

Product Number:	101580
Lot Number:	502629
Expiration:	1-Sep-2021
Platform(s):	HD-1/HD-x & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101582	032213
Detector Reagent	101581	027002
SBG Reagent	101583	027003
Sample Diluent	103735	027005
Calibrator Diluent	101585	024509
Calibrator Concentrate	100655	932607
RGP Reagent	103159	N/A ¹

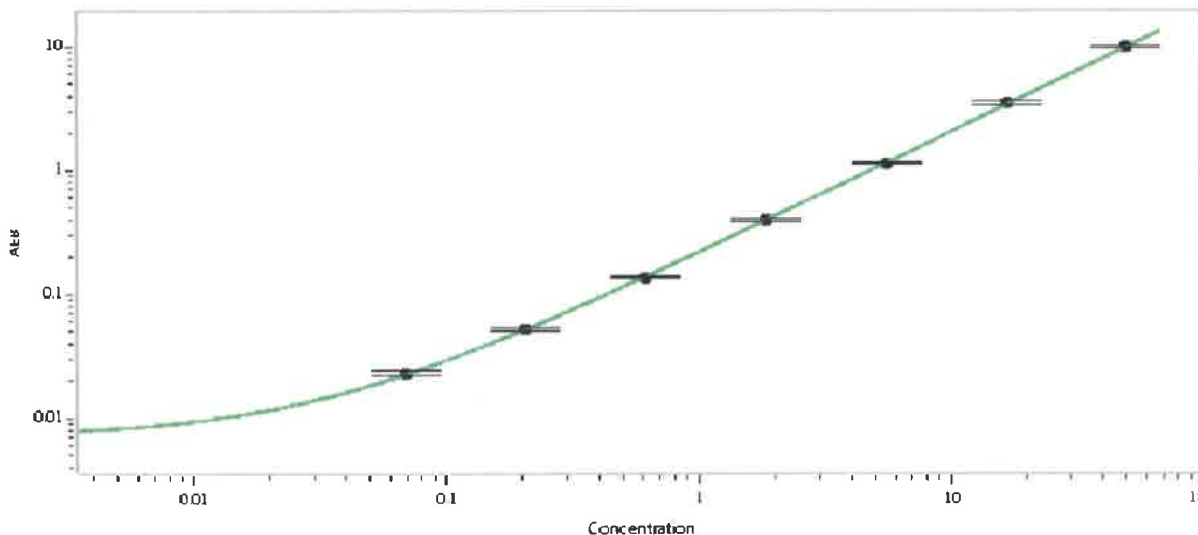
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-042817-01	3.07	2.27-3.41
Control 2	MF-051617-02	51.3	40.0-60.0
Panel 1	28301	30.3	22.3-33.5
Panel 2	28301	1.78	1.20-1.80
Panel 3	28301	2.47	1.70-2.55

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

Calibrator Concentrate Value Assignment (pg/mL)	1311
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.069
C	0.206
D	0.617
E	1.85
F	5.56
G	16.7
H	50.0

Review/Approval

Junpil Park

QC Supervisor

[Signature] 29 Jan 2021

Name

Title

Signature/Date

