

Release Materials*

Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	411006	19.0	15.1 - 22.7
Control 2	411006	129	106 - 158

Abeta 42

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	411006	6.84	5.45 - 8.18
Control 2	411006	31.4	25.4 - 38.1

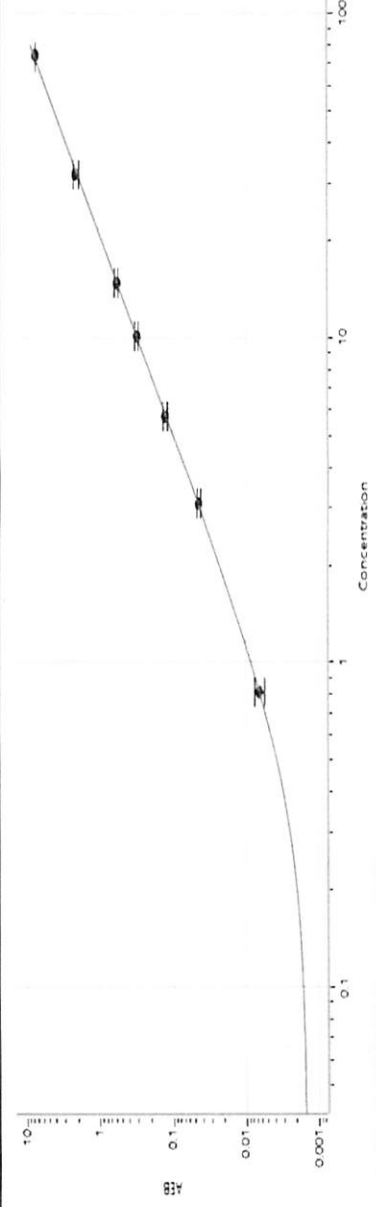
GFAP®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	411006	206	160 - 240
Control 2	411006	3174	2597 - 3895

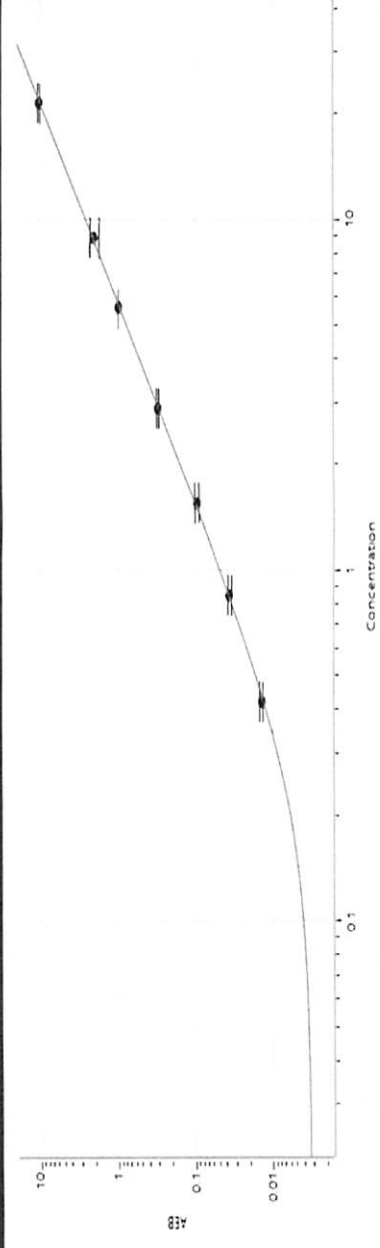
NF-Light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	411006	27.9	21.9 - 32.9
Control 2	411006	407	310 - 561

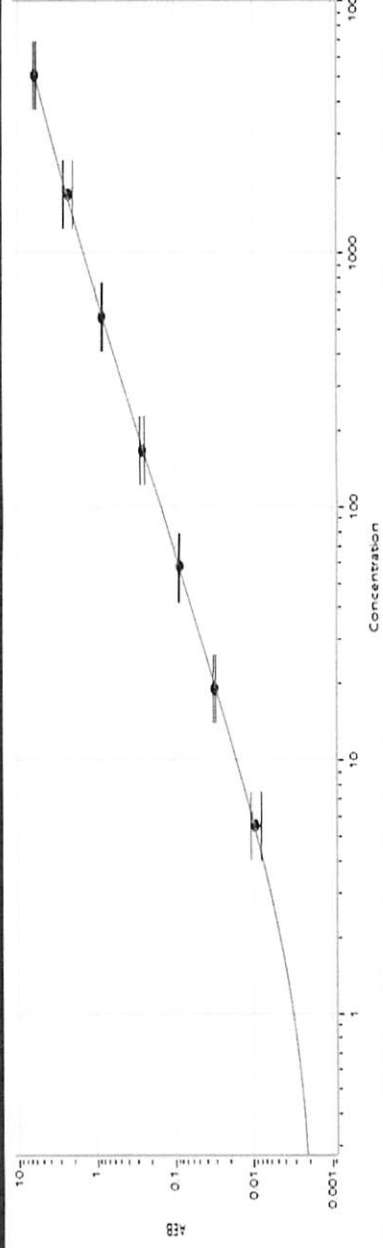
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Abeta 40 Calibrator Curve

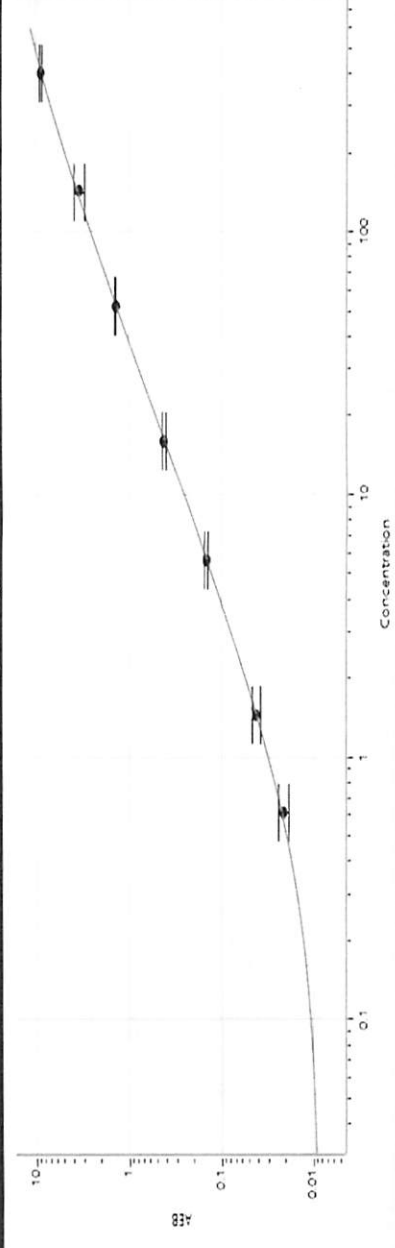
Calibrator	Concentration (pg/mL)
A	0.000
B	0.813
C	3.10
D	5.71
E	10.1
F	14.8
G	32.0
H	73.6

Abeta 42 Calibrator Curve

Calibrator	Concentration (pg/mL)
A	0.000
B	0.422
C	0.846
D	1.55
E	2.88
F	5.58
G	8.85
H	21.5

GFAP® Calibrator Curve

Calibrator	Concentration (pg/mL)
A	0.000
B	5.49
C	19.2
D	58.0
E	167
F	560
G	1728
H	5043

NF-Light® Calibrator Curve

Calibrator	Concentration (pg/mL)
A	0.000
B	0.611
C	1.44
D	5.58
E	15.8
F	52.0
G	142
H	400