

Product Number:	101599
Lot Number:	503336
Expiration:	22-Jan-2023
Platform(s):	HD-1/HD-X & SR-X



Component	Part Number	Lot Number
Bead Reagent	101601	204205
Detector Reagent	102012	204304
SBG Reagent	101602	204305
Sample Diluent	103886	204306
Calibrator Diluent	102014	202211
Calibrator Concentrate	100778	204206
RGP Reagent	103159	N/A <sup>1</sup>

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

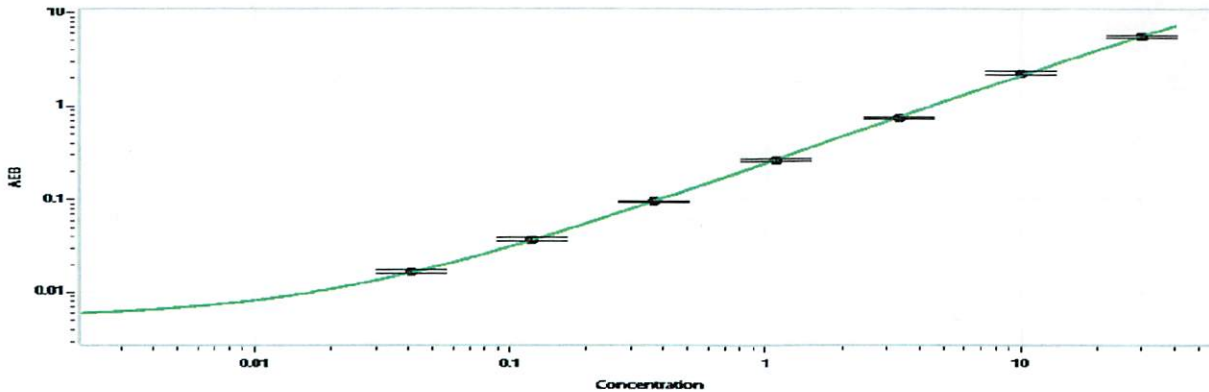
Data below represents results generated on the Simoa® HD-1 / HD-X & SR-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	209305	0.628	0.411-0.729
Control 2	209305	37.9	8.21-44.0
Panel 3	034908	11.1	8.12-12.5
Panel 4	034908	44.4	22.4-47.5

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

Calibrator Concentrate Value Assignment (pg/mL)	574
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 17 May 2022
	Name	Title	Signature/Date