

Product Number:	103345
Lot Number:	503762
Expiration:	28-Feb-2024
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	103338	305906	N/A
Detector Reagent	103339	306112	N/A
SBG Reagent	103340	306113	N/A
Sample Diluent	103341	306114	N/A
Calibrators	103342	314607	314607A
Control 1	103343	310209	310209A
Control 2	103344	310209	310209A
SBG Diluent (SR-X)†	103158	N/A ¹	N/A ¹
RGP Reagent	103159	N/A ¹	N/A ¹

¹ RGP is not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa™ Neuro 4-Plex B SR-X Kit Instructions (KI-0123).



Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials

Tau

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	310209	2.58	1.78 - 2.68
Control 2	310209	86.4	33.1 - 110.5
Panel 2	306108	2.64	2.00 - 3.30
Panel 3	306108	26.6	21.1 - 31.6
Panel 4	306108	97.3	78.1 - 117

NF-light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	310209	4.20	2.88 - 5.21
Control 2	310209	308	211 - 398
Panel 2	306108	11.8	9.90 - 14.9
Panel 3	306108	72.7	63.0 - 95.0
Panel 4	306108	355	304 - 456

GFAP®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	310209	154	100 - 167
Control 2	310209	12981	6303 - 16032
Panel 2	306108	85.0	64.5 - 96.8
Panel 3	306108	1079	754 - 1173
Panel 4	306108	6738	4626 - 7487

UCH-L1®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	310209	136	101 - 151
Control 2	310209	10130	7300 - 12748
Panel 2	306108	122	86.3 - 151
Panel 3	306108	1731	1557 - 2335
Panel 4	306108	12610	10263 - 15395

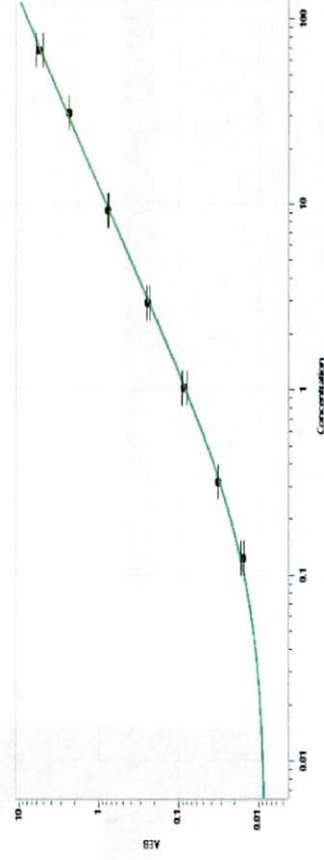
Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Review/Approval

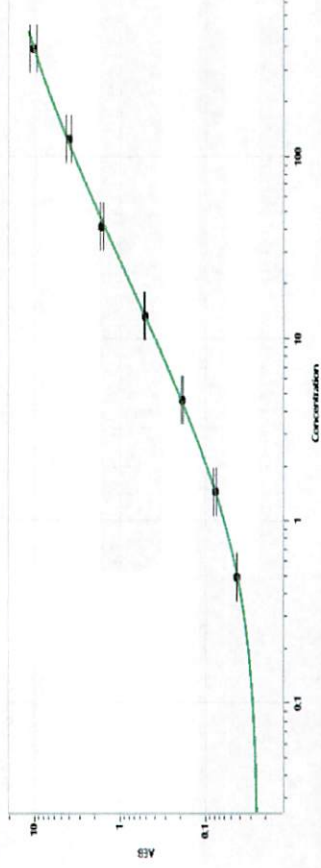
Sonia Khillan oc
Name Title

Sonia Kella 16 JUN 2023
Signature/Date

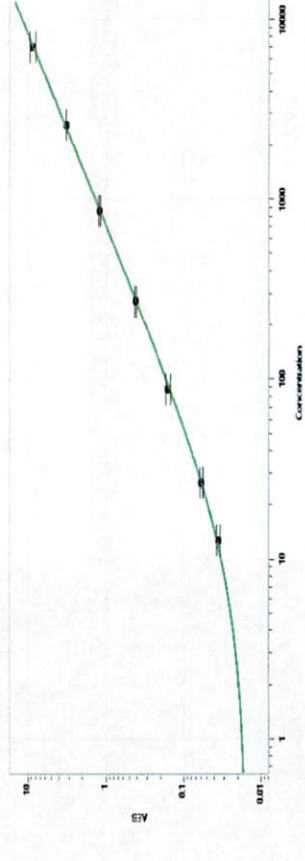
Tau Calibration Curve



NF-light® Calibration Curve



GFAP® Calibration Curve



UCHL1® Calibration Curve

