

Product Number:	103714
Lot Number:	502613
Expiration:	11-May-2021
Platform(s):	HD-1 / HD-X & SR-X



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

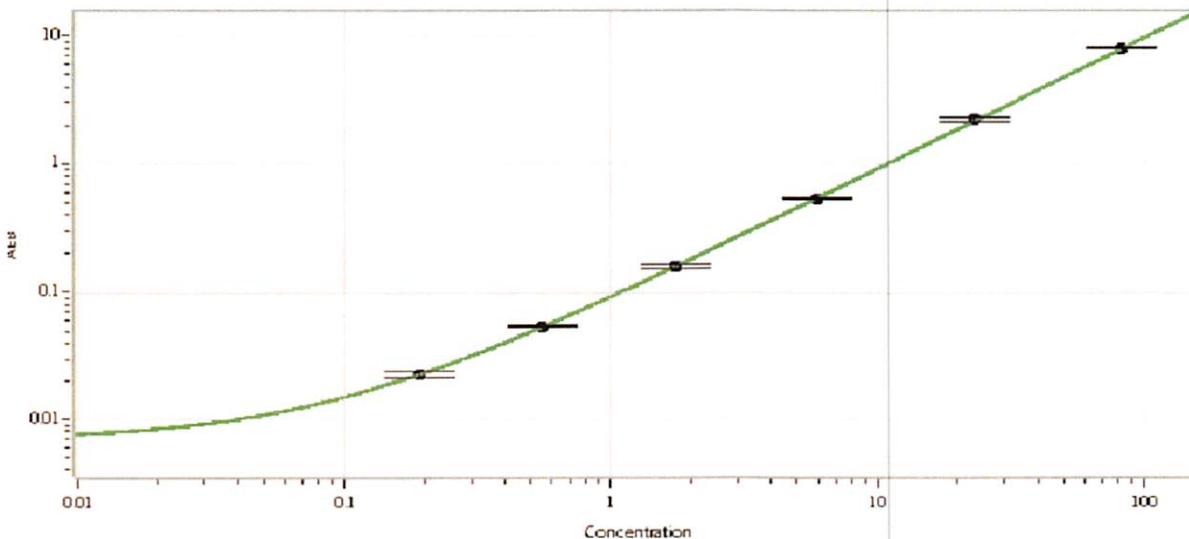
Component	Part Number	Lot Number
Bead Reagent	103698	031610
Detector Reagent	103700	031801
SBG Reagent	103702	031802
Sample Diluent	103704	031808
Reference Calibrator	103711	031601
Control 1	103712	031602
Control 2	103713	031602
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	031602	3.28	1.91-4.30
Control 2	031602	34.5	20.8-62.0
Panel 1	013903	0.355	0.230-1.08
Panel 2	013903	1.09	0.855-2.26
Panel 3	013903	2.50	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	0.193
C	0.552
D	1.74
E	5.96
F	23.3
G	82.3

Review/Approval

Shivani Goel

QC Manager

*Shivani Goel 03 Dec 2020*

Name

Title

Signature/Date