

Product Number:	103714
Lot Number:	502691
Expiration:	11-Jul-2021
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103698	101109
Detector Reagent	103700	101210
SBG Reagent	103702	101203
Sample Diluent	103704	101207
Reference Calibrator	103711	101201
Control 1	103712	101202
Control 2	103713	102805
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

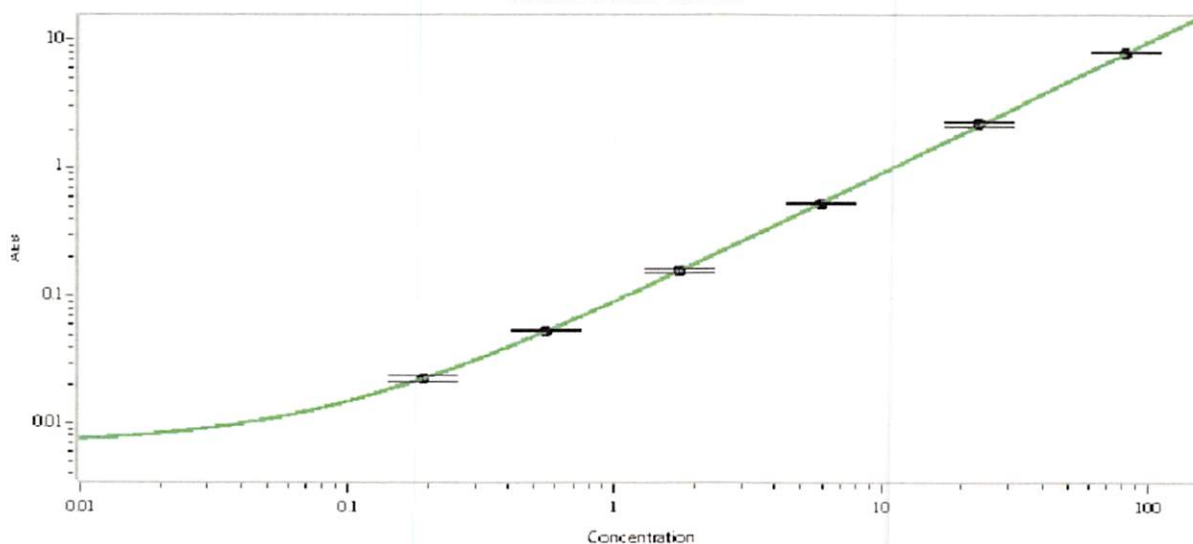


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	101202	3.34	2.57-3.89
Control 2	102805	80.6	50.5-93.8
Panel 1	013903	0.648	0.230-1.08
Panel 2	013903	1.54	0.855-2.26
Panel 3	013903	3.08	1.74-4.18

* 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	0.177
C	0.554
D	1.74
E	5.73
F	23.3
G	86.0

Review/Approval

Junpil Park

QC Supervisor

Name

Title

Signature/Date

 04 Feb 2021