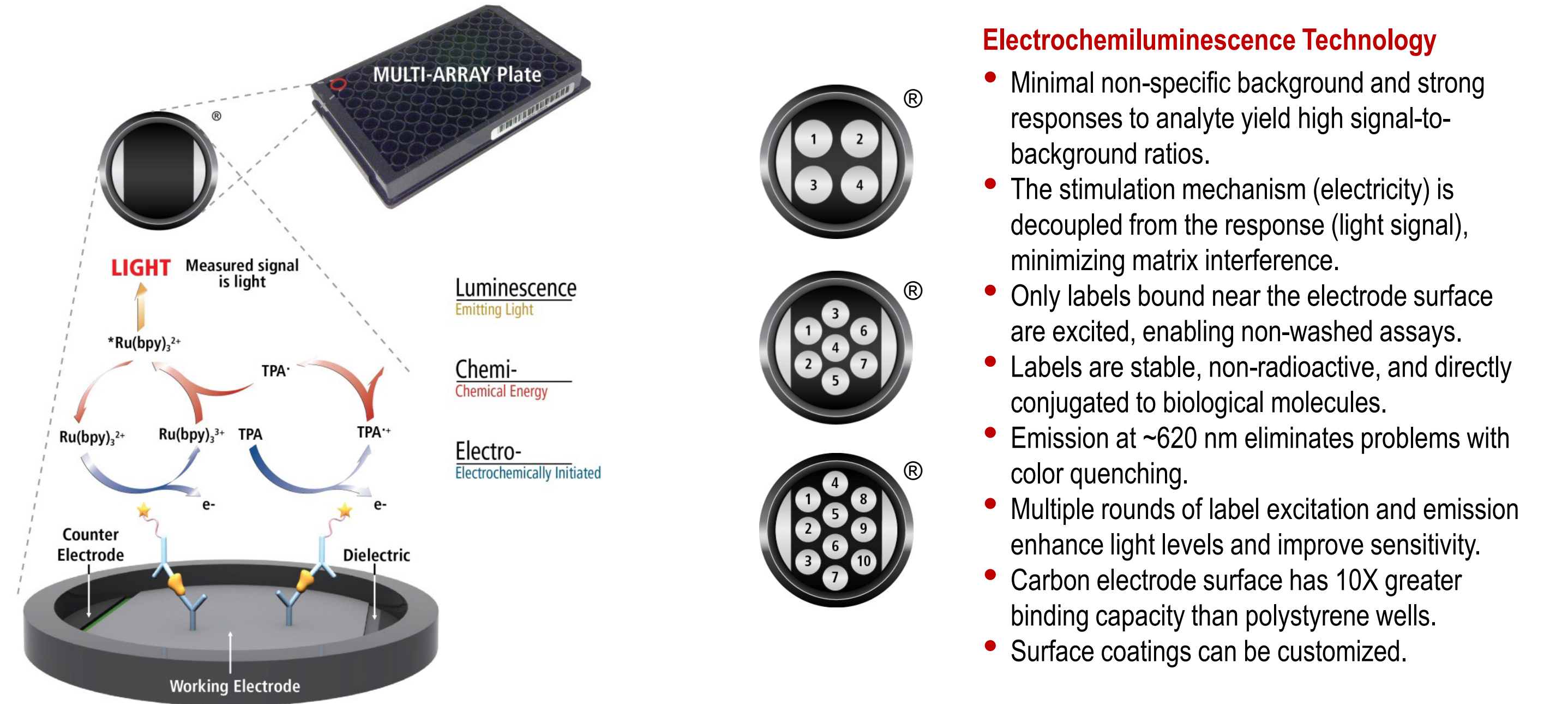
Catherine Demos, Daniel Romero, Nikhil Padmanabhan, Rachel Cohen, Jermaine Brown, Christopher Campbell, Martin Stengelin, Anu Mathew, George Sigal, and Jacob Wohlstadter
Meso Scale Diagnostics, LLC., Rockville, Maryland, USA

