

# Sensitive Serum/Plasma Neurofilament Light Immunoassay

Martin Stengelin, Pradeepthi Bathala, and Jacob N. Wohlstadter  
Meso Scale Discovery, Rockville, Maryland, USA

## 1 Abstract

Increased concentration of neurofilament light protein (NF-L) in CSF is a proven marker for neuronal injury. Obtaining a CSF sample is invasive; therefore measurement of this biomarker in blood would be preferred. However, use of blood is complicated by the extremely low NF-L concentrations and the presence of many potential assay interferents.

Using a next generation assay format based on MULTI-ARRAY® technology, MSD developed and characterized a highly sensitive immunoassay for NF-L in serum and plasma. Limit of blank (LOB) was determined by running 20 replicates of zero calibrator per plate and calculating the concentration corresponding to the average zero calibrator signal plus 2.5 standard deviations. Lower and upper limits of quantitation (LLOQ and ULOQ, respectively) were determined by running six plates on two days with four replicates per plate. LLOQ was defined as the lowest concentration with a total CV of 20% or less. The reference range in the normal population was established per CLSI EP28-A3c by testing serum samples from 80 apparently healthy individuals. Association of serum/plasma NF-L with neurocognitive status was evaluated by testing 120 samples from individuals with Alzheimer's, dementia, traumatic brain injury (TBI), or HIV with well-characterized neurocognitive status.

The LOB, LLOQ, and ULOQ for the assay were 0.7 pg/mL, 3.4 pg/mL, and 1,700 pg/mL, respectively. Total CVs of four QC samples (15 plates, 5 days) were 8%–14%. Serum and plasma samples diluted linearly. NF-L concentrations in the reference (healthy) sample set (5<sup>th</sup>, 25<sup>th</sup>, 50<sup>th</sup>, 75<sup>th</sup>, and 95<sup>th</sup> percentiles) were 10 pg/mL, 15 pg/mL, 19 pg/mL, 26 pg/mL, and 51 pg/mL, respectively. Median concentrations for mild and severe HIV-Associated Neurocognitive Disorder samples were 35 pg/mL and 56 pg/mL. Median concentration in Alzheimer's, dementia, and TBI samples were 220 pg/mL, 154 pg/mL, and 135 pg/mL, respectively. Matched serum and plasma concentrations correlated well ( $R^2 = 0.73$ ). NF-L concentrations in the normal population correlated with age ( $p=0.03$ ).

The assay enables quantitation of low NF-L levels in serum and plasma samples and should prove useful for studying neuronal injury caused by diseases such as Alzheimer's or HIV.

## 2 Methods

MSD's electrochemiluminescence (ECL) detection technology uses SULFO-TAG™ labels that emit light upon electrochemical stimulation initiated at the electrode surfaces of MULTI-ARRAY and MULTI-SPOT® microplates.

A next-generation MULTI-ARRAY technology with significantly higher sensitivity was used for the data presented in this poster.

