

Product Number:	102153
Lot Number:	503265
Expiration:	28-Oct-2022
Platform(s):	HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	102141	130201
Detector Reagent	102143	131908
SBG Reagent	102145	130107
Sample Diluent	102147	130108
Calibrators	102150	134311
Control 1	102151	134312
Control 2	102152	134312
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹



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

¹ SBG Diluent (SR-X) and RGP reagent are not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa™ Neuro 4-Plex A SR-X Kit Instructions (KI-0031).

Release Materials

Tau

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134312	2.40	1.47 - 3.08
Control 2	134312	84.6	71.9 - 108
Panel 1	111709	4.97	1.11 - 7.56
Panel 2	111709	4.62	1.34 - 7.79
Panel 3	111709	25.4	22.5 - 33.7

NF-light®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134312	4.92	2.69 - 7.55
Control 2	134312	462	370 - 555
Panel 1	111709	4.82	4.12 - 6.23
Panel 2	111709	9.05	7.75 - 11.6
Panel 3	111709	162	147 - 232

GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134312	27.8	17.5 - 39.5
Control 2	134312	1097	878 - 1317
Panel 1	111709	36.4	28.0 - 42.1
Panel 2	111709	104	85.8 - 129
Panel 3	111709	1537	1331 - 1996

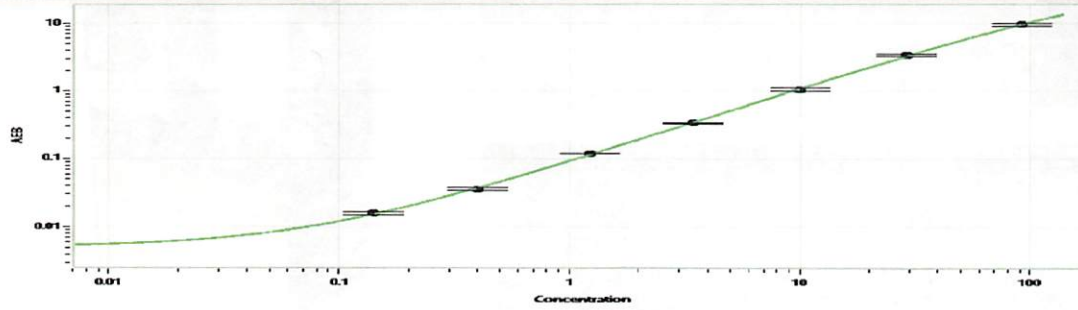
UCHL1®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134312	114	75.2 - 165
Control 2	134312	9330	7521 - 11281
Panel 1	111709	96.2	69.9 - 123
Panel 2	111709	116	83.9 - 136
Panel 3	111709	1594	1450 - 2176

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	09Mar2022
	Name	Title	Signature/Date	

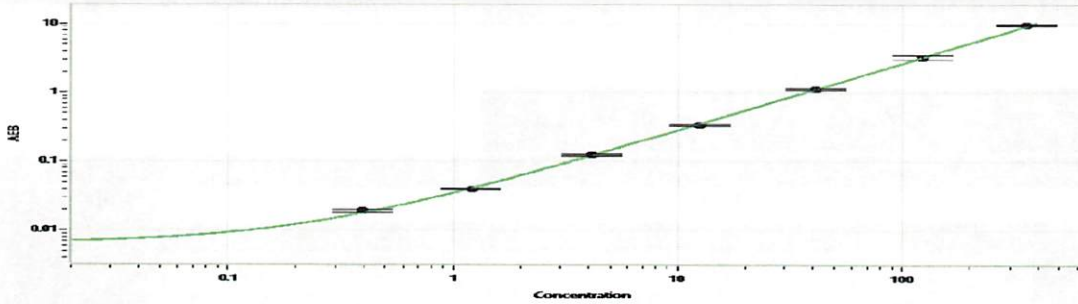
Tau Calibration Curve



Calibrator Levels
(pg/mL)

A	0.000
B	0.140
C	0.400
D	1.23
E	3.44
F	10.0
G	29.0
H	92.1

NF-light® Calibration Curve

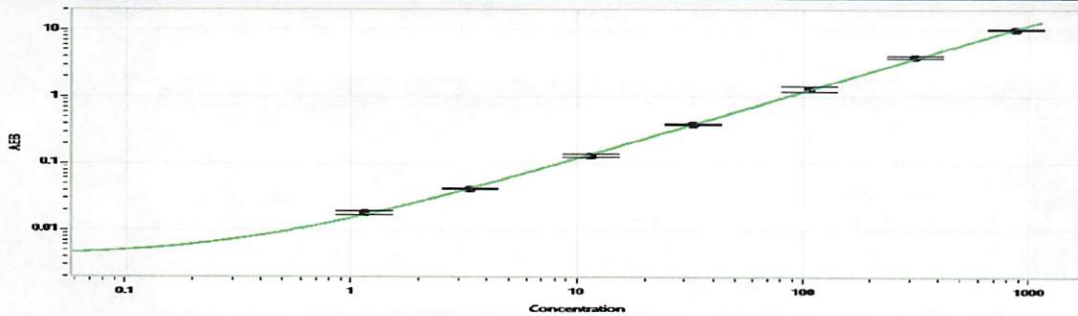


Calibrator Levels
(pg/mL)

A	0.000
B	0.396
C	1.21
D	4.15
E	12.6
F	41.2
G	122
H	356

84.6

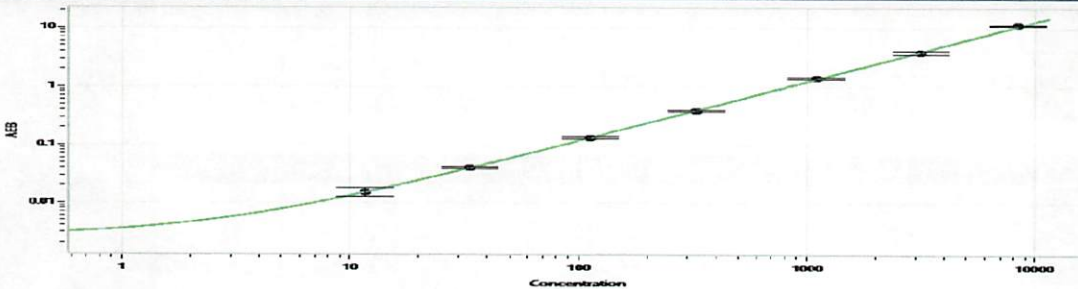
GFAP® Calibration Curve



Calibrator Levels
(pg/mL)

A	0.000
B	1.15
C	3.36
D	11.5
E	32.8
F	108
G	314
H	872

UCHL1® Calibration Curve



Calibrator Levels
(pg/mL)

A	0.000
B	11.6
C	33.3
D	112
E	326
F	1098
G	3136
H	8431

116.2945204

1594.346302

Shivani Goel

QC manager