

Product Number:	101605
Lot Number:	502948
Expiration:	7-May-2022
Platform(s):	HD-1 / HD-X



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

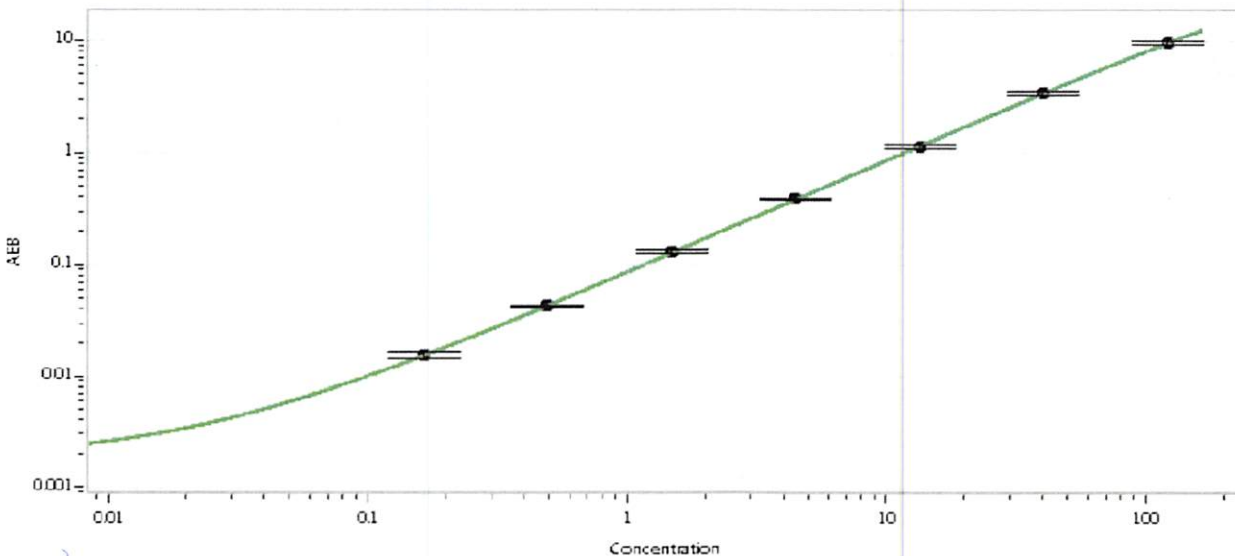
Component	Part Number	Lot Number
Bead Reagent	101607	114005
Detector Reagent	101606	117212
SBG Reagent	101608	117101
Sample Diluent	103200	117102
Calibrator Diluent	101610	112723
Calibrator Concentrate	100895	117610
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

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	117610	2.15	1.50-2.50	
Control 2	117610	128	90.0-150	
Panel 2	033901	0.549	0.252-0.618	
Panel 3	033901	29.5	19.1-31.1	
Panel 4	033901	79.1	54.0-83.1	

Calibrator Concentrate Value Assignment (pg/mL)	4195
---	------

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

 09 Jul 2024
Signature/Date