

Product Number:	101605
Lot Number:	502329
Expiration:	1-Apr-2021
Platform(s):	HD-1 / HD-X

Component	Part Number	Lot Number
Bead Reagent	101607	010507
Detector Reagent	101606	010702
SBG Reagent	101608	010904
Sample Diluent	103200	010703
Calibrator Diluent	101610	010705
Calibrator Concentrate	100895	010901
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



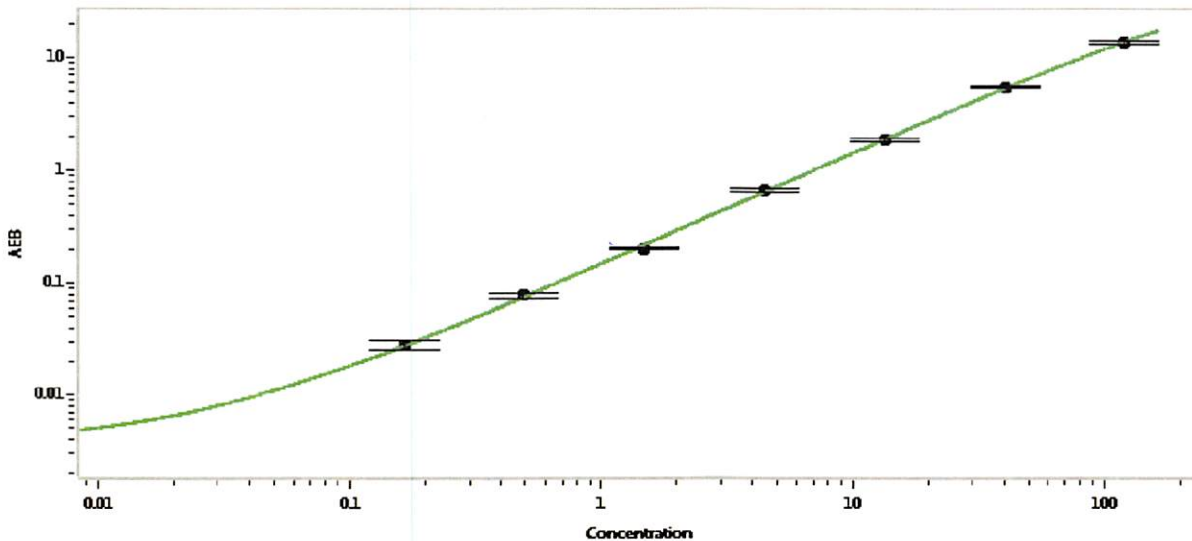
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	911406	2.29	1.09-3.60
Control 2	911406	150	87.5-204
Panel 1	829001	2.78	1.73-4.65
Panel 2	829002	4.19	2.946-9.02
Panel 3	829003	10.4	6.93-16.2
Panel 4	829004	49.6	38.4-64.9

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Value Assignment (pg/mL)	4866
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.165
C	0.494
D	1.48
E	4.44
F	13.3
G	40.0
H	120

Review/Approval

Shivani Goel QC Manager
Name Title

Shivani Goel 05 May 2020
Signature/Date