1 Abstract

Three blood-based biomarkers of neuroinjury—glial fibrillary acidic protein (GFAP), neurofilament light (NF-L), and total Tau (tTau)—have emerged as promising biomarkers for research regarding neurological disorders and neuroinjuries such as hypoxic-ischemic encephalopathy (HIE), traumatic brain injury (TBI) and Alzheimer's disease (AD). The low levels of GFAP, NF-L, and tTau in serum and plasma require highly sensitive assays to detect these important biomarkers. Here, we report on an ultrasensitive, electrochemiluminescence-based, multiplexed immunoassay for GFAP, NF-L, and tTau that has been analytically validated, providing a new tool for neuroinjury biomarker research. The MSD S-PLEX Neurology Panel 1 kit uses ultrasensitive S-PLEX assay technology to simultaneously measure GFAP, NF-L, and tTau in a 96-well plate format using standard liquid-handling techniques. Analytical validation was carried out in a series of studies based in part on Clinical and Laboratory Standards Institute guidelines EP05-A3, EP06-Ed2, EP07-Ed3, EP17-A2, EP25-A, and EP28-A3c to evaluate precision, dilution linearity, interference screening, detection capability, stability, spike recovery, and cross reactivity. The assay requires 5 µL of a sample for each replicate. The time to result is less than 6 hours and the quantifiable range of the assay is 2.4–4,230 pg/mL for GFAP, 7.65–10,000 pg/mL for NF-L, and 0.34–735 pg/mL for tTau. A set of 15 common potentially interfering substances were screened in serum and plasma, and none showed interference exceeding 18% difference in measurement compared to non-spiked samples. Within-laboratory precision was <11% CV for GFAP and NF-L and <19% CV for tTau for samples spanning the reportable range. Dilution up to 256-fold recovered within 80–120% of the expected value. MSD has developed and analytically validated a highly sensitive, multiplexed immunoassay for quantifying GFAP, NF-L, and tTau in plasma and serum. This assay is sufficiently sensitive to detect these analytes in normal serum and plasma and has a large dynamic range to accommodate elevated levels found in some neurological disorders.

2 Methods

MSD's electrochemiluminescence detection technology uses SULFO-TAG™ labels that emit light upon electrochemical stimulation initiated at the electrode surfaces of MULTI-ARRAY® and MULTI-SPOT® microplates. The MSD® S-PLEX Neurology Panel 1 is a three analyte ultrasensitive panel. The S-PLEX® platform uses ECL technology, retaining its well-known advantages and superior analytical performance. The improved sensitivity of S-PLEX assays is due to the new TURBO-TAG® and TURBO-BOOST® reagents. This assay was validated using a custom protocol designed to reduce time to results. Incubation times are reduced by 30% compared to the commercial kit protocol and are performed at +27°C.



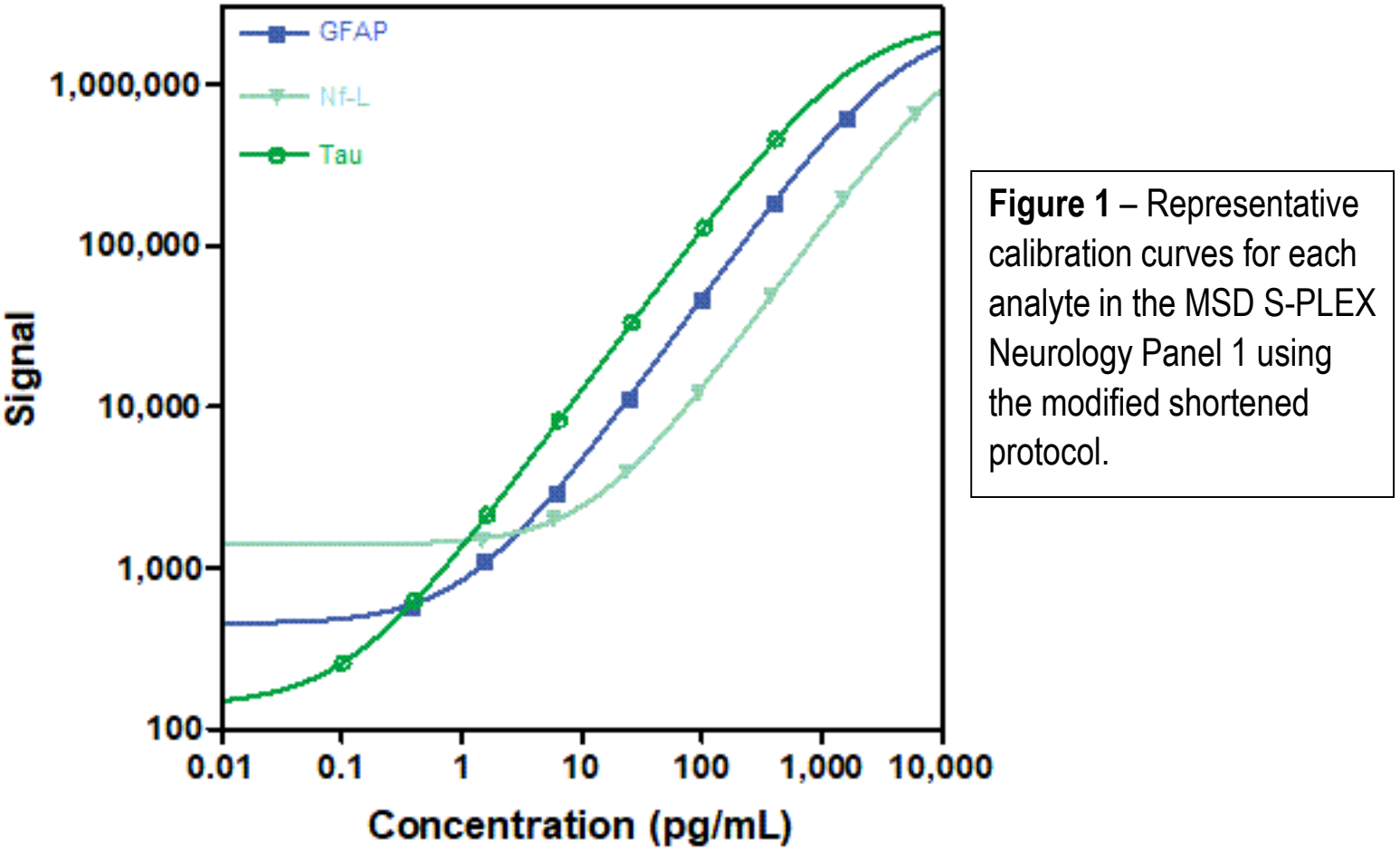
- Electrochemiluminescence Technology**
- Minimal non-specific background and strong responses to analyte yield high signal-to-background ratios.
 - The stimulation mechanism (electricity) is decoupled from the response (light signal), minimizing matrix interference.
 - Only labels bound near the electrode surface are excited, enabling non-washed assays.
 - Labels are stable, non-radioactive, and directly conjugated to biological molecules.
 - Emission at ~620 nm eliminates problems with color quenching.
 - Multiple rounds of label excitation and emission enhance light levels and improve sensitivity.
 - Carbon electrode surface has 10X greater binding capacity than polystyrene wells.
 - Surface coatings can be customized.

