

Product Number:	103670
Lot Number:	502721
Expiration:	12-Jan-2022
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103653	101208
Detector Reagent	103655	105415
SBG Reagent	103657	103313
Plasma Sample Diluent	103659	103312
CSF Sample Diluent	103727	103314
Calibrators	103667	103303
Control 1	103668	103304
Control 2	103669	103304
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials			
Abeta 40			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	103304	15.5	12.2-18.3
Control 2	103304	111	85.5-128
Panel 1	027301	32.7	23.3-35.0
Panel 2	018109	53.5	42.3-64.6
Panel 3	018109	134	100-163
Abeta 42			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	103304	5.11	3.97-5.96
Control 2	103304	28.2	21.1-31.7
Panel 1	027301	2.84	2.01-3.03
Panel 2	018109	22.9	16.6-24.8
Panel 3	018109	45.9	36.0-54.0
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	103304	217	161-242
Control 2	103304	4115	3085-4967
Panel 1	027301	52.6	37.8-56.7
Panel 2	018109	605	446-668
Panel 3	018109	2494	1916-2948
NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	103304	27.4	21.3-31.9
Control 2	103304	552	446-670
Panel 1	027301	10.1	8.37-12.6
Panel 2	018109	75.9	59.5-89.3
Panel 3	018109	399	310-474

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

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 31 Mar 2022
	Name	Title	Signature/Date