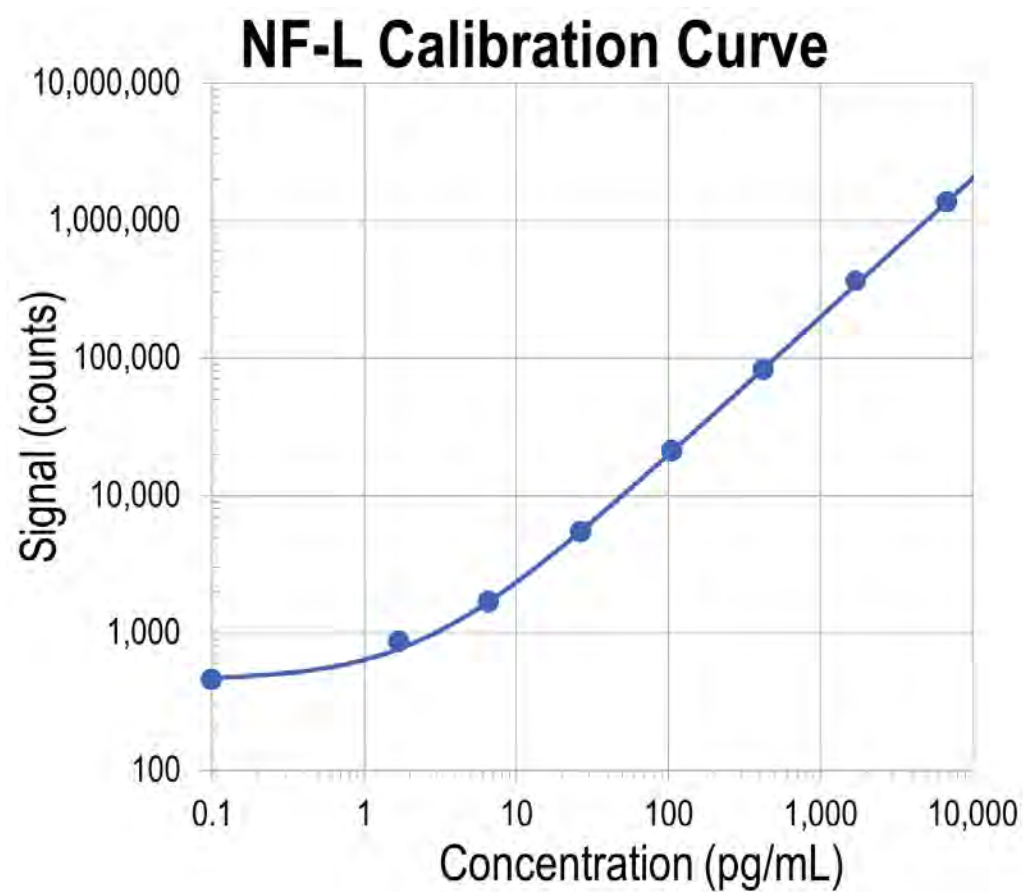


### Electrochemiluminescence Technology

- 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 Calibration Curve

Native purified NF-L from bovine spinal cord was used to calibrate the assay. The figure to the right shows a typical calibration curve. ECL technology enables a wide dynamic range (four orders of magnitude).



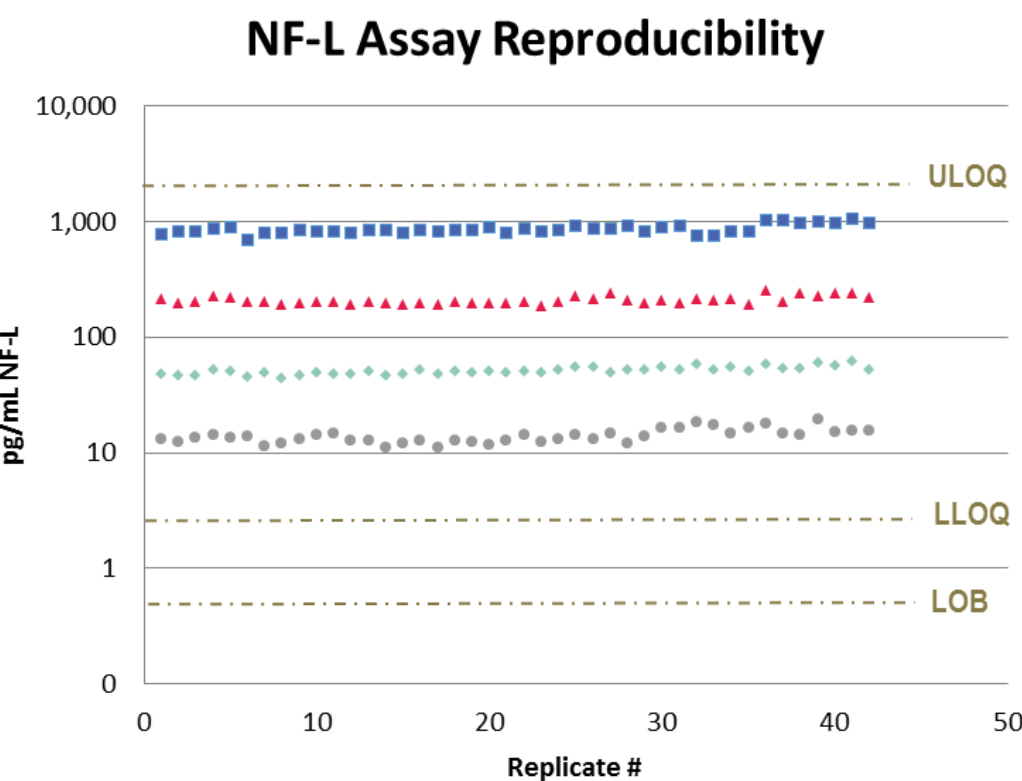
## 4 Assay Range

	Sample	Signal (counts)	pg/mL	Conc. %CV
ULOQ	ULOQ1	436,589	1,700	8
	ULOQ2	215,338	850	9
	ULOQ3	104,305	425	3
LLOQ	LLOQ1	25,689	106	7
	LLOQ2	13,419	53	9
	LLOQ3	7,021	27	7
	LLOQ4	3,804	13	7
	LLOQ5	2,268	7	8
	LLOQ6	1,342	3.4	13
	LLOQ7	962	1.7	29

