



**Release Materials\***

**Abeta 40**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Control 1</b>	329706	18.1	14.6 - 21.9
<b>Control 2</b>	329706	100	89.8 - 149

**Abeta 42**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Control 1</b>	329706	7.76	6.12 - 9.18
<b>Control 2</b>	329706	37.3	30.3 - 45.4

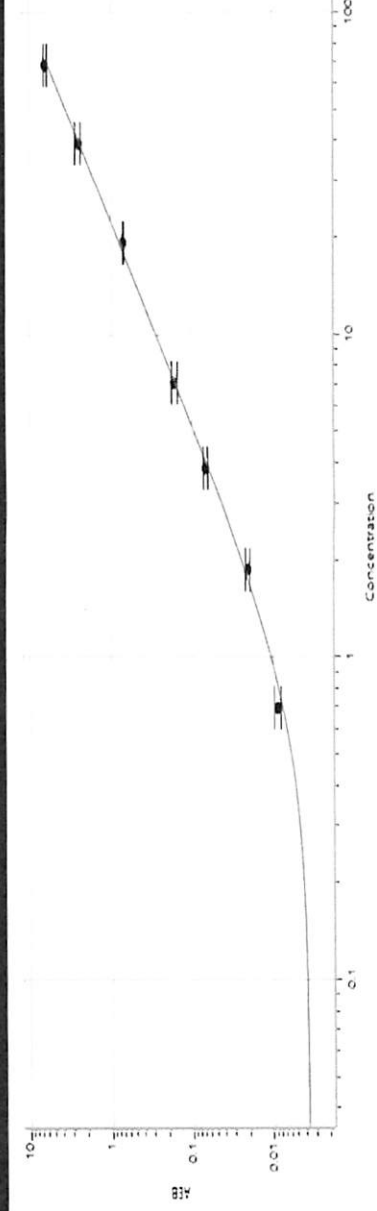
**GFAP®**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Control 1</b>	329706	202	161 - 242
<b>Control 2</b>	329706	4375	3450 - 5174

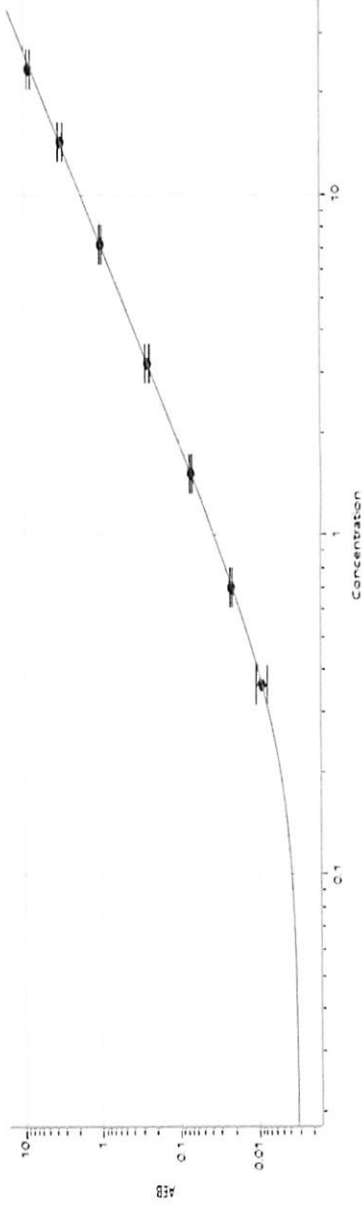
**NF-Light®**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Control 1</b>	329706	24.8	19.6 - 29.5
<b>Control 2</b>	329706	559	459 - 691

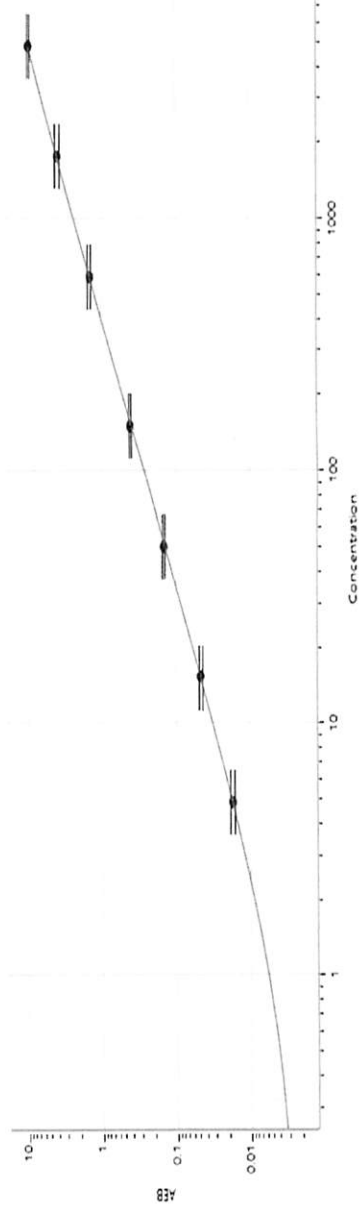
\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

**Abeta 40 Calibrator Curve**Calibrator Levels  
(pg/mL)

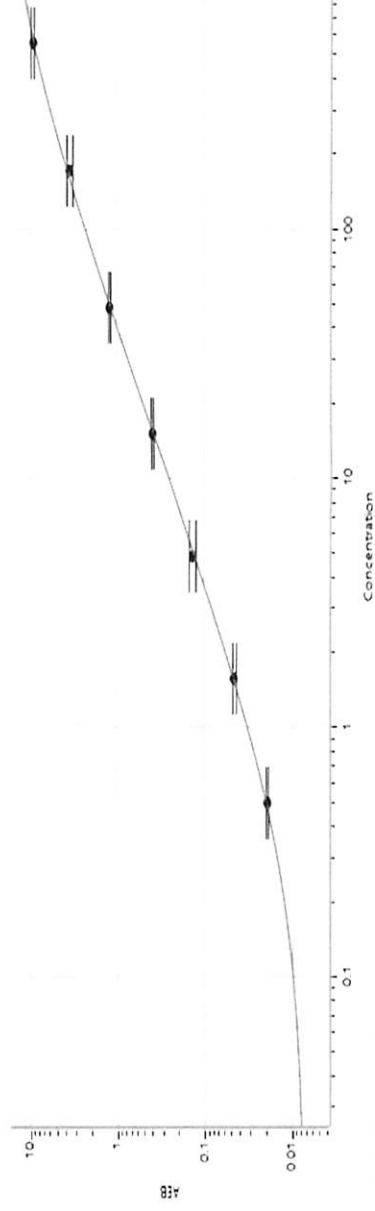
A	0.000
B	0.693
C	1.87
D	3.83
E	7.07
F	19.2
G	39.4
H	68.3

**Abeta 42 Calibrator Curve**Calibrator Levels  
(pg/mL)

A	0.000
B	0.360
C	0.700
D	1.50
E	3.20
F	7.11
G	14.3
H	23.5

**GFAP® Calibrator Curve**Calibrator Levels  
(pg/mL)

A	0.000
B	4.88
C	15.2
D	49.8
E	151
F	591
G	1757
H	4803

**NF-Light® Calibrator Curve**Calibrator Levels  
(pg/mL)

A	0.000
B	0.498
C	1.56
D	4.82
E	15.0
F	48.3
G	170
H	558