

Product Number:	101605
Lot Number:	502916
Expiration:	5-Mar-2022
Platform(s):	HD-1 / HD-X

Component	Part Number	Lot Number
Bead Reagent	101607	107606
Detector Reagent	101606	112613
SBG Reagent	101608	107718
Sample Diluent	103200	107717
Calibrator Diluent	101610	106420
Calibrator Concentrate	100895	108212
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

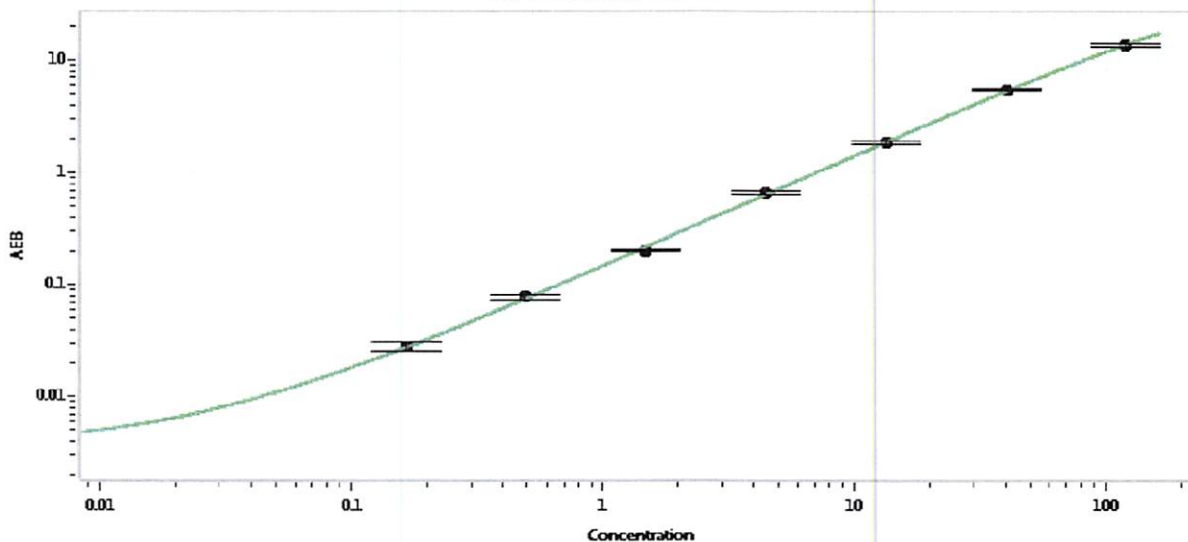


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*				* 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	108212	1.85	1.50-2.50	
Control 2	108212	119	90.0-150	
Panel 1	033901	4.20	0.659-6.76	
Panel 2	033901	0.347	0.264-0.576	
Panel 3	033901	20.7	19.6-29.4	
Panel 4	033901	61.4	53.7-80.6	

Calibrator Concentrate Value Assignment (pg/mL)	5874
---	------

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

Junpil Park
Name

QC Supervisor
Title

 23 Jun 2021
Signature/Date