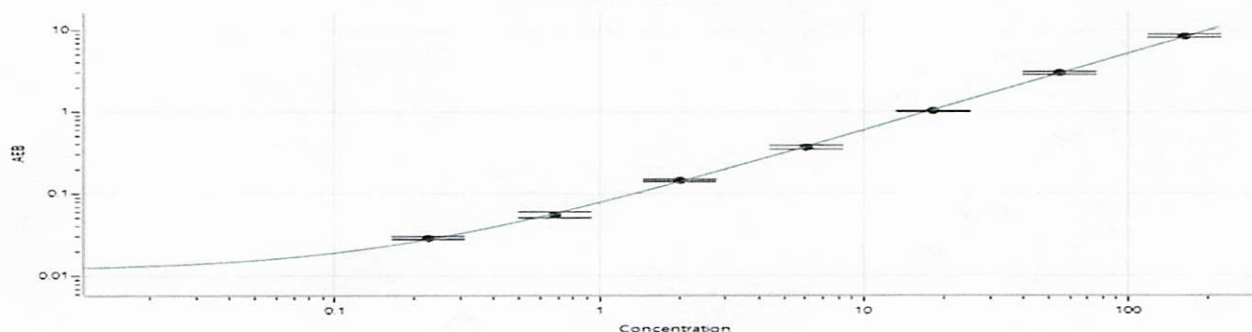


Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	430516	7.52	5.81 - 10.8
Control 2 Concentrate	430517	128	91.6 - 170

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

Calibrator Concentrate Reconstitution Volume (uL)	676
Control 1 Concentrate Reconstitution Volume (uL)	676
Control 2 Concentrate Reconstitution Volume (uL)	676



A	0.000
B	0.225
C	0.675
D	2.02
E	6.07
F	18.2
G	54.7
H	164

 Name Title Signature/Date