

**CERTIFICATE OF ANALYSIS**


Product Description: S-PLEX Neurology Panel 1 (human) Calibrator Blend
Catalog Number: C01639-2
Lot Number: A00S0133
Expiration Date: 31 OCT 2027
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
GFAP	18120000	fg/mL
Neurofilament L	91400000	fg/mL
Tau (total)	3220000	fg/mL

Dilution to Top of Curve: 20 X

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		11 JUN 2025