

Product Number:	101622
Lot Number:	503559
Expiration:	23-May-2023
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	221404
Detector Reagent	101623	221408
SBG Reagent	101625	221409
Sample Diluent	103732	220913
Calibrator Diluent	101627	218621
Calibrator Concentrate	100652	126802
RGP Reagent	103159	N/A ¹

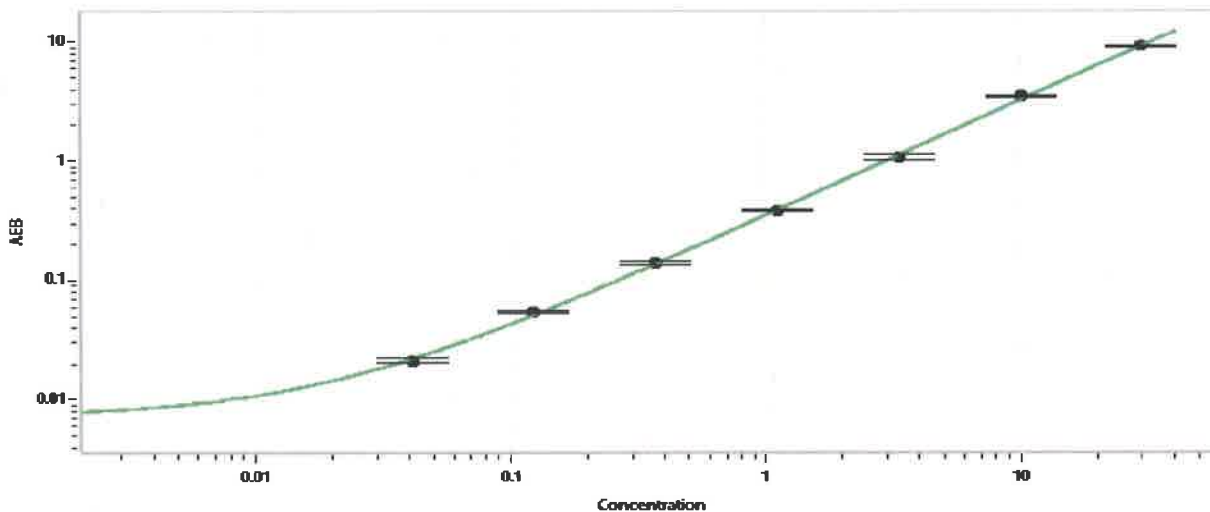
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	108406	1.43	1.23-1.92
Control 2	108406	24.6	17.5-28.9
Panel 1	029301	46.3	39.8-59.7
Panel 2	029401	5.19	4.35-6.53
Panel 3	029301	1.53	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)	541
---	-----

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: Sonia Khillan Name: Sonia Khillan Title: QC Signature/Date: Sonia Khillan / 01SEP2022