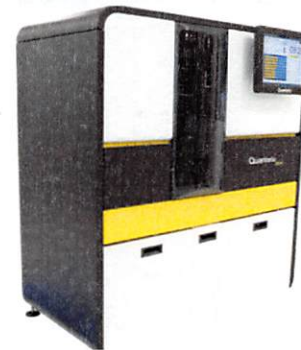


Product Number:	102153
Lot Number:	502023
Expiration:	24-Jun-2020
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	102141	925505
Detector Reagent	102143	925506
SBG Reagent	102145	917502
Sample Diluent	102147	924601
Calibrators	102150	920401
Controls	102151-102152	920402
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	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 A SR-X Kit Instructions (KI-0031).



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

Release Materials

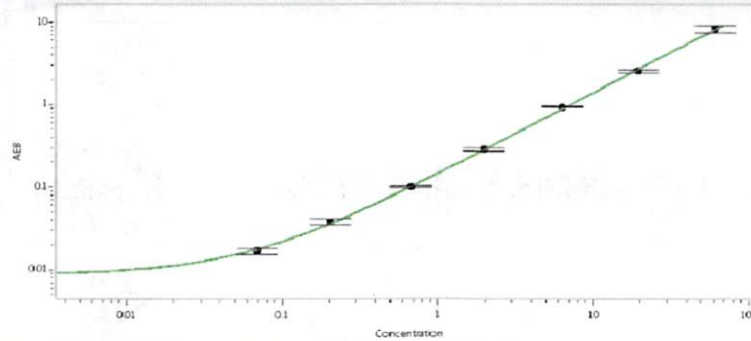
Tau			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	920402	1.61	1.10-1.71
Control 2	920402	72.7	57.1-85.6
Panel 1	MF-101617-01	0.930	0.605-1.92
Panel 2	MF-101617-02	33.1	22.7-44.5
Panel 3	MF-101617-03	137	96.8-194
NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	920402	5.70	4.23-6.34
Control 2	920402	462	371-556
Panel 1	MF-101617-01	5.11	3.08-6.38
Panel 2	MF-101617-02	36.7	16.7-50.4
Panel 3	MF-101617-03	394	234-541
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	920402	22.7	16.5-24.7
Control 2	920402	1036	809-1213
Panel 1	MF-101617-01	61.4	40.9-62.1
Panel 2	MF-101617-02	214	111-227
Panel 3	MF-101617-03	3149	2085-3303
UCHL1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	920402	121	78.3-138
Control 2	920402	10710	7808-11713
Panel 1	MF-101617-01	70.1	42.9-87.8
Panel 2	MF-101617-02	473	249-503
Panel 3	MF-101617-03	6174	4232-6348

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

Review/Approval: Patrizia Stadler Assoc. Director, QC P. Shodler 18 Sep 2019

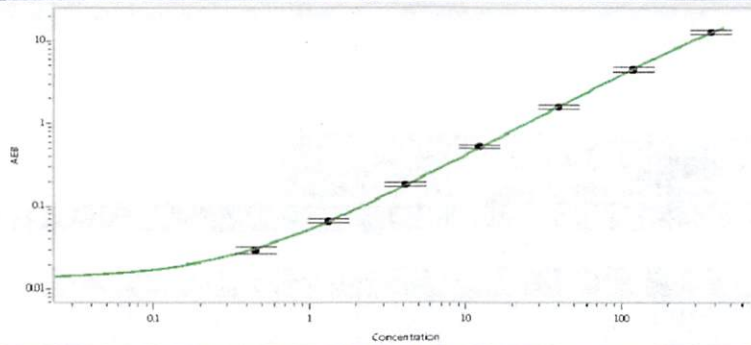
Name Title Signature/Date

Tau Calibration Curve



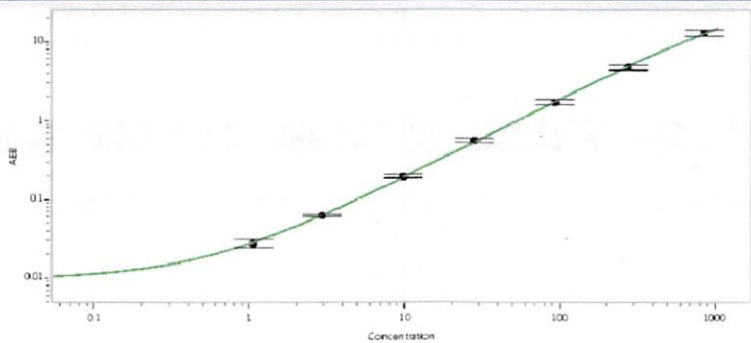
Calibrator Levels (pg/mL)	
A	0.000
B	0.070
C	0.204
D	0.679
E	1.99
F	6.34
G	19.7
H	61.3

NF-light® Calibration Curve



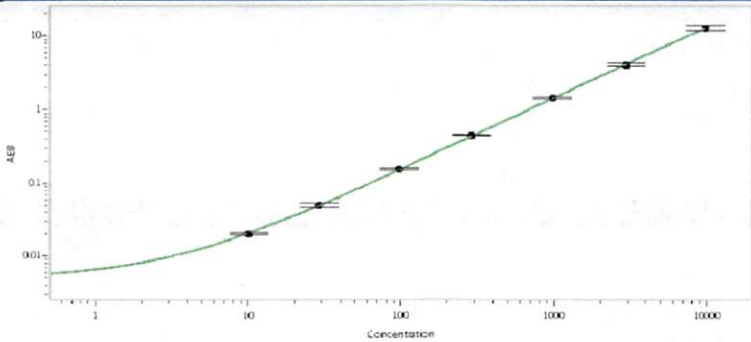
Calibrator Levels (pg/mL)	
A	0.000
B	0.455
C	1.31
D	4.12
E	12.2
F	40.0
G	120
H	376

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.06
C	2.93
D	9.77
E	28.2
F	93.6
G	277
H	855

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	10.1
C	28.9
D	97.8
E	288
F	981
G	2981
H	9782