The LOB was determined by running 20 replicates of zero calibrator per plate and calculating the concentration corresponding to the average zero calibrator signal plus 2.5 standard deviations. The LLOQ and ULOQ were determined by running six plates on two days with four replicates per plate. LLOQ was defined as the lowest concentration with a total CV of 20% or less. ULOQ was defined as the highest concentration with a total CV of 20% or less.

LOB	0.7 pg/mL
LLOQ	3.4 pg/mL
ULOQ	1,700 pg/mL

## 5 Reproducibility



	QC 01	QC 02	QC 03	QC 04
NF-L (pg/mL)	862	207	51	14
CV (21 plates, 7 days)	9%	8%	8%	14%

The reproducibility of the NF-L assay was determined by including two replicates each of four QC samples spanning the assay range in each run. The assay had good reproducibility: Total CVs over 21 plates run on 7 days were between 8 and 14%.

## 6 Clinical Samples

**Reference range samples:** 80 serum samples from apparently healthy individuals were tested to determine the NF-L reference range. Samples reflected a representative adult U.S. population with respect to race, age, and gender.

**CHARTER samples:** 120 well-characterized serum and EDTA plasma samples from the CNS HIV Anti-Retroviral Therapy Effects Research (CHARTER) cohort were used for this study. This sample set included 11 individuals with HIV-Associated Dementia (HAD), 25 individuals with mild neurocognitive disorder (MND), and 30 individuals with no or asymptomatic neurocognitive impairment (ANI).

**Remnant serum, EDTA plasma, and heparin plasma samples** from apparently healthy individuals and from individuals with Alzheimer's, dementia, and TBI were obtained from commercial sample vendors. Only limited information such as age and gender was provided for most of these samples.

