

Product Number:	101622
Lot Number:	502970
Expiration:	9-May-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	115407
Detector Reagent	101623	116109
SBG Reagent	101625	115602
Sample Diluent	103732	113508
Calibrator Diluent	101627	112912
Calibrator Concentrate	100652	115704
RGP Reagent	103159	N/A ¹

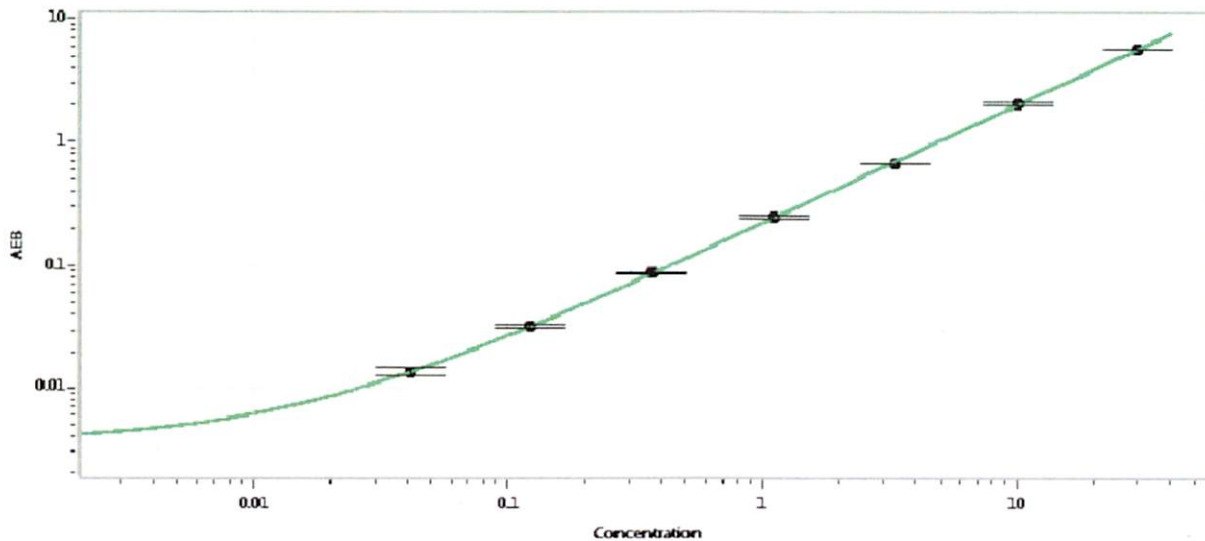
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108406	1.52	1.12-2.04
Control 2	108406	21.1	17.2-25.9
Panel 1	029301	53.9	39.8-59.7
Panel 2	029401	5.65	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
Name

QC Supervisor
Title

 29 Jun 2024
Signature/Date