3 Results

A. Detection Capability

Table 1 – Limits of detection and quantification for the S-PLEX Neurology Panel 1 assay using the modified shortened protocol.

Limits of Detection and Quantification				
Conc. (pg/mL)	LOB	LOD	LLOQ	ULOQ
GFAP	1.72	2.05	2.30	4,230
NF-L	3.23	4.99	7.65	10,030
tTau	0.255	0.324	0.687	735



Quantification of the analytes was determined by reference to an eight point calibration curve (Figure 1) run on each plate and fit to a four parameter logistic (4PL) model. The limits of quantification were determined according to CLSI guidelines EP17-A2 by measuring four high level, eight low level, and four blank samples in quintuplicate on three plates each for two reagent lots. The total error (TE) was calculated via the root mean squared model for each sample on each reagent lot and the TE cutoff value was ≤ 25%.

- Lower limit of quantification (LLOQ) was determined as the lowest concentration sample where the TE ≤ 25%, and the higher value between the two reagent lots.
- Upper limit of quantification (ULOQ) was determined similarly as the highest concentration with TE ≤ 25%, and the lower value between the reagent lots.
- The limit of blank (LOB) was determined by the measurement of four different blank samples on two reagent lots. The measurements of the blank samples were rank-ordered and LOB was determined as the 95th percentile of the calculated concentrations.
- The limit of detection (LOD) was determined by the measurement of four different blank samples and four different low level samples on two reagent lots. LOD is the minimum concentration such that 95% of measurements of a sample with concentration equal to LOD will exceed the LOB.

B. Clinical Range – Alzheimer's Diseased Samples

- To demonstrate the ability of the assays to measure the relevant concentration ranges of the target analytes in plasma and CSF, healthy and diseased matched plasma and CSF samples were tested on the MSD S-PLEX Neurology Panel 1.
- All samples were within the limits of quantification. Data are shown in Figure 2.

