

Product Number:	102153
Lot Number:	501157
Expiration:	23-Aug-2018
Platform(s):	HD-1 and SR-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102141	80762
Detector Reagent	102143	80760
SBG Reagent (HD-1)	102145	80544
Sample Diluent	102147	80543
Calibrators	102150	73200
Controls	102151-102152	73201
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

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

Release Materials

Tau

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	73201	2.73	1.65-3.24
Control 2	73201	105	86.6-130
Panel 1	MF-101617-01	1.11	0.532-2.26
Panel 2	MF-101617-02	27.0	23.4-35.1
Panel 3	MF-101617-03	131	107-161

NF-light®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	73201	6.12	4.92-7.38
Control 2	73201	423	364-566
Panel 1	MF-101617-01	6.11	3.00-7.55
Panel 2	MF-101617-02	40.4	27.5-48.4
Panel 3	MF-101617-03	455	363-544

GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	73201	20.8	12.5-24.4
Control 2	73201	832	656-984
Panel 1	MF-101617-01	61.1	40.6-69.8
Panel 2	MF-101617-02	173	145-252
Panel 3	MF-101617-03	2989	2397-3595

UCHL1®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	73201	138	100-171
Control 2	73201	11430	8573-12859
Panel 1	MF-101617-01	69.5	49.2-90.2
Panel 2	MF-101617-02	434	361-542
Panel 3	MF-101617-03	5240	4612-6919

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

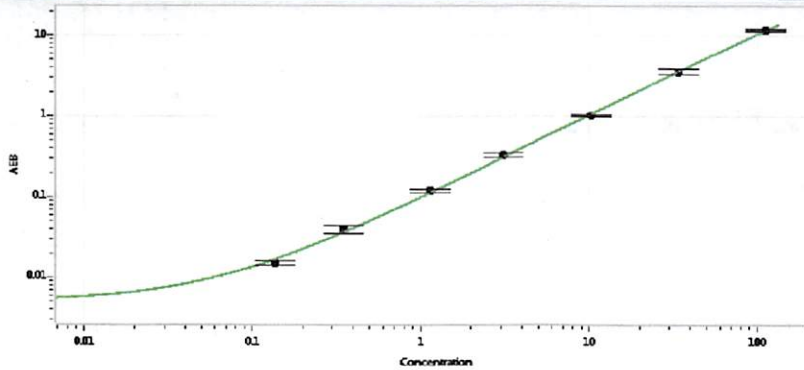
Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
22 MAR 2018
Signature/Date

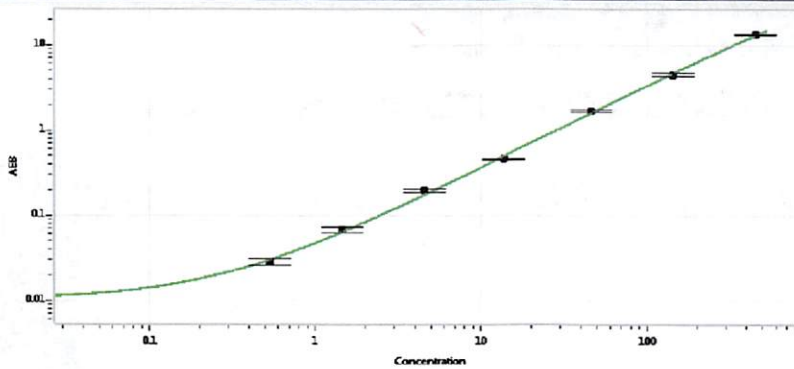
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.136
C	0.348
D	1.14
E	3.11
F	10.3
G	33.9
H	112

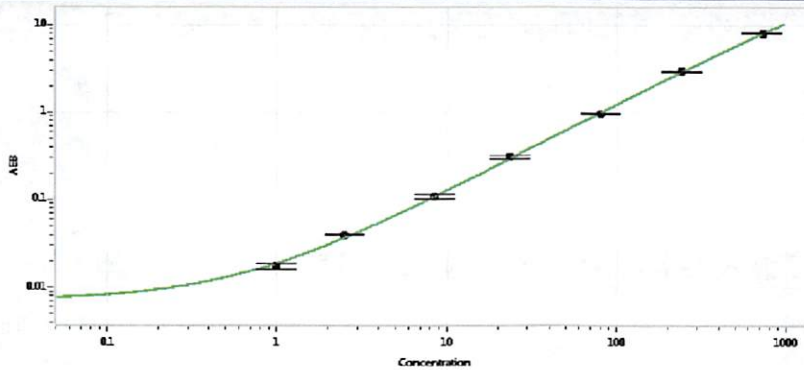
NF-light® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.533
C	1.46
D	4.57
E	13.6
F	45.8
G	143
H	453

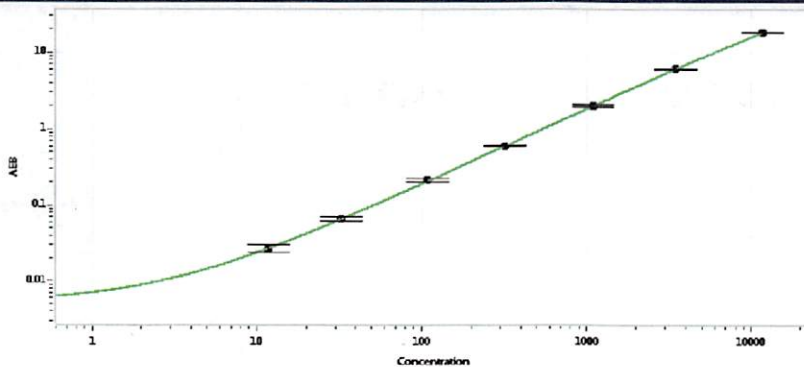
GFAP® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.987
C	2.52
D	8.49
E	23.7
F	80.5
G	243
H	725

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	11.9
C	32.8
D	108
E	322
F	1106
G	3492
H	11636