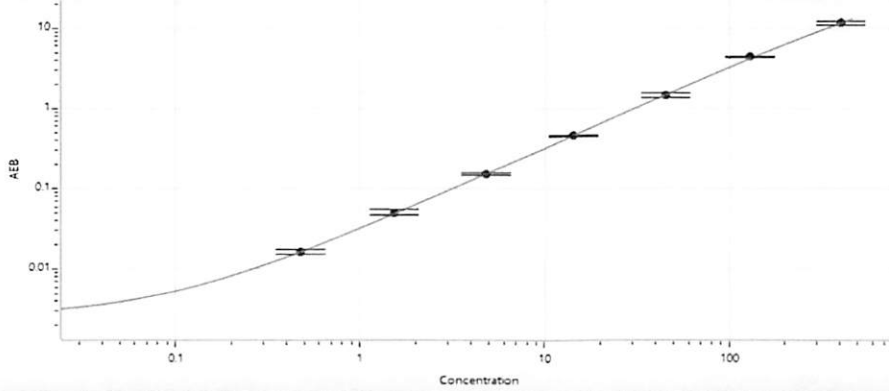


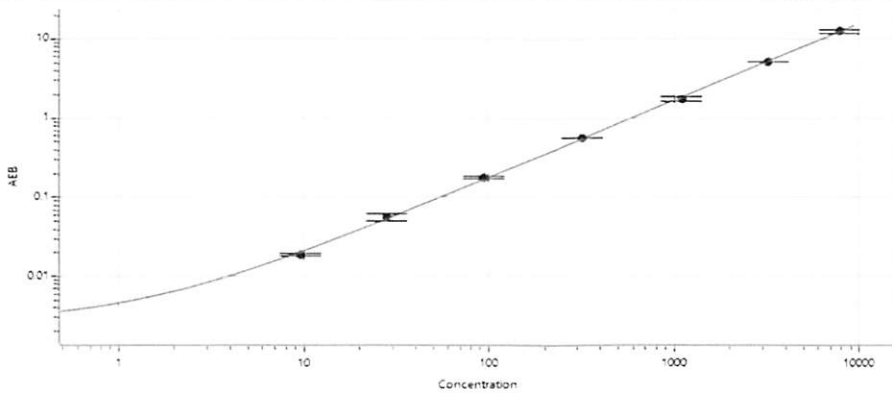
NF-Light® Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.476
C	1.54
D	4.86
E	14.3
F	45.4
G	131
H	405

GFAP® Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	9.54
C	28.1
D	94.1
E	317
F	1097
G	3219
H	7838