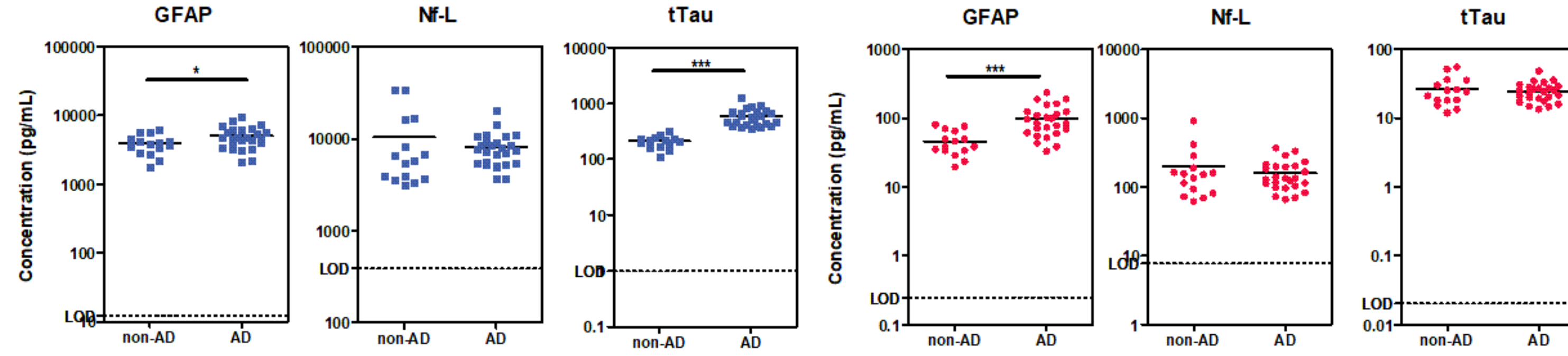


Figure 2 – Sample concentrations of AD (n=25) vs non-AD (n=15) CSF (left, blue) and EDTA plasma (right, red). Samples were tested blinded and in duplicate. Mean intra-plate CV was <5% for all analytes. Dashed lines indicate dilution-adjusted LOD (100-fold for CSF and 2-fold for EDTA Plasma). Differences of means were calculated by student's t-test. * = p-value ≤ 0.05, *** = p-value ≤ 0.0001

C. Dilution Linearity

Dilution linearity of the MSD S-PLEX Neurology Panel 1 was determined in three stages adapted from the CLSI guidelines within EP06-A2.

- Linearity was shown to be within 80–120% over a 256-fold dilution range by testing serially diluted pooled adult human serum and plasma that had been spiked with calibrator at the upper end of the quantitative range. Representative data are shown in Figure 3.
- Dilution linearity was verified using 64 individual neonatal samples diluted 5- (recommended dilution), 10- and 20-fold. Further dilutions were not tested due to the low sample volume available. Summary NF-L results for all donors are shown in Figure 4. GFAP, NF-L, and tTau are all observed to dilute linearly within 20%.
- Dilution linearity of neonatal serum samples was tested by preparing "high" and "low" pools of HIE samples into a 5-sample panel using equal-volume mixing. All analytes showed linearity within 20%. Representative data are shown in Figure 5.