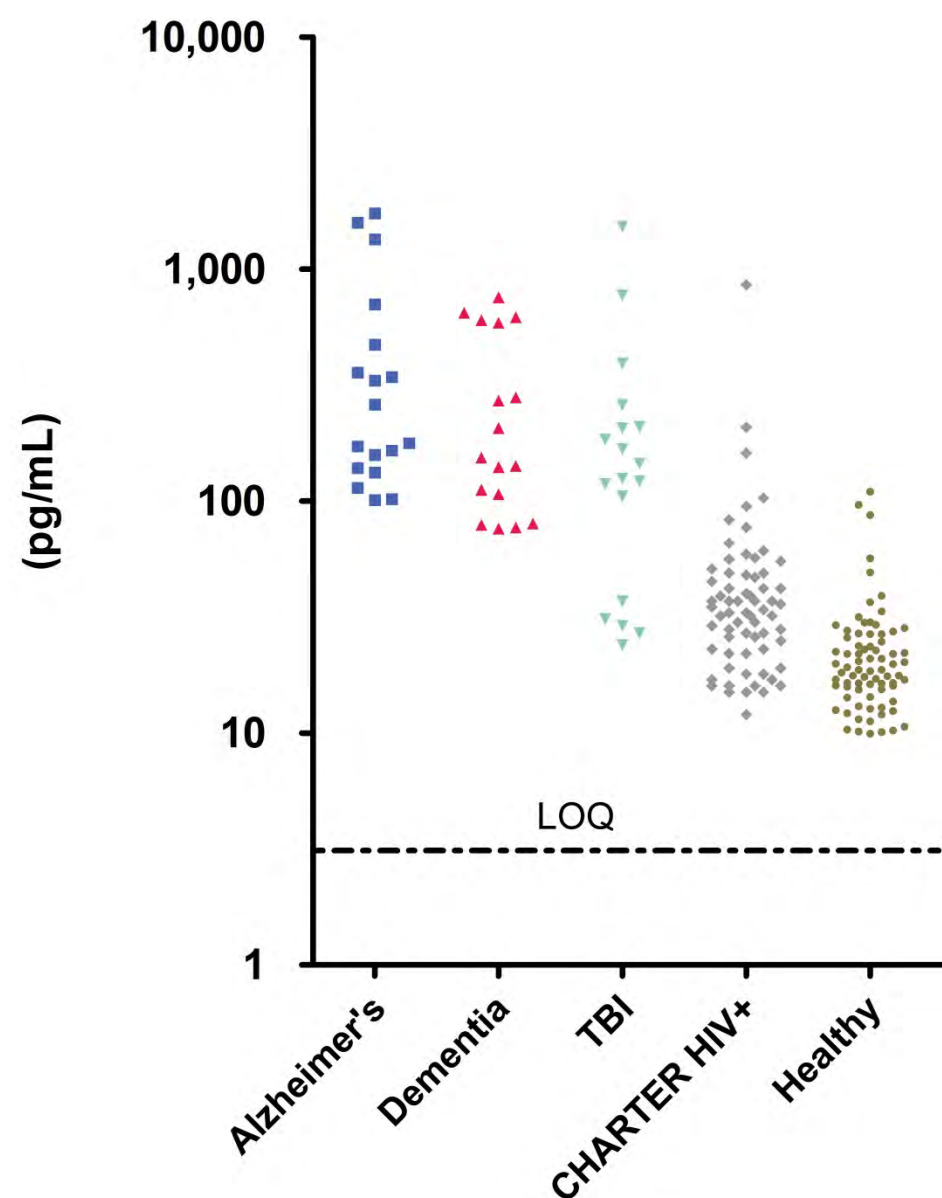
### Acknowledgements

Samples from the CHARTER cohort were used for this study. This publication was made possible from NIH funding through the NIMH and NINDS institutes by the following grants: Manhattan HIV Brain Bank (MHBB) U24MH100931; Texas NeuroAIDS Research Center (TNRC) U24MH100930; National Neurological AIDS Bank (NNAB) U24MH100929; California NeuroAIDS Tissue Network (CNTN) U24MH100928; Data Coordination Center (DCC) U24MH100925.

Research reported in this publication was supported by the National Institute Of Allergy And Infectious Diseases, the National Institute of Aging, and the National Institute of Mental Health, Division of AIDS, of the National Institutes of Health under Award Number U24AI118663.

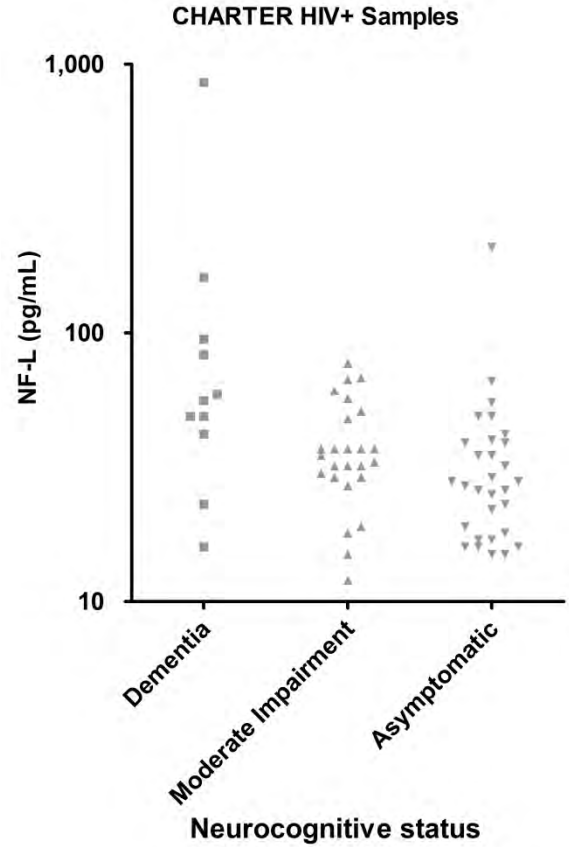
The content is solely the responsibility of the authors and does not necessarily represent the official views of the National Institutes of Health or NNTC.

## 7 NF-L Concentrations in Serum and Plasma Samples



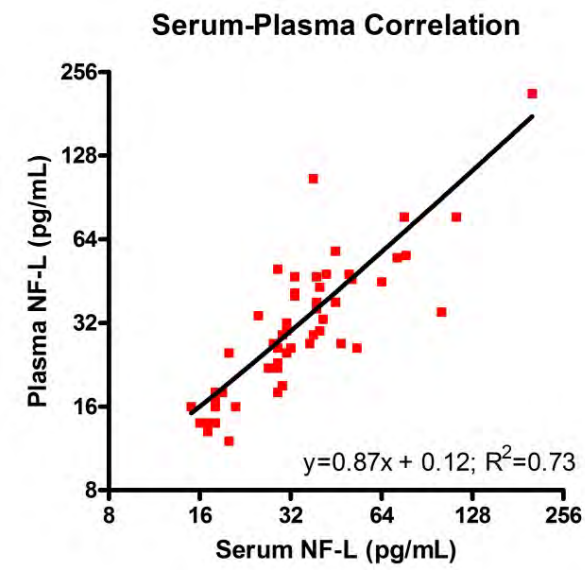
The figure to the right shows dependence of NF-L serum or plasma concentrations on neurocognitive status in samples from the CHARTER cohort.

