

Product Number:	101580
Lot Number:	503112
Expiration:	2-Aug-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	123013
Detector Reagent	101581	123401
SBG Reagent	101583	123402
Sample Diluent	103735	123308
Calibrator Diluent	101585	123403
Calibrator Concentrate	100655	123509
RGP Reagent	103159	N/A ¹

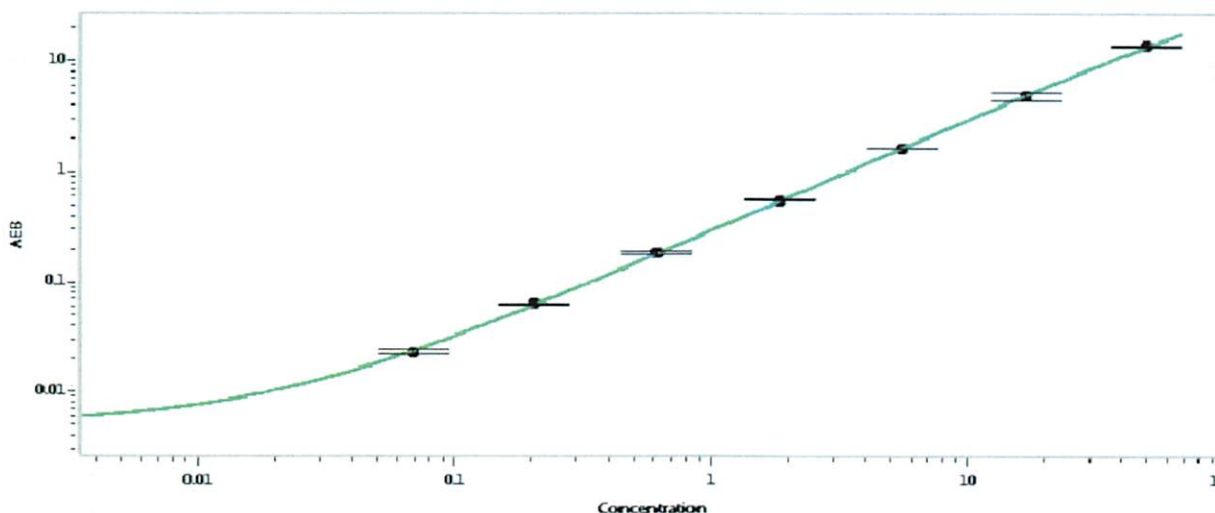
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	024603	1.04	0.594 - 2.32
Control 2	024603	17.6	13.9 - 36.9
Panel 2	115317	2.41	1.37 - 4.10
Panel 3	116301	20.3	13.1 - 41.8

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

Calibrator Concentrate Value Assignment (pg/mL)	999
---	-----

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 31 Aug 2021