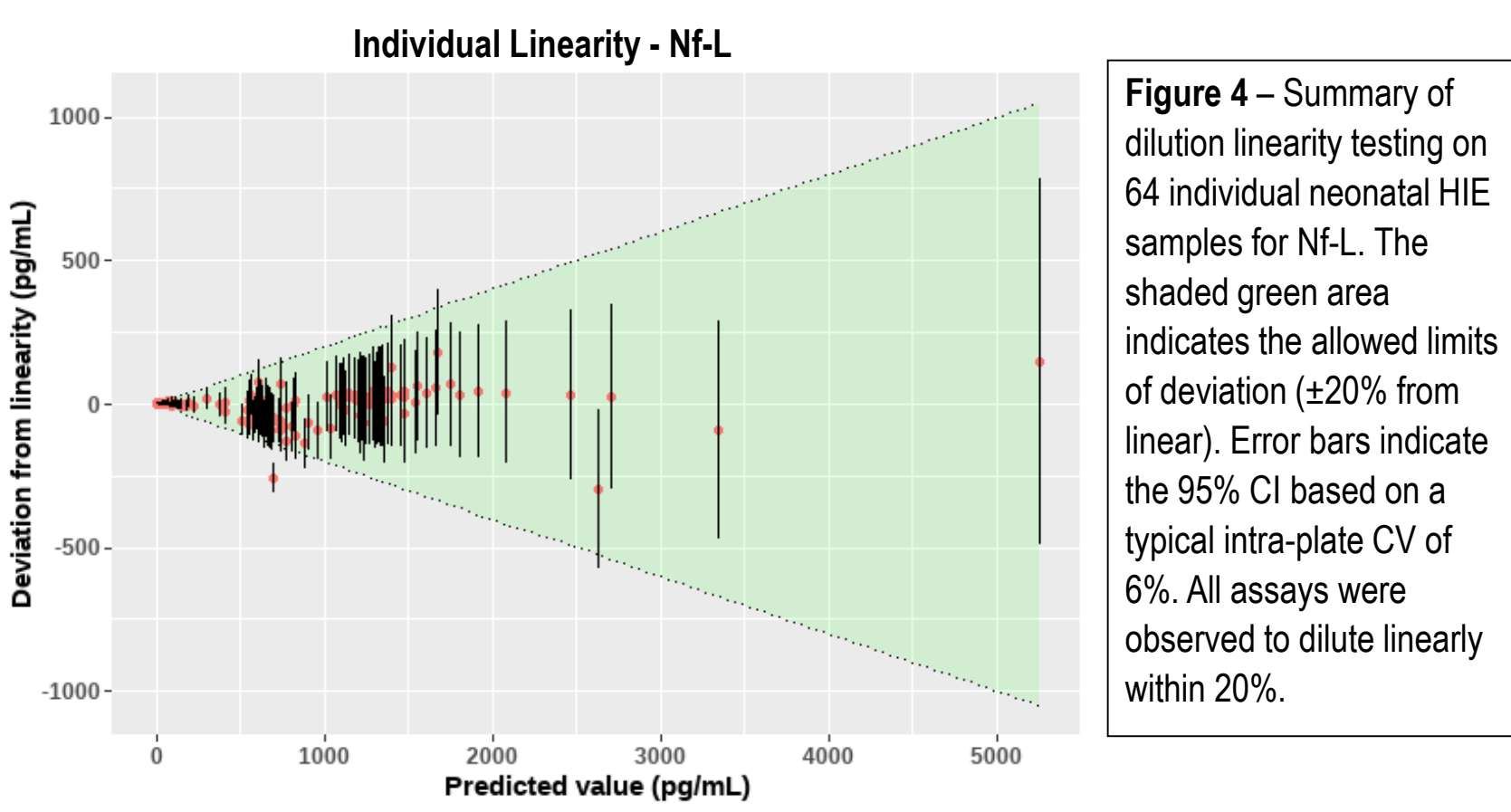
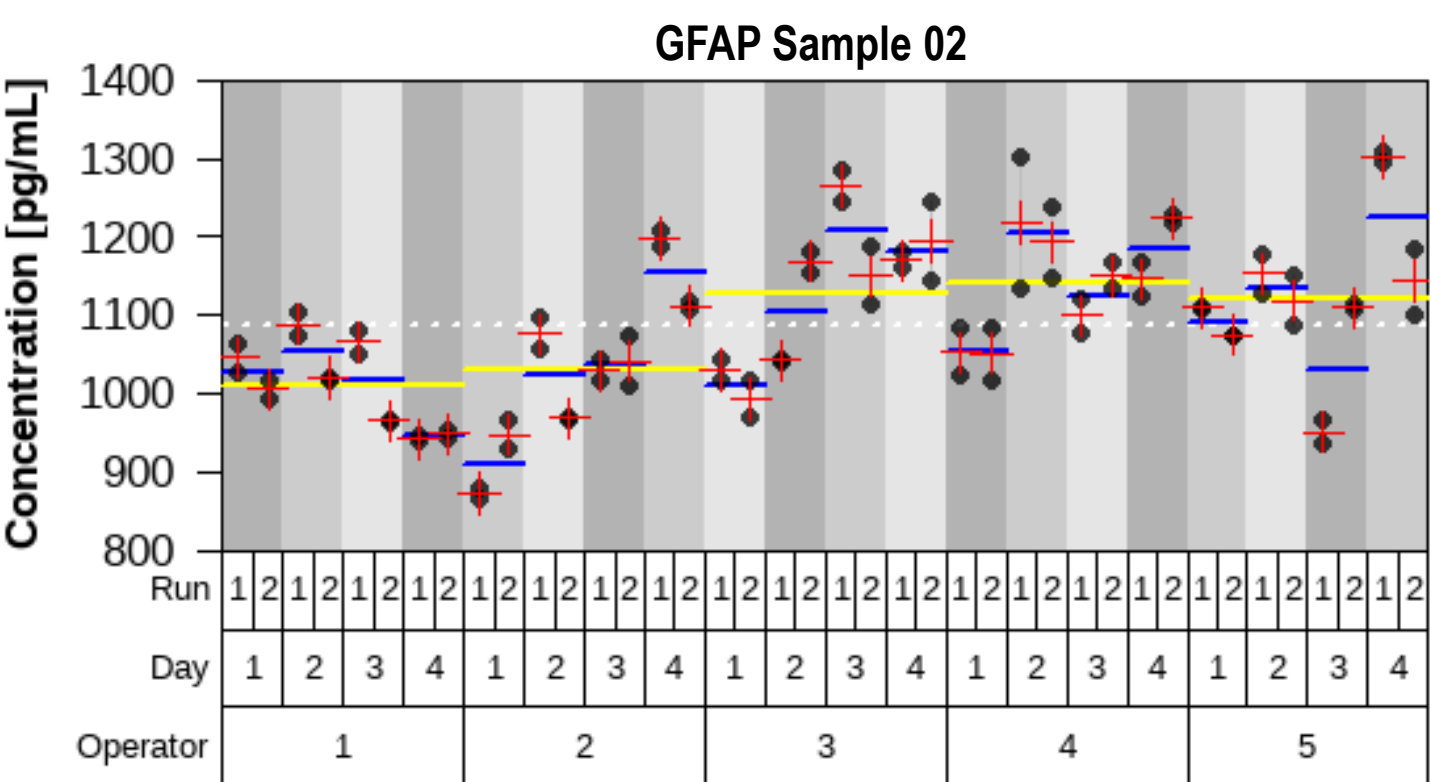


