

Product Number:	102860
Lot Number:	503122
Expiration:	2-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	102852	124506
Detector Reagent	102854	124508
SBG Reagent	102855	124509
Sample Diluent	102856	124602
Calibrator Diluent	102859	124603
Calibrator Concentrate	102857	124212
RGP Reagent	103159	N/A ¹

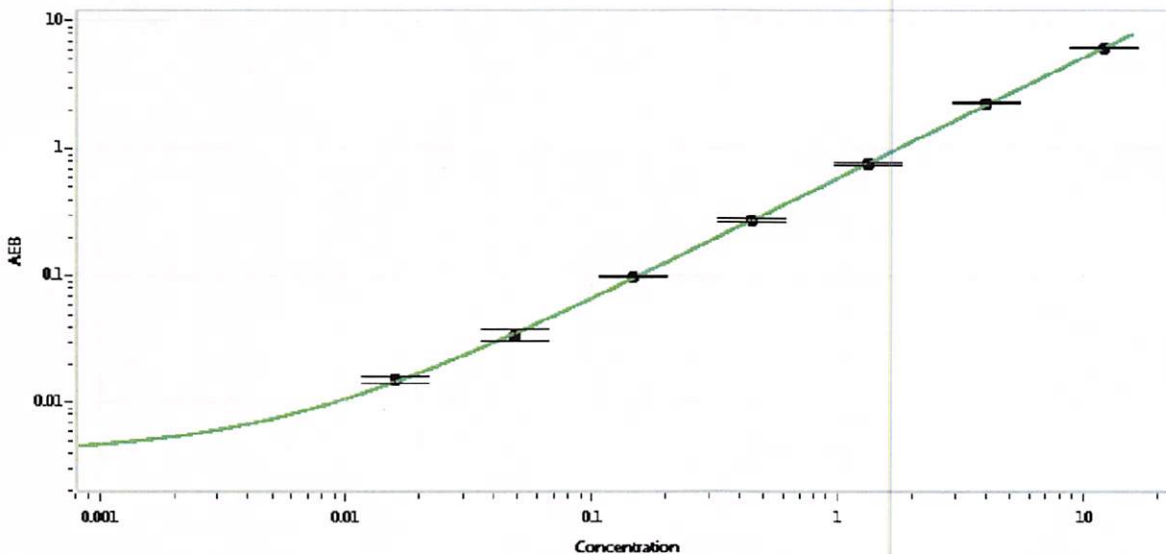
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	121506	0.286	0.229 - 0.344
Control 2	121506	5.14	3.93 - 5.90
Panel 1	033802	0.141	0.109 - 0.242
Panel 2	033802	2.05	1.70 - 2.61
Panel 3	033802	10.6	8.09 - 12.1

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

Calibrator Concentrate Value Assignment (pg/mL)	403
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.016
C	0.049
D	0.148
E	0.444
F	1.33
G	4.00
H	12.0

Review/Approval

Junpil Park

QC Supervisor

[Signature] 1/5/2021

Name

Title

Signature/Date