

Release Materials*

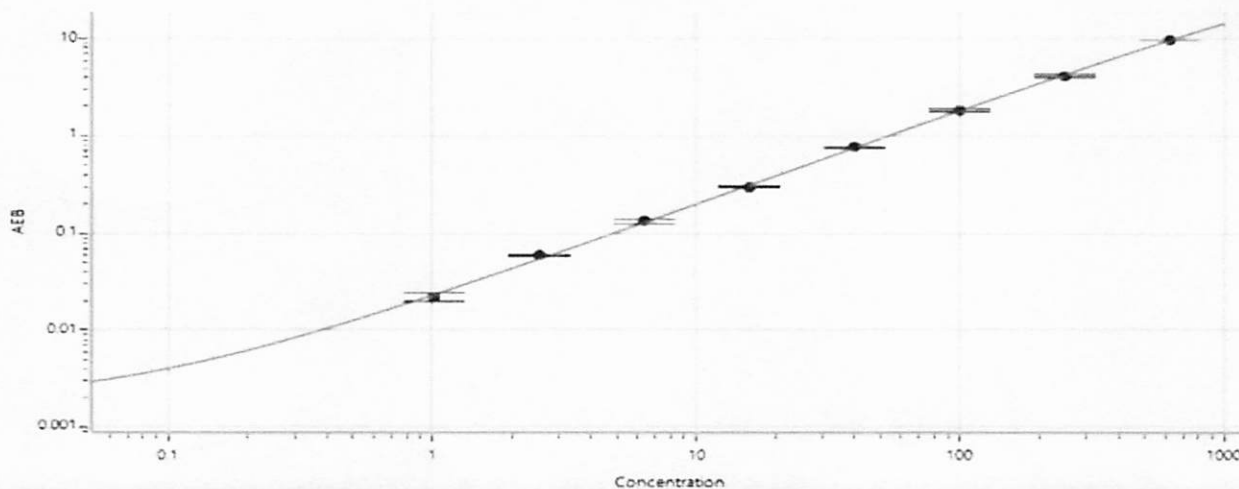
GFAP®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	402949	87.3	59.6 - 111
Control 2	402950	1097	776 - 1441

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

Calibrator Concentrate Reconstitution Volume (uL)	1144
Control 1 Concentrate Reconstitution Volume (uL)	1144
Control 2 Concentrate Reconstitution Volume (uL)	1144

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	1.02
C	2.55
D	6.35
E	16.0
F	39.7
G	99.0
H	248
I	621

Release Materials*

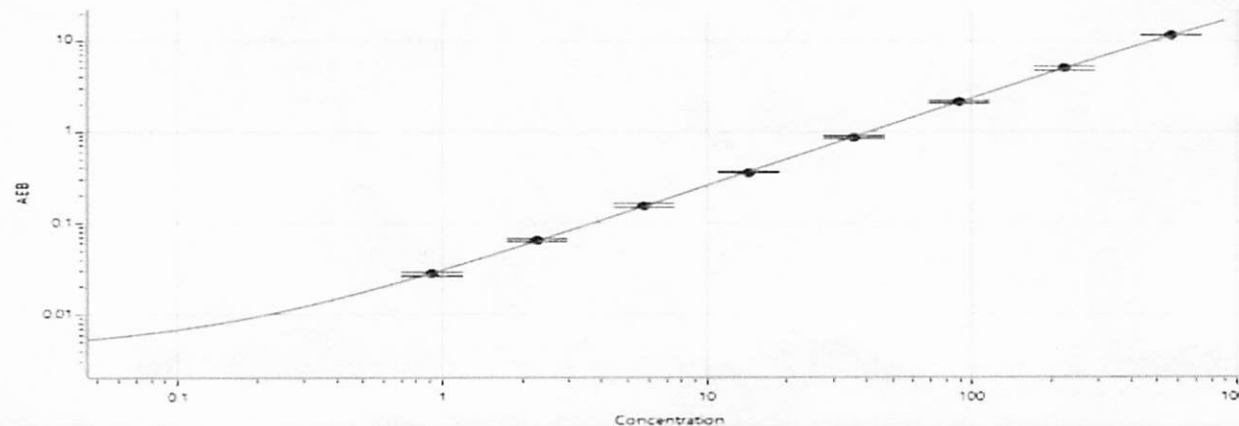
NF-Light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	402949	84.1	56.5 - 105
Control 2	402950	1048	734 - 1362

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

Calibrator Concentrate Reconstitution Volume (uL)	1144
Control 1 Concentrate Reconstitution Volume (uL)	1144
Control 2 Concentrate Reconstitution Volume (uL)	1144

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.915
C	2.29
D	5.72
E	14.3
F	35.7
G	89.4
H	223
I	559