

Product Number:	103277
Lot Number:	503423
Expiration:	24-Nov-2022
Platform(s):	HD-1/ HDX



Component	Part Number	Lot Number
Bead Reagent	103265	210917
Detector Reagent	103268	211004
SBG Reagent	103269	211007
Sample Diluent	103270	211008
Calibrator Diluent	103266	211009
Calibrator Concentrate	103272	032905
RGP Reagent	103159	N/A <sup>1</sup>

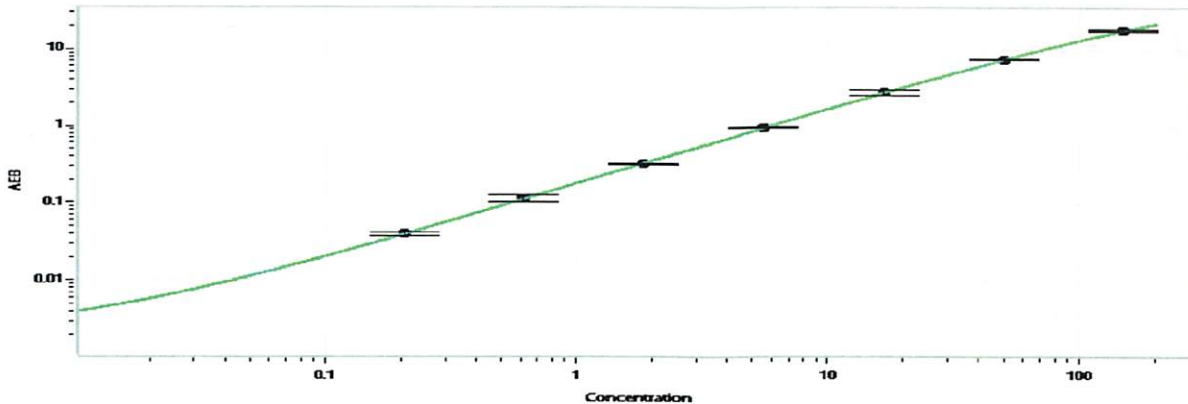
Data below represents results generated on the Simoa® HD-1 /HD-X Analyzer.

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	203709	3.54	2.82-4.22	
Control 2	203709	84.8	66.8-114	
Panel 1	205507	1.90	1.35-2.47	
Panel 2	205507	31.4	22.4-46.5	
Panel 3	205507	77.8	61.1-112	

Calibrator Concentrate Value Assignment (pg/mL)	5358
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.206
C	0.617
D	1.85
E	5.56
F	16.7
G	50.0
H	150

Review/Approval

Richa Shrestha  
Name

QC Supervisor  
Title

*[Signature]*  
03 May 2022  
Signature/Date