

Product Number:	100860
Lot Number:	504041
Expiration:	27-Nov-2024
Platform(s):	HD-1/HD-X



Component	Part Number	Lot Number
Bead Reagent	100865	333403
Detector Reagent	100862	333512
SBG Reagent	100868	333603
Sample Diluent	100870	333107
Calibrators	101111	331808
Control 1	100875	331809
Control 2	100876	331809
RGP Reagent	103159	N/A <sup>1</sup>

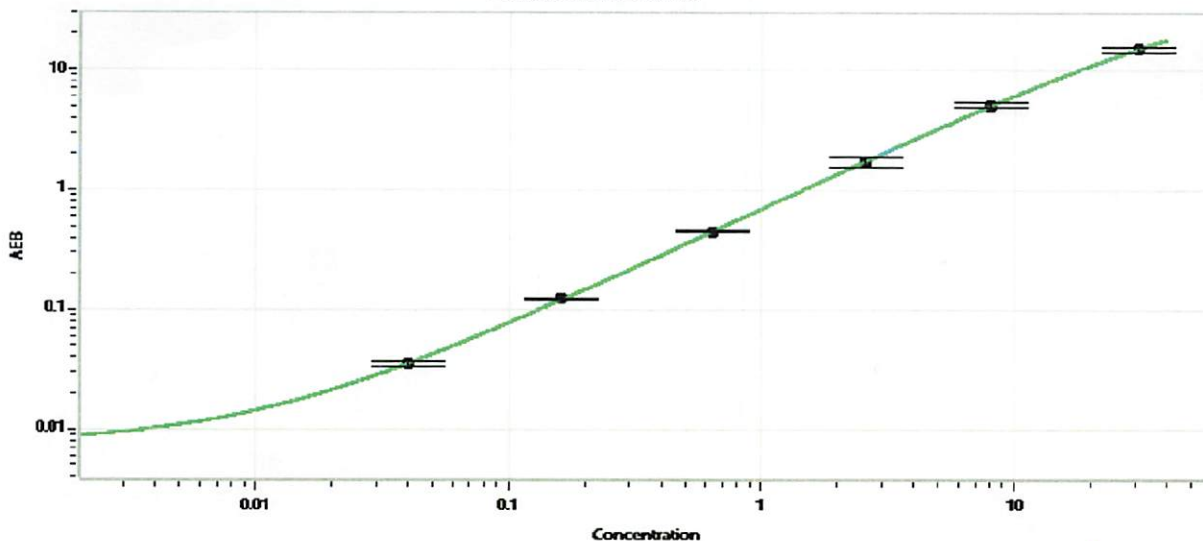
<sup>1</sup> Reagents are not Kit Lot Specific

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

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	331809	6.00	4.50 - 6.75
Control 2	331809	0.125	0.097 - 0.145
Panel 1	306704	0.688	0.534 - 1.26
Panel 2	306704	2.81	1.37 - 7.72
Panel 3	306704	17.9	11.2 - 24.4

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.040
C	0.161
D	0.634
E	2.55
F	8.06
G	30.5

Review/Approval

ROBERT SOMOGYI

QC MANAGER

*[Signature]* 22DEC2023

Name

Title

Signature/Date