

Product Number:	101580
Lot Number:	502938
Expiration:	3-Feb-2022
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	103601
Detector Reagent	101581	113712
SBG Reagent	101583	106314
Sample Diluent	103735	103405
Calibrator Diluent	101585	103802
Calibrator Concentrate	100655	104006
RGP Reagent	103159	N/A ¹

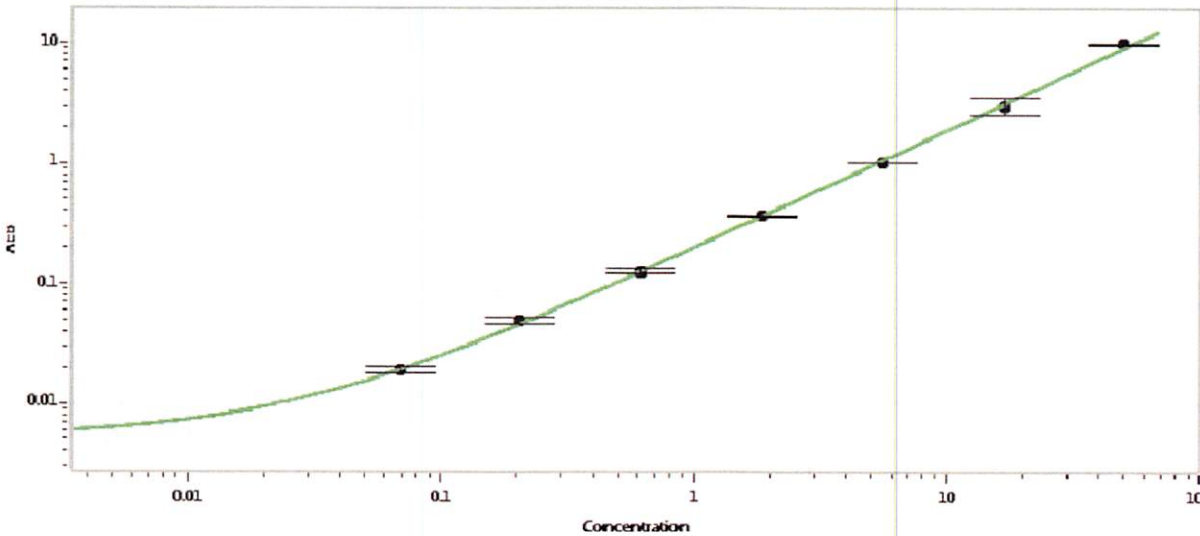
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	024603	1.49	0.606-2.36
Control 2	024603	27.1	17.1-35.5
Panel 1	116301	1.54	0.410-2.22
Panel 2	115317	4.01	2.64-4.92
Panel 3	116301	36.8	27.2-40.7

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

Calibrator Concentrate Value Assignment (pg/mL)	1090
---	------

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
Name

QC Supervisor
Title

[Signature] 17 Jun 2021
Signature/Date