

Product Number:	101580
Lot Number:	503061
Expiration:	14-Jul-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	119506
Detector Reagent	101581	119714
SBG Reagent	101583	119710
Sample Diluent	103735	119713
Calibrator Diluent	101585	119702
Calibrator Concentrate	100655	119711
RGP Reagent	103159	N/A ¹

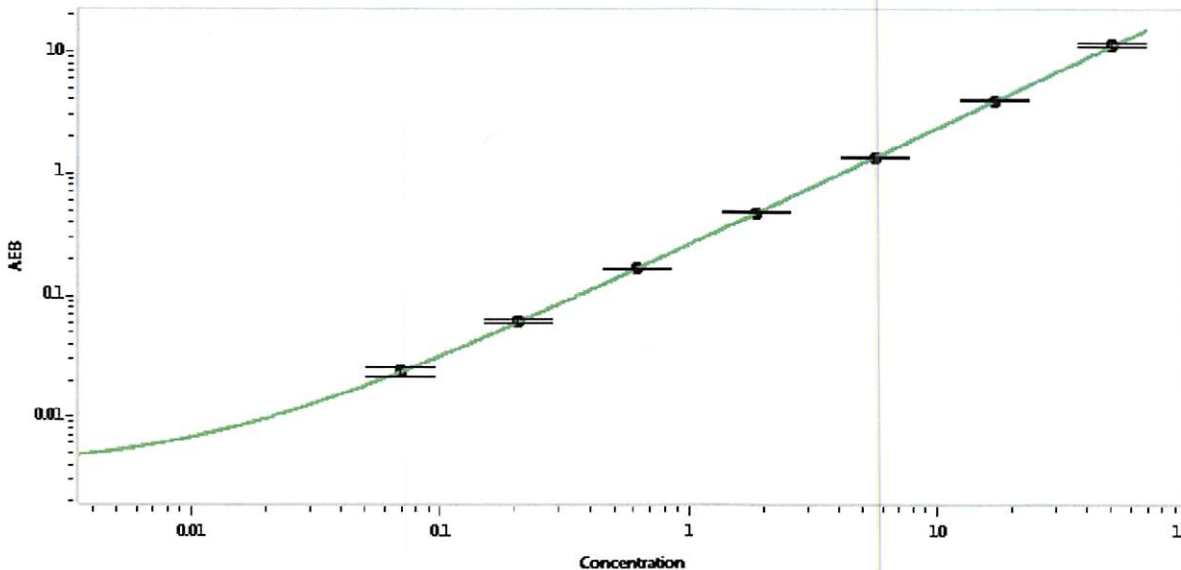
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	024603	1.40	0.762-2.33
Control 2	024603	27.8	19.1-34.9
Panel 2	115317	4.83	0.929-5.69
Panel 3	116301	42.7	16.0-47.2

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

Calibrator Concentrate Value Assignment (pg/mL)	1143
---	------

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/Date 28 Jul 2021