Figure 4 – Summary of dilution linearity testing on 64 individual neonatal HIE samples for NF-L. The shaded green area indicates the allowed limits of deviation (±20% from linear). Error bars indicate the 95% CI based on a typical intra-plate CV of 6%. All assays were observed to dilute linearly within 20%.

D. Precision

Table 2 – Precision estimates for components of potential variability.

%CV	Repeatability (Within-Run/Residual Error)	Between-Operator	Between-Day	Between-Run	Within Laboratory Precision (Total)
GFAP	4.0%	5.2%	3.6%	7.3%	10.9%
NF-L	5.9%	2.1%	3.8%	6.0%	10.2%
tTau	4.8%	11.4%	3.1%	13.5%	19.3%



- A precision study was designed based on CLSI Guideline EP05-A3 and modified to assess precision within laboratory (total) and between operator, day, run, and replicate (repeatability). Briefly, 10 samples (five EDTA plasma, five serum) were spiked with a blended calibrator preparation at five levels spanning the expected range and vialied in single-use frozen aliquots. Five analytical controls spanning the expected range were lyophilized in single-use vials. Five operators each performed four days of testing with two simultaneous, vial-independent runs per operator per day, spanning a total of 20 calendar days, resulting in 80 measurements per sample overall.
- For all samples within the limits of quantification, all assays achieved a total %CV ≤ 20%. Detailed results for a representative sample are shown in Figure 6, and precision estimates for the entire quantitative range are listed in Table 2. The residual error (error between replicates) is <6% CV for all analytes, and the total within-laboratory precision was 11% for GFAP, 10% for NF-L, and 19% for tTau.

Figure 6 – Representative measurements of one sample from the precision study. The mean concentration between replicates (red cross), between runs per day (blue line) and between operators (yellow line) are plotted against individual measurements. Variance component analysis was completed with R using the package 'VCA.'

E. Interference Screening

Fifteen common laboratory and medical substances listed below were screened for interference with the MSD S-PLEX Neurology Panel 1. Samples spiked with the maximum clinically relevant concentration or a mock spike were compared, and none showed interference exceeding 18% in paired-difference testing. An example summary of interferent screening for the ampicillin effect on GFAP is shown in Table 3.

- Ampicillin
- Morphine
- Benzodiazepine
- Phenobarbital
- Morphine
- Erythropoietin
- Phenytoin
- Piperacillin
- Bilirubin (unconjugated)
- Tazobactam
- Gentamicin
- Topiramate
- Levetiracetam
- Vancomycin
- Melatonin
- Bilirubin (conjugated)

Table 3 – Summary of interferent screening for ampicillin on GFAP. In all cases, the point estimate for the percent difference was below 18%, which led us to conclude that ampicillin is not a significant cause of interference for GFAP at high or low concentrations. In most cases, the 95% confidence interval crossed 0, which indicates that the null hypothesis (i.e., ampicillin does not alter analyte measurements) could not be rejected.

Analyte	Sample Matrix	Analyte Level	Control Concn. (pg/mL)	Test Concn. (pg/mL)	Percent Difference (%)	Interference	95% CI
GFAP	EDTA Plasma	Spiked	1,070	1,020	-4.7	Not Detected	[-9.8, 0.47]
	EDTA Plasma	Unspiked	9.90	10.2	3.2	Not Detected	[-5.5, 12]
	Serum	Spiked	982	960	-2.2	Not Detected	[-6.9, 2.4]
	Serum	Unspiked	17.9	16.9	-5.3	Not Detected	[-13, 2.8]

F. Spike Recovery

Spike recovery was determined first by spiking individual adult serum and plasma samples with three levels of calibrator, then diluting 2-fold (recommended dilution for adult samples). These GFAP, NF-L, and tTau assays are shown to recover the spiked calibrator within 80%–120% at 2x dilution in serum and EDTA plasma, as shown in Table 4. Next, individual neonatal HIE samples (n=64) were used to verify spike recovery. Samples were spiked with <5% overall volume of calibrator and diluted 5- (recommended dilution for neonatal samples), 10- and 20-fold, as shown in Figure 7.

