

Product Number:	102860
Lot Number:	503958
Expiration:	22-Mar-2024
Platform(s):	HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	102852	330501
Detector Reagent	102854	326804
SBG Reagent	102855	327705
Sample Diluent	102856	326506
Calibrator Diluent	102859	326512
Calibrator Concentrate	102857	331005
RGP Reagent	103159	N/A ¹

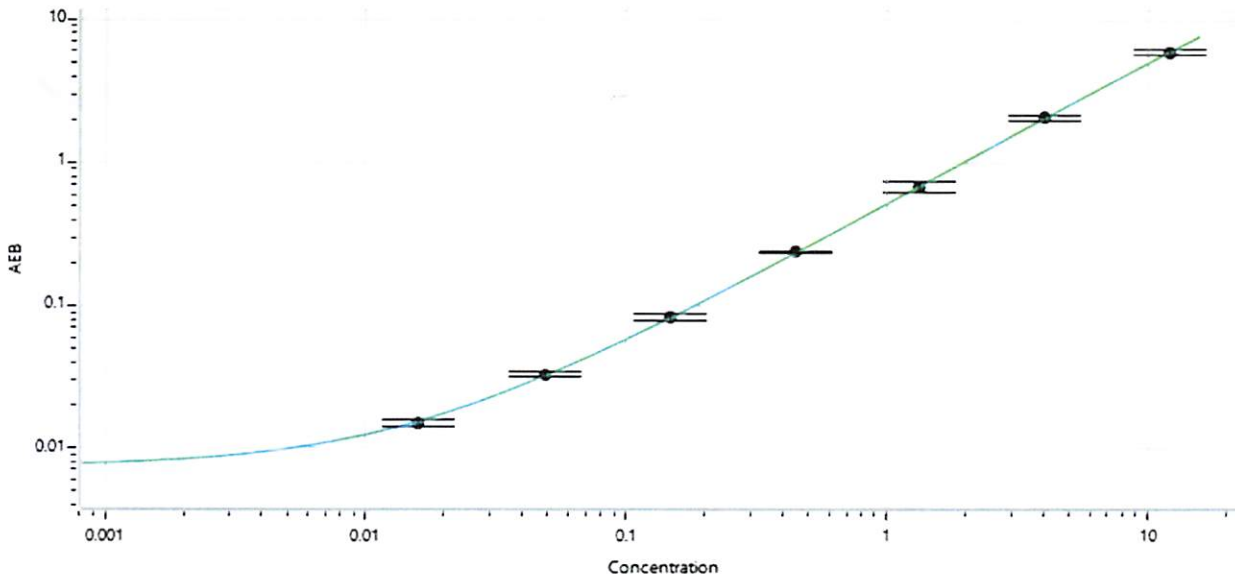
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	331005	0.201	0.180 - 0.300
Control 2	331005	4.21	3.60 - 6.00
Panel 1	231306	0.265	0.223 - 0.335
Panel 2	234108	3.04	2.40 - 3.75
Panel 3	234108	10.4	8.02 - 13.5

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

Calibrator Concentrate Value Assignment (pg/mL)	295
---	-----

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: Sonia Khillan Name, QC Title, [Signature] Signature/Date, 17 NOV 2023 Date