



S-PLEX[®] Neurology Panel 1 Kits

For use with cerebral spinal fluid (CSF), serum, plasma, and cell culture supernatants

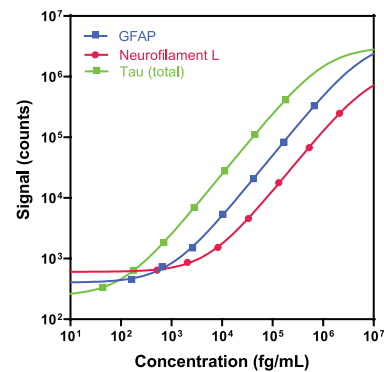
S-PLEX[®] from MESO SCALE DISCOVERY[®] TrueSensitivity[®], Simple Execution

- Ultrasensitive singleplex and multiplex kits
- Superior sample matrix performance
- Require minimal sample volume
- Run on existing MSD instruments
- Reliable performance and quality

Sensitivity you can trust

S-PLEX Neurology Panel 1				
Analyte	Median LLOD (fg/mL)	LLOD range (fg/mL)	LLOQ (fg/mL)	ULOQ (fg/mL)
GFAP	520	290–600	1,340	450,000
Neurofilament L	1,700	1,400–2,100	8,480	1,400,000
Tau (total)	75	62–120	180	120,000

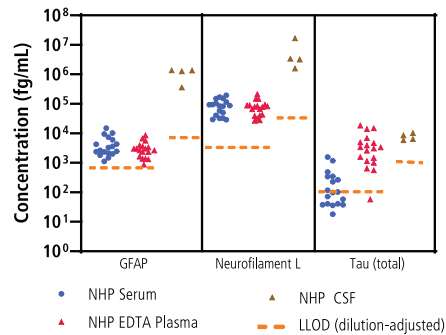
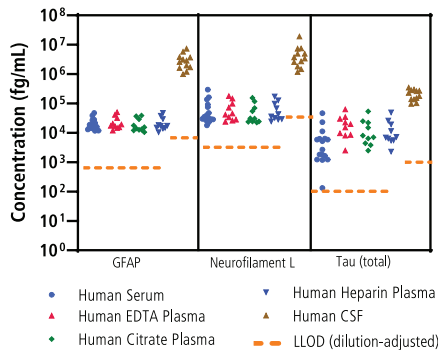
Wide dynamic range



Numerous studies have confirmed GFAP, Neurofilament L, and Tau as critical biomarkers in CSF and blood-based samples for neurological research. Ultrasensitive immunoassays are required to reliably measure these biomarkers in blood samples because of their low concentrations. MSD has developed an ultrasensitive multiplex panel using the S-PLEX platform under rigorous design control and according to "Fit for Purpose" principles¹ to provide the analytical performance required for reliable measurement of these biomarkers in human and NHP samples.

Native Level Detection

The S-PLEX Neurology Panel 1 enables researchers to measure neurological biomarkers in serum, plasma, and CSF from humans and NHPs.

