

Product Number:	103186
Lot Number:	502554
Expiration:	14-Sep-2021
Platform(s):	HD-1/ HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	102246	026804
Detector Reagent	102248	026904
SBG Reagent	102250	026908
Sample Diluent	102252	026905
Calibrators	102255	027308
Control 1	102256	027309
Control 2	102257	027309
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

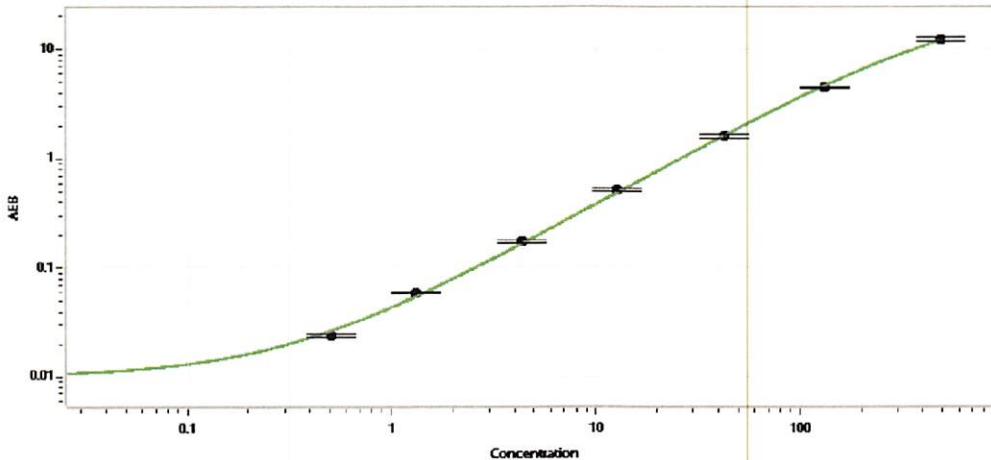


Data below represents results generated on the Simoa HD-1® Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	027309	3.79	2.93-4.39
Control 2	027309	145	120-180
Panel 1	926601	18.1	14.1-21.1
Panel 2	922602	6.69	4.94-7.41
Panel 3	922603	39.7	28.4-42.6
Panel 4	922604	150	129-204

* 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	0.505
C	1.31
D	4.39
E	12.7
F	42.5
G	132
H	487

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivam Goel 13 Oct 2020
Signature/Date