Table 4 – Spike recovery range of each analyte in 2-fold diluted serum and EDTA plasma.

	Spike Level (pg/mL)	Serum (n=6) 2x Diluted		EDTA Plasma (n=6) 2x Diluted	
		Average % Recovery	% Recovery Range	Average % Recovery	% Recovery Range
GFAP	12,500	88%	86–91%	83%	74–98%
	2,500	84%	81–87%	88%	65–109%
	625	84%	80–90%	92%	68–121%
NF-L	3333	129%	113–139%	109%	90–130%
	667	118%	113–123%	110%	99–121%
	167	112%	81–127%	88%	67–116%
tTau	250	96%	92–102%	94%	85–102%
	50	94%	89–97%	91%	80–100%
	12.5	96%	94–98%	90%	92–108%

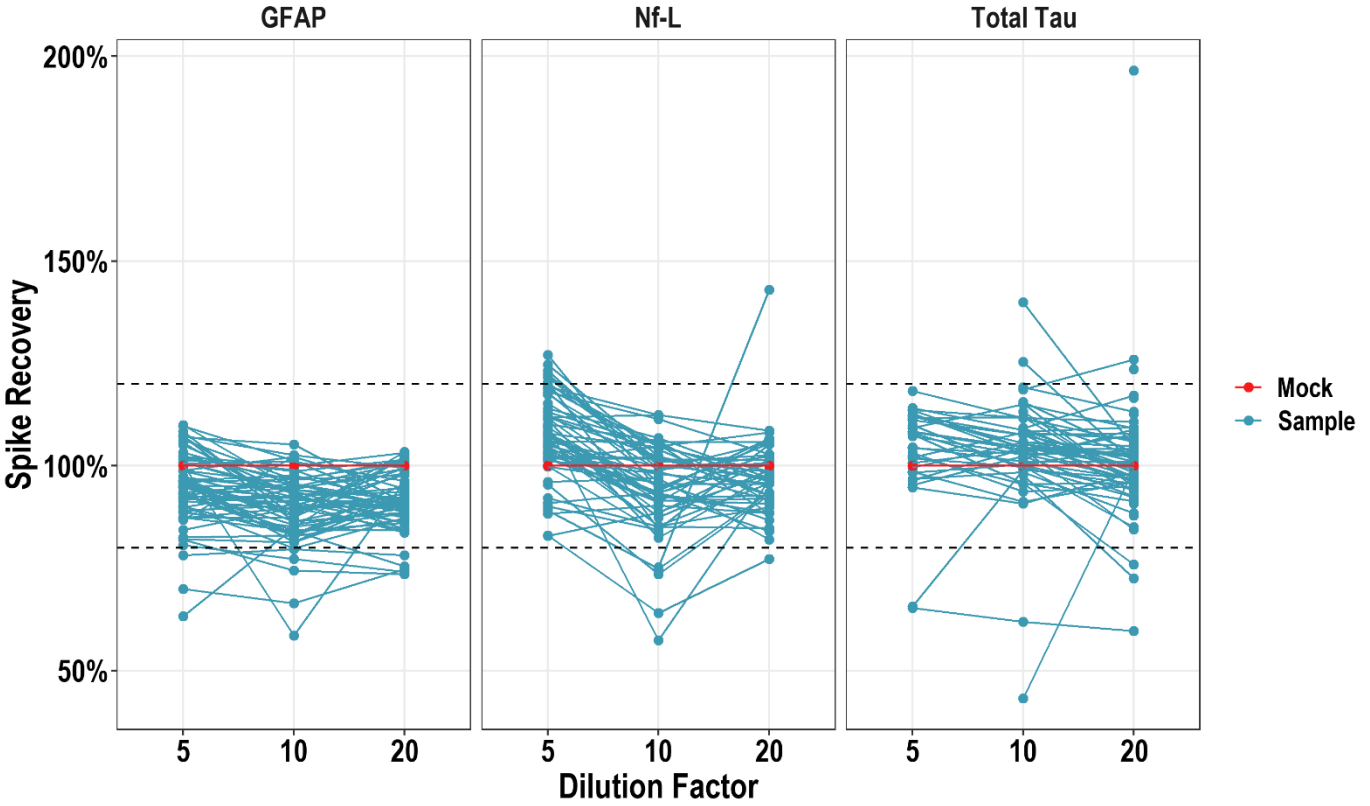


Figure 7 – Spike recovery testing of 64 individual neonatal HIE samples (cord blood) showing recovery of GFAP, NF-L, and tTau in most samples remains within ±20% (dashed lines) at 5-, 10-, and 20-fold dilution.

