

Product Number:	103186
Lot Number:	502742
Expiration:	15-Jan-2022
Platform(s):	HD-1/ HD-X

Component	Part Number	Lot Number
Bead Reagent	102246	104109
Detector Reagent	102248	104101
SBG Reagent	102250	104015
Sample Diluent	102252	104001
Calibrators	102255	104002
Control 1	102256	104003
Control 2	102257	105001
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

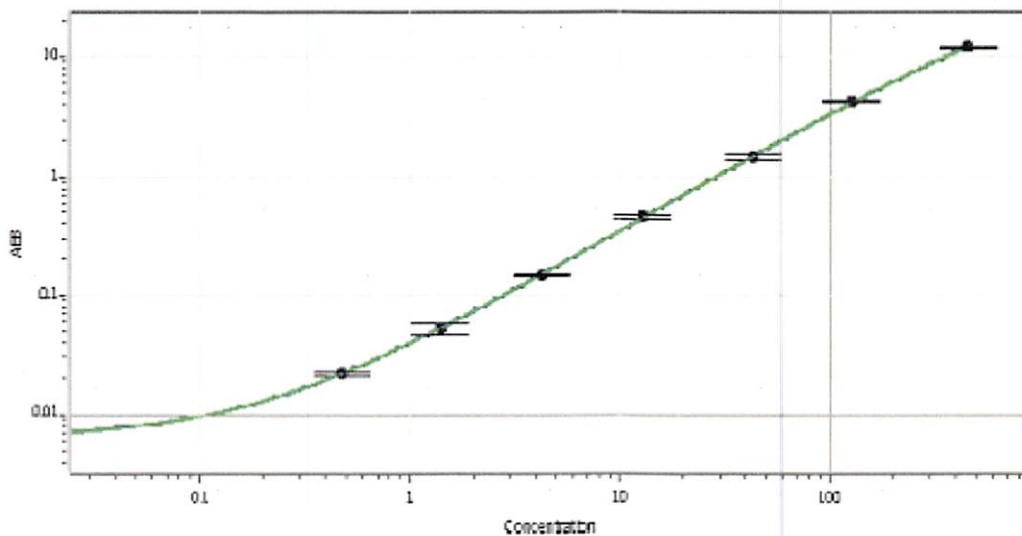


Data below represents results generated on the Simoa HD-1® Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	104003	3.72	2.41-4.12
Control 2	105001	165	138-207
Panel 1	922601	19.1	14.3 - 21.4
Panel 2	922602	7.10	4.97 - 7.45
Panel 3	922603	41.6	28.8 - 43.2
Panel 4	922604	185	133 - 201

* 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	0.479
C	1.40
D	4.27
E	12.9
F	43.3
G	128
H	451

Review/Approval

Junpil Park

QC Supervisor

Name

Title

Signature/Date

[Signature] 26 Feb 2021