

Product Number:	103836
Lot Number:	504171
Expiration:	10-Apr-2025
Platform(s):	HD-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	103820	328906	N/A
Detector Reagent	103822	329003	N/A
SBG Reagent	103824	329004	329004A
Sample Diluent	103826	328407	N/A
Calibrator Diluent	103829	328315	N/A
Control Diluent	104435	328316	N/A
Calibrator Concentrate	104443	335302	N/A
Control 1 Concentrate	104445	335303	N/A
Control 2 Concentrate	104446	335304	N/A
RGP 2.0 Reagent	104321	N/A <sup>1</sup>	N/A <sup>1</sup>
Activation Buffer Reagent	104324	N/A <sup>1</sup>	N/A <sup>1</sup>

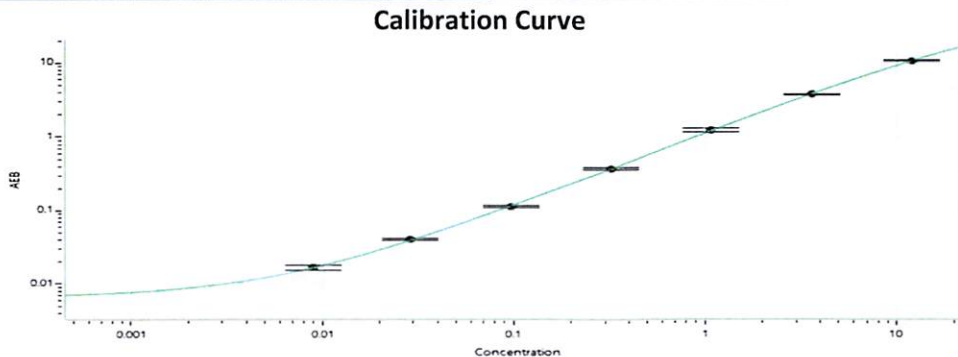
<sup>1</sup> Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	335303	0.174	0.074 - 0.251
Control 2 Concentrate	335304	3.50	2.15 - 4.00
Panel 1	204807	0.047	0.032 - 0.060
Panel 2	204807	0.039	0.025 - 0.055
Panel 3	204807	1.23	0.932 - 1.73
Panel 4	204807	2.77	1.75 - 3.64

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

Calibrator Concentrate Reconstitution Volume (uL)	1045
Control 1 Concentrate Reconstitution Volume (uL)	1045
Control 2 Concentrate Reconstitution Volume (uL)	1045

Calibrator Levels (pg/mL)	
A	0
B	0.009
C	0.029
D	0.098
E	0.325
F	1.08
G	3.60
H	12.0



Review/Approval

Cody Carlson                      QC Manager  
 \_\_\_\_\_  
 Name                                      Title

*Cody Carlson* 12 MAR 2024  
 \_\_\_\_\_  
 Signature/Date