

Product Number:	101622
Lot Number:	503029
Expiration:	19-Jul-2022
Platform(s):	HD-1 / HD-X & SR-X



Data below represents results generated on the Simoa™ HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101624	122408
Detector Reagent	101623	122707
SBG Reagent	101625	122708
Sample Diluent	103732	120032
Calibrator Diluent	101627	121307
Calibrator Concentrate	100652	115704
RGP Reagent	103159	N/A ¹

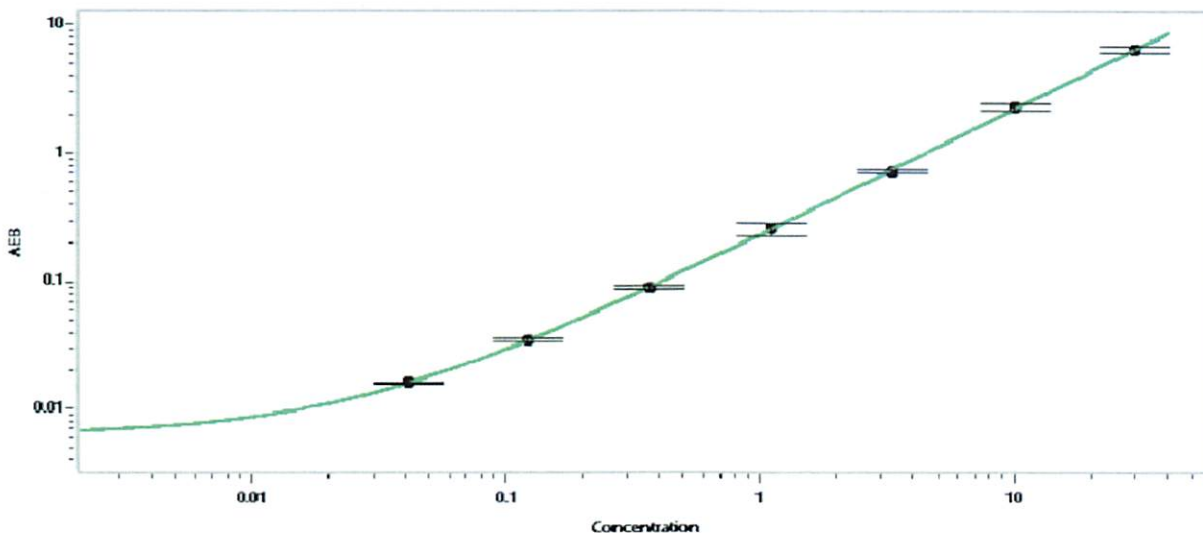
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108406	1.60	1.19-1.99
Control 2	108406	26.6	16.4-28.8
Panel 1	029301	51.2	39.8-59.7
Panel 2	029401	5.86	4.35-6.53
Panel 3	029301	1.56	1.22-1.82

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

Calibrator Concentrate Value Assignment (pg/mL)	657
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Junpil Park

QC Supervisor

 31 Aug 2021

Name

Title

Signature/Date