



Available Now!

Validated Analytical Testing Service for Ultrasensitive Tau (pT217) and (pT181) Measurements

MSD is pleased to introduce a validated analytical testing service for measuring phosphorylated Tau at threonine 217 (pT217) and threonine 181 (pT181) in human samples through our Bioanalytical Lab (BAL). The pT217 and pT181 singleplex S-PLEX[®] assays achieve fg/mL levels of detection and enable the interrogation of these biomarkers in both normal and diseased samples.

MSD's Bioanalytical Lab provides comprehensive services to analyze your samples. The pT217 and pT181 assays have undergone rigorous analytical method validation, allowing MSD to offer RUO, GLP, and GCLP sample testing.

Accelerate your neurology biomarker research program using our new testing service for pT217 and pT181 biomarkers.

Key Benefits

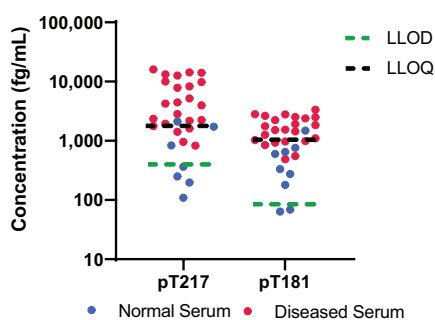
- ✓ Assays developed using MSD's ultrasensitive S-PLEX platform
- ✓ Assays validated for serum, EDTA plasma, and CSF
- ✓ Assays optimized for accuracy, precision, sensitivity, specificity, linearity, and recovery

TrueSensitivity[®] You can Trust

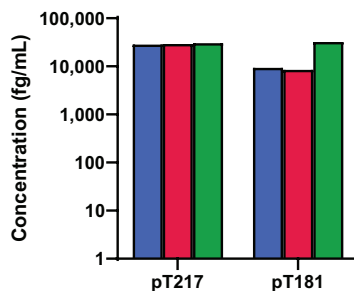
Analyte	LLOD (fg/mL)	LLOQ (fg/mL)	ULOQ (fg/mL)
pT217	395	1,836	3,760,000
pT181	84.7	1,079	2,210,000

Tau (pT217) and (pT181): Key Biomarkers in Neurological Disorders

Recent studies have shown that pT217 and pT181 are relevant biomarkers for the study of various neurological disorders, including Alzheimer's disease (AD). Meaningful measurements of pT217 and pT181 in blood matrices in addition to traditional CSF samples require improved assay sensitivity. MSD's new assays address this unmet need.



Representative BAL data above show the biomarker profile between normal serum (•N=7) and diseased serum from AD patients (•N=23). Dashes show the limit of detection (--), and lower limit of quantification (--).

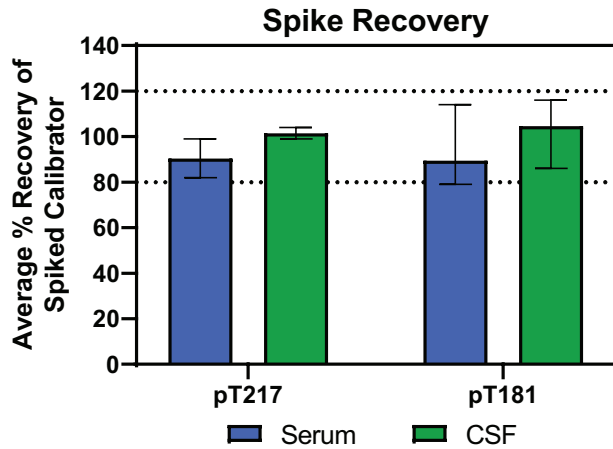


- Plasma spiked with Assay Specific Calibrator
- Plasma spiked with Assay Specific Calibrator+non-Phosphorylated Tau
- Plasma spiked with Assay Specific Calibrator+Mutated pT217

MSD's pT217 and pT181 assays are specific to the target molecule. Assay specificity was evaluated by MSD's BAL in plasma samples spiked with assay-specific calibrators in the presence of either non-phosphorylated Tau (400 pg/mL) or mutated phosphorylated Tau217 proteins (400 pg/mL).

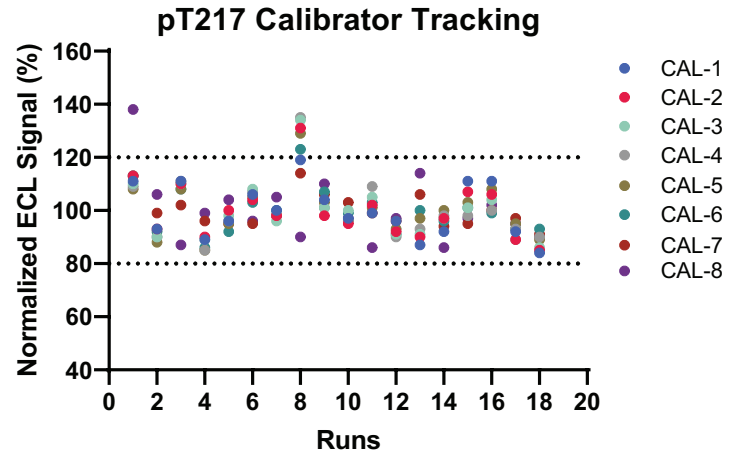
Analytical Method Validation in Multiple Sample Matrices

Matrix studies such as dilution linearity and spike recovery measurements in different sample types are part of the rigorous analytical method validation process in MSD's Bioanalytical Lab. Representative percent recovery data for pT217 and pT181 assays are shown below as an example of the high level of accuracy observed in key validated sample matrices.



Superior Performance Between Runs

During validation, calibrator and quality control sample reproducibilities were tested. The graph below shows the normalized ECL signal tracking for each calibrator level across n=18 runs, which varied by less than $\pm 20\%$ in most runs. The inter-run calibrator and control signal %CVs were $< 15\%$ (data not shown).



Contact us to discuss your sample testing service needs.

Step 1

Contact MSD or Account Manager to receive a quote, or request a quote online.

Step 2

Place your order referencing your quote number.

Step 3

Ship samples to MSD with appropriate documentation.

Step 4

Receive a detailed report with analytical results.

Phone: 1-240-314-2793 . Fax: 1-301-990-2776

Email: bioanalytical.services@mesoscale.com
customerservice@mesoscale.com

Meso Scale Diagnostics, MSD, mesoscale.com, www.mesoscale.com, TrueSensitivity, S-PLEX, MSD (design), 96 WELL 10-SPOT (design), and Spot the Difference are trademarks and/or service marks owned by Meso Scale Diagnostics, LLC.

©2023 Meso Scale Diagnostics, LLC. All rights reserved.



VIEW MORE ONLINE



MK-DS-1171-v1-2023Jul

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



www.mesoscale.com