G. Cross-Reactivity

Cross-reactivity between GFAP, NF-L, and tTau was calculated for both calibrator and native samples by using individual calibrators and individual detection antibodies. Some cross-reactivity of the GFAP calibrator with the NF-L (3.5%) and tTau (2.7%) assays was observed, presumably due to non-specific binding of the recombinant GFAP calibrator, shown in Table 5. This cross-reactivity was not observed for native samples spiked at the CAL 01 level or at 5-times the CAL 01 level. For all target and assay combinations, cross-reactivity was <0.1% in native matrices and should not affect sample quantitation, as shown in the representative data in Table 6.

Table 5 – Summary cross-reactivity data for individual calibrators and detection antibodies.

Calibrator & Detect		GFAP	NF-L	tTau
Spot	Sample Used	Cross-reactivity (%)		
GFAP	Cal 01	—	0.51	0.00
	Cal 02	—	0.53	0.00
NF-L	Cal 01	3.50	—	0.01
	Cal 02	2.52	—	0.00
tTau	Cal 01	2.74	0.24	—
	Cal 02	2.07	0.28	—

Table 6 – Representative cross-reactivity data in native samples.

Plasma 1 (5x CAL01 Spike)		Cross reactivity for GFAP	Cross reactivity for NF-L	Cross reactivity for tTau
	GFAP	—	0.029	0.001
	NF-L	0.043	—	0.004
Plasma 2 (5x CAL01 Spike)	Tau	0.03	0.029	—
		Cross reactivity for GFAP	Cross reactivity for NF-L	Cross reactivity for tTau
	GFAP	—	0.041	0.002
Plasma 3 (CAL01 Spike)	NF-L	0.029	—	0.004
	Tau	0.018	0.03	—
		Cross reactivity for GFAP	Cross reactivity for NF-L	Cross reactivity for tTau
	GFAP	—	0.025	0.001
Plasma 4 (CAL01 Spike)	NF-L	0.014	—	0.003
	Tau	0.008	0.018	—
		Cross reactivity for GFAP	Cross reactivity for NF-L	Cross reactivity for tTau
	GFAP	—	0.031	0.001
	NF-L	0.014	—	0.003
	Tau	0.01	0.025	—

4 Conclusion

MSD has developed and analytically validated a highly sensitive, multiplexed immunoassay for quantifying GFAP, NF-L, and tTau in plasma and serum. This assay is sufficiently sensitive to detect these analytes in normal serum and plasma and has a large dynamic range to accommodate elevated levels found in some neurological disorders, making it suitable for associated research.

5 Funding

Research reported in this publication was supported by the National Institute Of Neurological Disorders And Stroke of the National Institutes of Health under Award Number U01NS114144. The content is solely the responsibility of the authors and does not necessarily represent the official views of the National Institutes of Health.

DOWNLOAD POSTER



MSD® products are for research use only. Not for use in diagnostic procedures. MESO SCALE DISCOVERY, Meso Scale Diagnostics, MSD, mesoscale.com, www.mesoscale.com, methodicalmind.com, www.methodicalmind.com, DISCOVERY WORKBENCH, InstrumentLink, MESO, MesoSphere, Methodical Mind, MSD GOLD, MULTI-ARRAY, MULTI-SPOT, QuickPlex, ProductLink, SECTOR, SECTOR HTS, SECTOR PR, SULFO-TAG, TeamLink, TrueSensitivity, TURBO-BOOST, TURBO-TAG, N-PLEX, R-PLEX, S-PLEX, T-PLEX, U-PLEX, V-PLEX, MSD (design), MSD (luminescent design), Methodical Mind (head logo), 96 WELL SMALL-SPOT (design), 96 WELL 1-, 4-, 7-, 9-, & 10-SPOT (designs), 384 WELL 1- & 4-SPOT (designs), N-PLEX (design), R-PLEX (design), S-PLEX (design), T-PLEX (design), U-PLEX (design), V-PLEX (design), It's All About U, Spot the Difference, The Biomarker Company, and The Methodical Mind Experience are trademarks and/or service marks owned by or licensed to Meso Scale Diagnostics, LLC. All other trademarks and service marks are the property of their respective owners. ©2023 Meso Scale Diagnostics, LLC. All rights reserved.