

Product Number:	103186
Lot Number:	502855
Expiration:	8-Mar-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	102246	106912
Detector Reagent	102248	106911
SBG Reagent	102250	106716
Sample Diluent	102252	106717
Calibrators	102255	106801
Control 1	102256	106818
Control 2	102257	109005
RGP Reagent	103159	N/A ¹

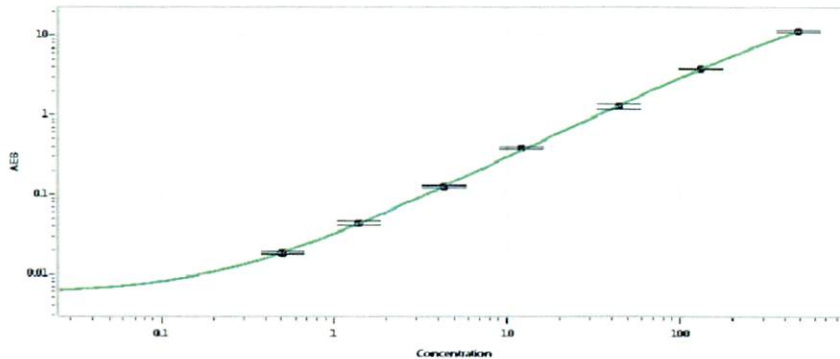
¹ RGP is not Kit Lot Specific

Release Materials*					
	Lot	Previously reported		Corrected **	
		Result (pg/mL)	Mean Range (pg/mL)	Result (pg/mL)	Mean Range (pg/mL)
Control 1	106818	3.90	2.73-5.70	3.14	2.27-4.32
Control 2	109005	235	180-270	188	144-250
Panel 1	008501	16.1	9.26-19.5	12.8	8.59-20.5
Panel 2	008501	13.7	7.42-14.5	10.9	7.39-15.1
Panel 3	008501	75.1	46.0-88.2	59.8	43.8-91.8
Panel 4	009401	216	144-253	172	128-302

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

** Values and ranges are updated as of 26Oct2021. Corrected result and calibrator levels should be used for reference.

Calibration Curve



	Calibrator Levels (pg/mL)	
	Previously reported	Corrected**
A	0.000	0.000
B	0.576	0.479
C	1.43	1.12
D	4.75	3.77
E	13.9	10.8
F	44.8	36.4
G	134	108
H	424	350

Review/Approval Shivani Goel QC Manager Shivani Goel 05Apr2022
Name Title Signature/Date