

Product Number:	103714
Lot Number:	502923
Expiration:	9-Mar-2022
Platform(s):	HD-1 / HD-X & SR-X



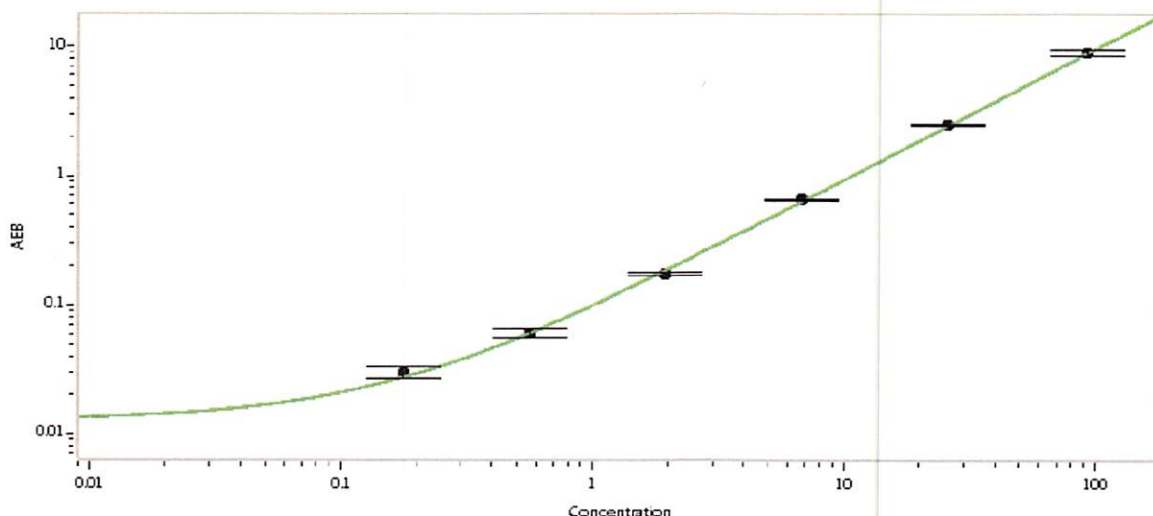
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	103698	112903
Detector Reagent	103700	112904
SBG Reagent	103702	112905
Sample Diluent	103704	113013
Reference Calibrator	103711	115405
Control 1	103712	115406
Control 2	103713	115406
RGP Reagent	103159	N/A ¹

¹ 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	115406	3.98	2.92-4.37	
Control 2	115406	125	105-157	
Panel 1	013903	0.722	0.258-1.11	
Panel 2	013903	1.86	0.841-2.44	
Panel 3	013903	3.07	1.79-4.31	

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.179
C	0.570
D	1.93
E	6.80
F	25.8
G	92.6

Review/Approval

Junpil Park

QC Supervisor

Name

Title

Signature/Date

[Signature] 10 Jun 2021