

Product Number:	103520
Lot Number:	503170
Expiration:	4-Oct-2022
Platform(s):	HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103510	127706
Detector Reagent	103512	127802
SBG Reagent	103514	127703
Sample Diluent	103516	127704
Calibrators	103517	134901
Control 1	103518	134902
Control 2	103519	134902
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 2-Plex B SR-X Kit Instructions (KI-0062).



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

**Release Materials**

**NF-light®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134902	5.15	3.22 - 5.89
Control 2	134902	440	339 - 509
Panel 1	128902	14.9	7.33 - 16.3
Panel 2	128902	14.3	8.91 - 15.3
Panel 3	128902	138	101 - 152
Panel 4	128902	376	261 - 436

**GFAP®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134902	111	80.0 - 143
Control 2	134902	12576	9784 - 15179
Panel 1	128902	87.9	44.4 - 101
Panel 2	128902	67.1	45.0 - 79.6
Panel 3	128902	5093	3319 - 6109
Panel 4	128902	23252	17135 - 25703

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

Review/Approval

Shivani Goel  
Name

QC

Title

Shivani Goel

12 Jun 2022

Signature/Date



