

Release Materials*

Tau

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	408602	2.13	1.45 - 2.65
Control 2	408602	90.8	74.0 - 112

NF-Light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	408602	6.13	4.84 - 7.49
Control 2	408602	482	414 - 621

GFAP®

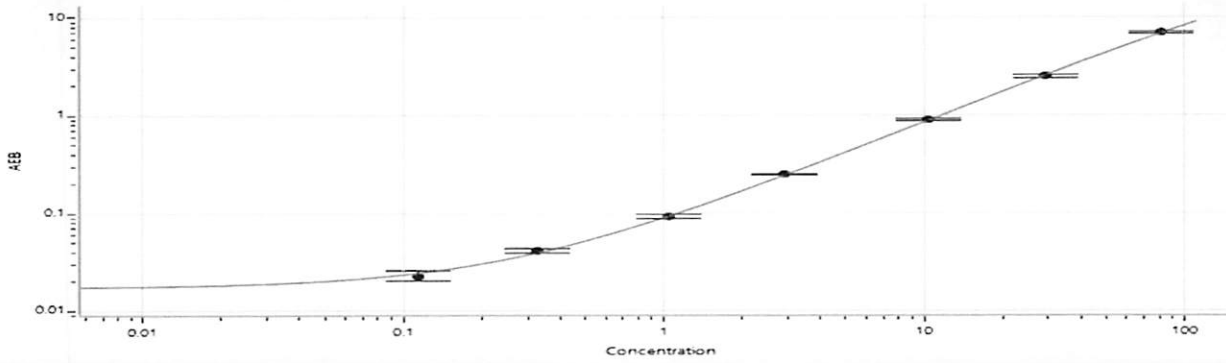
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	408602	93.4	72.9 - 113
Control 2	408602	7741	6697 - 10348

UCHL1®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	408602	115	98.8 - 151
Control 2	408602	9778	8446 - 12669

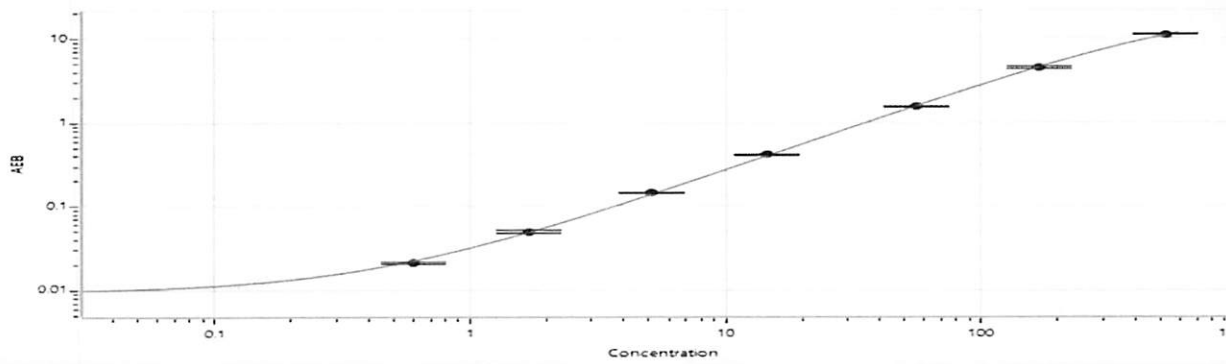
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Tau Calibrator Curve



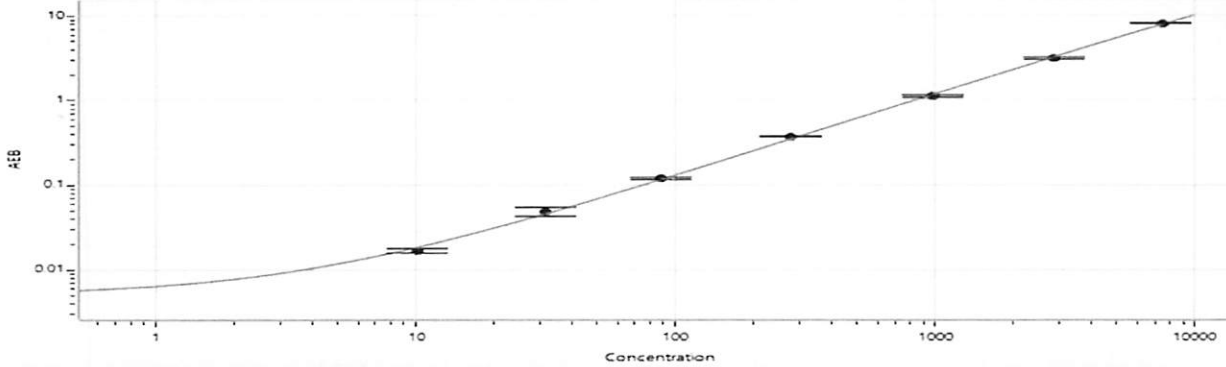
Calibrator Levels (pg/mL)	
A	0.000
B	0.114
C	0.328
D	1.05
E	2.91
F	10.4
G	29.3
H	82.1

NF-Light® Calibrator Curve



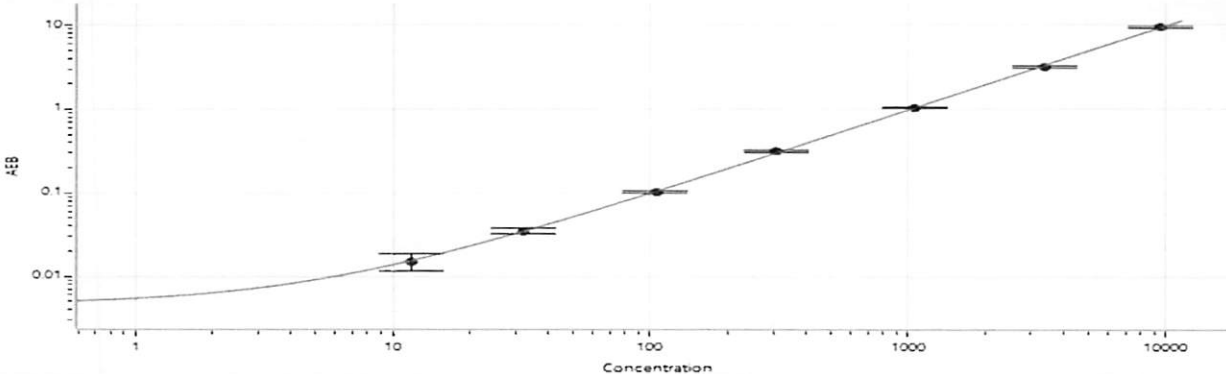
Calibrator Levels (pg/mL)	
A	0.000
B	0.605
C	1.70
D	5.16
E	14.6
F	55.6
G	170
H	529

GFAP® Calibrator Curve



Calibrator Levels (pg/mL)	
A	0.000
B	10.1
C	31.4
D	87.6
E	279
F	977
G	2844
H	7418

UCHL1® Calibrator Curve



Calibrator Levels (pg/mL)	
A	0.000
B	11.8
C	31.9
D	105
E	309
F	1059
G	3375
H	9490