

Product Number:	100196
Lot Number:	504250
Expiration:	18-Jan-2025
Platform(s):	HD-1/HD-X



Component	Part Number	Lot Number	Lot Number
Bead Reagent	100391	324306	N/A
Detector Reagent	100315	324805	N/A
SBG Reagent	100316	324803	N/A
Sample Diluent	100317	319916	319916A
Calibrator Diluent	100665	324009	N/A
Calibrator Concentrate	100664	320709	N/A
RGP Reagent	103159	N/A ¹	N/A ¹

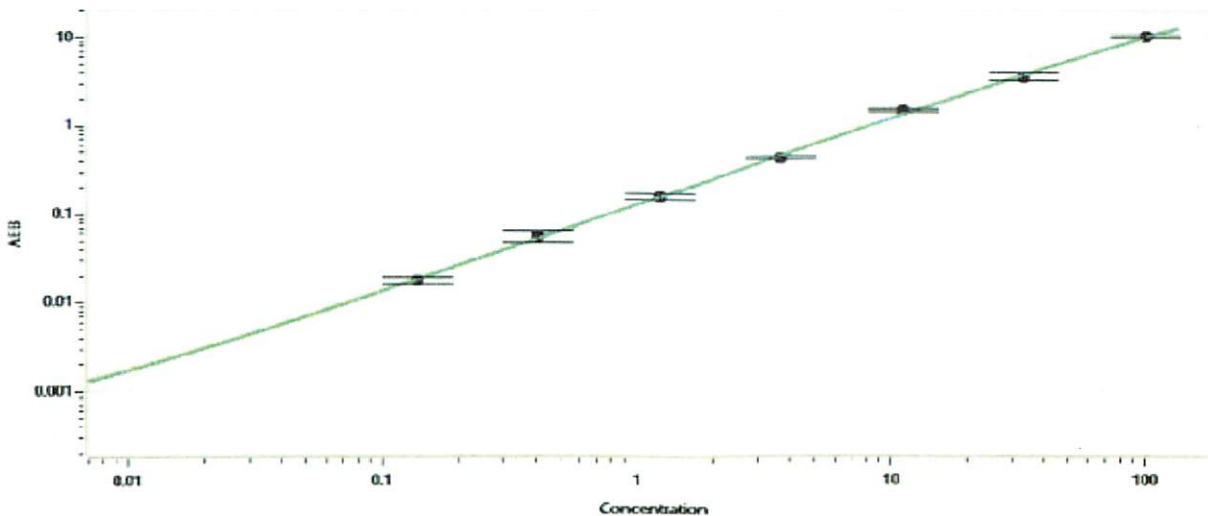
¹ Reagents are not Kit Lot Specific

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

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	123410	8.30	5.34 - 13.4
Control 2	123410	83.7	51.9 - 143
Panel 1	029002	1.61	1.19 - 2.82
Panel 2	029002	7.42	5.13 - 14.0
Panel 3	029002	81.1	46.8 - 134

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

Calibrator Concentrate Value Assignment (pg/mL)	1910
--------------------------------------------------------	------



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval

Binh Vu
Name

QC Supervisor
Title

[Signature] 01MAY2024
Signature/Date