

Product Number:	103345
Lot Number:	502652
Expiration:	20-Nov-2021
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103338	034307
Detector Reagent	103339	034408
SBG Reagent	103340	034412
Sample Diluent	103341	034410
Calibrators	103342	033705
Control 1	103343	034501
Control 2	103344	034501
SBG Diluent (SR-X) †	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> 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)	Mean Range (pg/mL)
Control 1	034501	1.75	1.46-2.27
Control 2	034501	73.9	64.2-96.2
Panel 2	921102	1.23	0.312 - 2.93
Panel 3	921103	27.2	20.3 - 34.0
Panel 4	921104	43.0	29.9 - 61.1

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	034501	7.44	6.18-9.51
Control 2	034501	646	541-837
Panel 2	921102	6.24	5.75 - 8.63
Panel 3	921103	168	147 - 221
Panel 4	921104	575	555 - 832

GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	034501	144	109-184
Control 2	034501	9563	7644-12407
Panel 2	921102	31.5	27.8 - 53.1
Panel 3	921103	847	638 - 1195
Panel 4	921104	5026	3584 - 6789

UCH-L1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	034501	134	112-168
Control 2	034501	10585	9114-13671
Panel 2	921102	104	84.5 - 127
Panel 3	921103	3115	2398 - 4156
Panel 4	921104	7881	6436 - 10728

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

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 28 Dec 2020  
Signature/Date