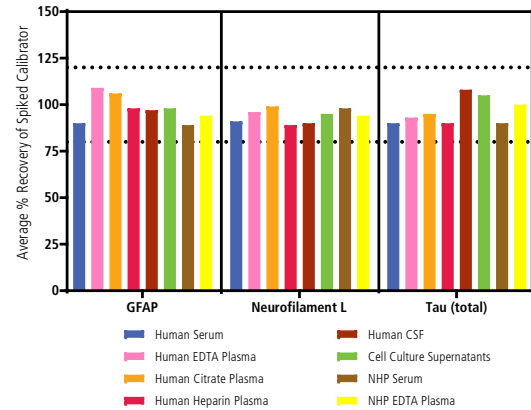
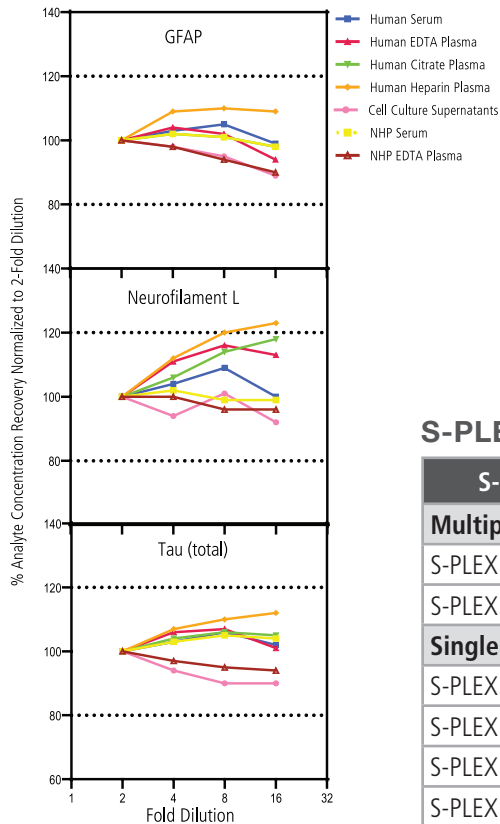


1. Lee JW, et al. Fit-for-purpose method development and validation for successful biomarker measurement. Pharm Res. 2006; 23:3 12-28.

For Research Use Only. Not for use in diagnostic procedures.

Overcoming Matrix Challenges and Recovery

Interference from the sample matrix presents a challenge when measuring neurological biomarkers. The dilution linearity and spike recovery profiles shown below exemplify the excellent matrix tolerance of assays in the S-PLEX Neurology Panel 1.



S-PLEX Neurology Products

S-PLEX Neurology	Assays	Catalog Numbers
Multiplex and Custom Panels		
S-PLEX Neurology Panel 1 Kits	GFAP, Neurofilament L, Tau (total)	K15639S (Hu) and K15640S (NHP)
S-PLEX Custom Neurology	Customized panel	K151ANV (Hu) and K156ANV (NHP)
Singleplex Kits		
S-PLEX GFAP Kits	GFAP	K151AJXS (Hu) and K156AJXS (NHP)
S-PLEX Neurofilament L Kits	Neurofilament L	K151AKGS (Hu) and K156AKGS (NHP)
S-PLEX Tau (total) Kits	Tau (total)	K151AGPS (Hu) and K156AGPS (NHP)
S-PLEX Tau (pT181) Kits	Tau (pT181)	K151AGMS (Hu) and K156AGMS (NHP)
S-PLEX Tau (pT231) Kits	Tau (pT231)	K151AGNS (Hu) and K156AGNS (NHP)

Related MSD Neurology Products

In addition to the S-PLEX platform, MSD provides solutions to support and simplify your neuroinflammation and neurodegeneration research needs. We offer singleplex and multiplex assays to measure biomarkers such as A β 38, A β 40, A β 42 peptides and α -Synuclein under various MSD's platforms. All assays ship in 1 to 3 days.

- V-PLEX[®] and V-PLEX Plus A β Peptide Panel 1 (6E10) Kit
- V-PLEX Neuroinflammatory Panel 1 (human) Kit
- V-PLEX and V-PLEX Plus A β Peptide Panel 1 (4G8) Kit
- U-PLEX[®] and U-PLEX Plus Human α -Synuclein Kit

We are continually expanding our portfolio by developing new and improved assays for neurology-focused biomarkers across various assay platforms. For the most-up-to-date information and a complete listing of our products, visit www.mesoscale.com/en/applications and click "Neurobiology/Neurodegeneration."

Phone: 1-240-314-2795 . Fax: 1-301-990-2776
 Email: customerservice@mesoscale.com

For Research Use Only. Not for use in diagnostic procedures.

MESO SCALE DISCOVERY, MSD, www.mesoscale.com, TrueSensitivity, S-PLEX, U-PLEX, V-PLEX, MSD (design), 96 WELL 10-SPOT (design), S-PLEX (design), and Spot the Difference are trademarks and/or service marks owned by 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.

VIEW MORE ONLINE



MK-DS-1166-v1-Feb2023