The figure to the left summarizes testing of commercial samples from individuals with Alzheimer's, dementia, and TBI; from HIV positive individuals from the CHARTER cohort, and from apparently healthy individuals. NF-L concentrations in the reference (healthy) sample set (5<sup>th</sup>, 25<sup>th</sup>, 50<sup>th</sup>, 75<sup>th</sup>, and 95<sup>th</sup> percentiles) were 10 pg/mL, 15 pg/mL, 19 pg/mL, 26 pg/mL, and 51 pg/mL, respectively. Median concentrations for mild and severe HIV-Associated Neurocognitive Disorder samples were 35 pg/mL and 56 pg/mL. Median concentration in Alzheimer's, dementia, and TBI samples were 220 pg/mL, 154 pg/mL, and 135 pg/mL, respectively.



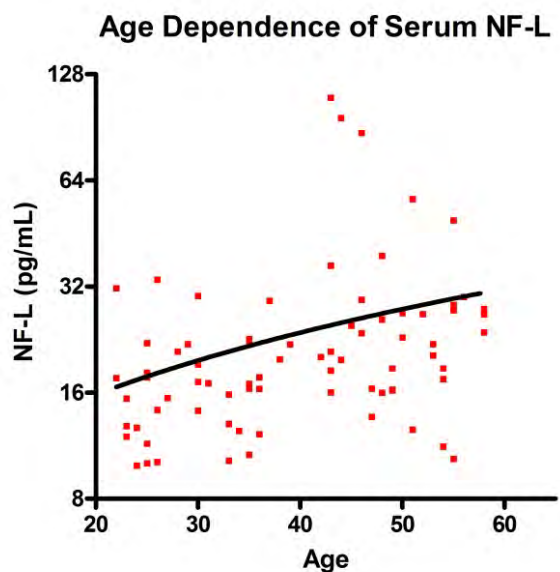
## 8 Serum Plasma Correlation

The CHARTER cohort included matched serum and EDTA plasma samples. Plasma and serum correlated well.



## 9 Dependence on Age and Gender

Serum/plasma NF-L concentrations in apparently healthy individuals did not depend on gender, but increased with age.



## 10 Dilution Linearity; Parallelism

	Dilution Factor	EDTA Plasma				Dilution Factor	Serum		
		Signal (counts)	pg/mL (dilution corrected)	Recovery			Signal (counts)	pg/mL (dilution corrected)	Recovery
Normal Donor 1	neat	4,627	11.3	100%	Normal Donor 9	neat	10,477	27.9	100%
	2	3,106	13.8	122%		2	5,502	27.6	99%
	4	1,844	12.8	114%		4	3,352	30.4	109%
Normal Donor 2	8	1,360	14.2	126%	Normal Donor 10	8	2,037	30.2	108%
	neat	n/a	n/a	n/a		neat	6,442	16.5	100%
	2	7,116	36.8	100%		2	3,775	17.6	107%
Normal Donor 3	4	4,249	40.7	111%	Normal Donor 11	4	2,374	19	116%
	8	2,525	41.6	113%		8	1,479	17	103%
	neat	6,314	16.3	100%		neat	7,569	19.9	100%
Normal Donor 4	2	3,096	14.2	87%	Normal Donor 12	2	4,083	19.9	100%
	4	1,917	14.8	91%		4	2,622	22.9	115%
	8	1,336	16.3	100%		8	1,580	21.9	110%
TBI Donor 5	neat	3,846	9.25	100%	TBI Donor 13	neat	24,565	68.7	100%
	2	2,311	9.68	105%		2	14,172	77.8	113%
	4	1,388	8.75	95%		4	7,375	77.5	113%
TBI Donor 6	8	999	8.56	93%	TBI Donor 14	8	4,286	84.1	122%
	neat	21,081	57.6	100%		neat	24,707	67.7	100%
	2	10,941	58.4	101%		2	13,946	75.3	111%
TBI Donor 7	4	5,934	60.1	104%	TBI Donor 15	4	7,062	72.9	108%
	8	3,398	61.9	107%		8	4,363	84.1	124%
	neat	22,263	60.9	100%		neat	63,231	174	100%
TBI Donor 8	2	12,579	67.7	111%	TBI Donor 16	2	30,801	169	97%
	4	6,911	71.2	117%		4	15,686	170	98%
	8	3,562	65.6	108%		8	8,032	168	97%
TBI Donor 9	neat	46,606	132	100%	TBI Donor 17	neat	10,255	27.6	100%
	2	24,679	138	105%		2	5,613	28.6	104%
	4	14,752	162	123%		4	3,046	27.8	101%
TBI Donor 10	8	7,563	159	121%	TBI Donor 18	8	1,948	30.4	110%
	neat	38,141	108	100%		neat	67,595	192	100%
	2	19,826	110	102%		2	35,272	199	103%
TBI Donor 11	4	10,539	114	106%	TBI Donor 19	4	16,925	187	97%
	8	5,510	112	104%		8	9,096	195	101%

