

Product Number: 100860
Lot Number: 503051
Expiration: 1-Jun-2022
Platform(s): HD-1/ HD-X

Component	Part Number	Lot Number
Bead Reagent	100865	119503
Detector Reagent	100862	119509
SBG Reagent	100868	119606
Sample Diluent	100870	119206
Calibrators	101111	116101
Control 1	100875	116102
Control 2	100876	116102
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

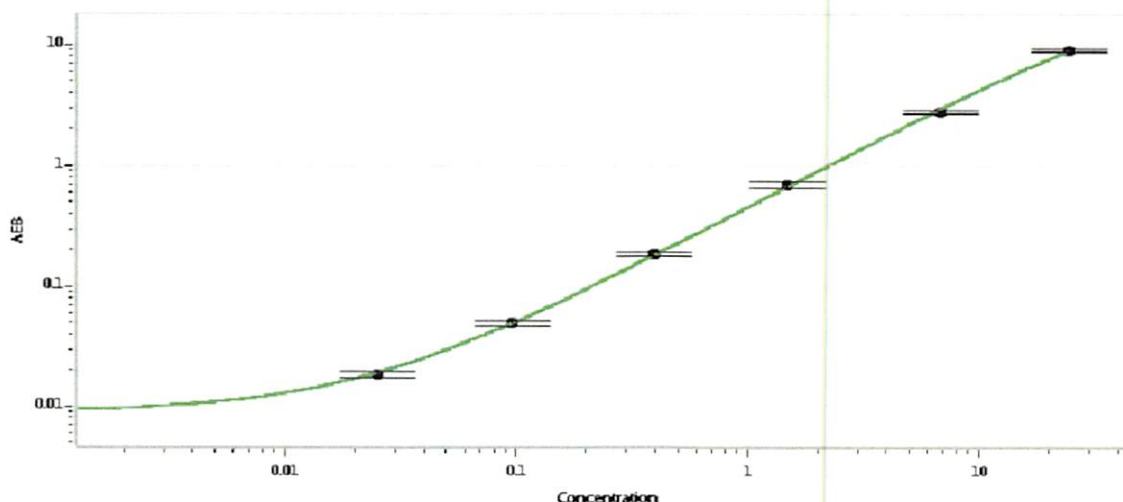


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	116102	4.19	3.47-5.43
Control 2	116102	0.081	0.051-0.101
Panel 1	115916	1.22	0.912-1.37
Panel 2	115916	4.82	3.92-9.41
Panel 3	115916	33.6	26.6-49.9

* 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 0.025
C 0.096
D 0.395
E 1.48
F 6.73
G 24.2

Review/Approval

Junpil Park

QC Supervisor

Name

Title

Signature/Date

 23 Jul 2021