

Product Number:	101478
Lot Number:	503369
Expiration:	14-Feb-2023
Platform(s):	HD-1/HD-X & SR-X



Component	Part Number	Lot Number
Bead Reagent	101480	209605
Detector Reagent	101481	209807
SBG Reagent	101482	209808
Sample Diluent	101483	209810
Calibrators	100758-100765	213502
Control 1	100768	213503
Control 2	100769	213503
RGP Reagent	103159	N/A ¹

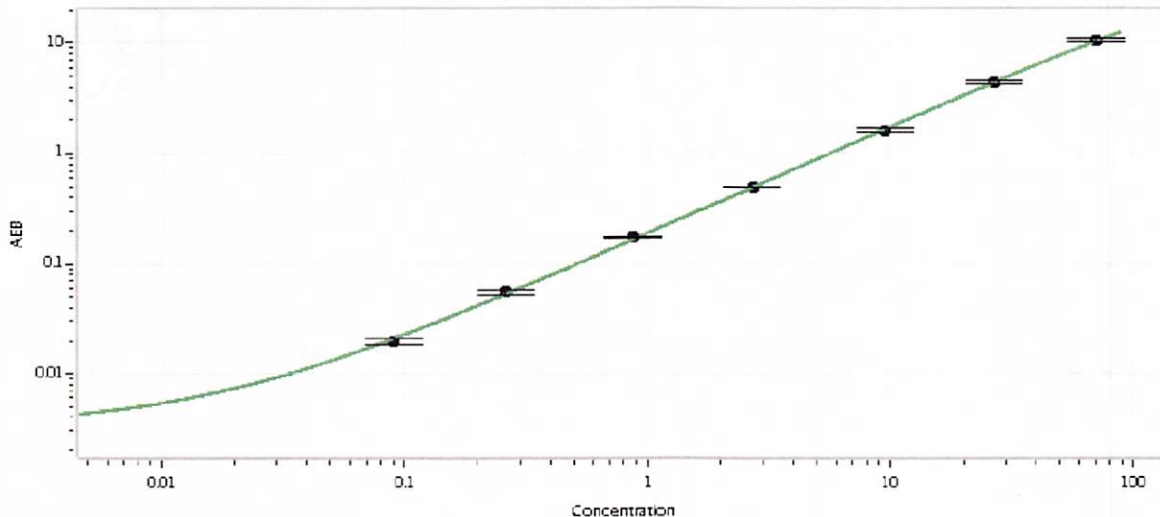
Data below represents results generated on the Simoa® HD-1 / HD-X & SR-X Analyzer.

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	213503	2.78	2.12-3.17
Control 2	213503	59.6	46.5-69.8
Panel 1	213804	1.03	0.433-1.27
Panel 2	122912	4.97	3.65-5.91
Panel 3	122912	54.7	41.5-68.9

* 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	0.090
C	0.260
D	0.878
E	2.71
F	9.56
G	26.9
H	70.0

Review/Approval Matt Smart Director, QC Matt Smart 09 Jun 2022
 Name Title Signature/Date