## 11 Spike Recovery

	Spike Conc. (pg/mL)	EDTA Plasma				Spike Conc. (pg/mL)	Heparin Plasma				Spike Conc. (pg/mL)	Serum		
		Signal (counts)	Conc. (pg/mL)	Recovery (%)			Signal (counts)	Conc. (pg/mL)	Recovery (%)			Signal (counts)	Conc. (pg/mL)	Recovery (%)
Donor 1	5,000	149,113	5,525	109	Donor 4	5,000	91,586	3,466	68	Donor 7	5,000	106,977	4,022	79
	2,500	61,072	2,350	91		2,500	36,722	1,441	55		2,500	47,451	1,845	71
	1,250	33,533	1,320	99		1,250	22,359	892	66		1,250	28,779	1,139	85
Donor 2	0	2,208	81	Ave. = 100	Donor 5	0	1,852	65	Ave. = 63	Donor 8	0	1,986	71	Ave. = 78
	5,000	129,315	4,822	96		5,000	85,490	3,245	64		5,000	123,716	4,621	92
	2,500	55,282	2,137	84		2,500	39,001	1,528	60		2,500	51,649	2,001	78
Donor 3	1,250	34,122	1,343	104	Donor 6	1,250	22,113	882	68	Donor 9	1,250	27,029	1,072	82
	0	1,278	39	Ave. = 95		0	1,054	29	Ave. = 64		0	1,403	45	Ave. = 84
	5,000	123,001	4,596	91		5,000	97,547	3,682	73		5,000	103,581	3,899	77
Control Spike (Calibr. Diluent)	2,500	49,047	1,904	75	Control Spike (Calibr. Diluent)	2,500	43,238	1,687	66	Control Spike (Calibr. Diluent)	2,500	50,438	1,956	77
	1,250	26,876	1,066	82		1,250	25,066	996	76		1,250	26,491	1,051	82
	0	1,263	39	Ave. = 83		0	1,332	42	Ave. = 72		0	1,090	31	Ave. = 79

## 12 Conclusion

Using a next generation assay format based on MULTI-ARRAY technology, MSD developed and characterized a highly sensitive immunoassay for NF-L in serum and plasma. The sensitivity of this assay is an order of magnitude better than typical concentrations in normal human serum or plasma. The assay should prove useful for studying neuronal injury caused by diseases such as Alzheimer's or HIV.

MSD products are For Research Use Only. Not for use in diagnostic procedures.

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, MESO, [www.mesoscale.com](http://www.mesoscale.com), DISCOVERY WORKBENCH, MSD GOLD, MULTI-ARRAY, MULTI-SPOT, QUICKPLEX, SECTOR, SECTOR PR, SECTOR HTS, STREPTAVIDIN GOLD, SULFO-TAG, TrueSensitivity, TURBO-BOOST, TURBO-TAG, R-PLEX, S-PLEX, T-PLEX, U-PLEX, V-PLEX, MSD (design), 96 WELL SMALL SPOT (design), 96 WELL 1-, 4-, 7-, 9-, & 10-SPOT (designs), 384 WELL 1- & 4-SPOT (designs), R-PLEX (design), S-PLEX (design), T-PLEX (design), U-PLEX (design), V-PLEX (design), It's All About U, and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC. ©2019 Meso Scale Diagnostics, LLC. All rights reserved.

DOWNLOAD POSTER

