

Product Number:	101635
Lot Number:	503035
Expiration:	13-Apr-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	101637	110324
Detector Reagent	101636	110508
SBG Reagent	101638	118306
Sample Diluent	101639	110411
Calibrator Diluent	101640	110412
Calibrator Concentrate	100661	111014
RGP Reagent	103159	N/A ¹

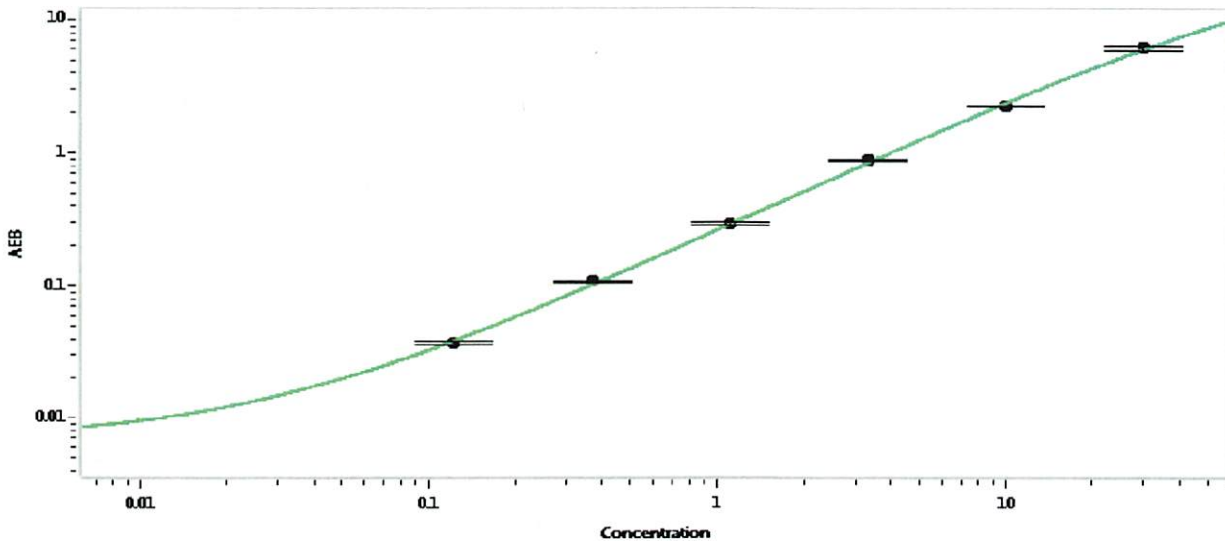
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	111014	0.565	0.450-0.750
Control 2	111014	55.2	45.0-75.0
Panel 1	035301	0.362	0.262-0.456
Panel 2	035301	4.12	3.25-4.88
Panel 3	035301	46.1	35.5-53.2

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

Calibrator Concentrate Value Assignment (pg/mL)	2283
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.123
C	0.370
D	1.11
E	3.33
F	10.0
G	30.0
H	

Review/Approval

Junpil Park
Name

Supervisor, QC
Title

 15 Jul 2021
Signature/Date