

Product Number:	103277
Lot Number:	503890
Expiration:	17-Jul-2024
Platform(s):	HD-1/ HDX



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

Component	Part Number	Lot Number
Bead Reagent	103265	319811 ¹
Detector Reagent	103268	323404
SBG Reagent	103269	323405
Sample Diluent	103270	320108
Calibrator Diluent	103266	320107
Calibrator Concentrate	103272	323508
RGP Reagent	103159	N/A ²

¹ Control number for Bead reagent lot is 319811A

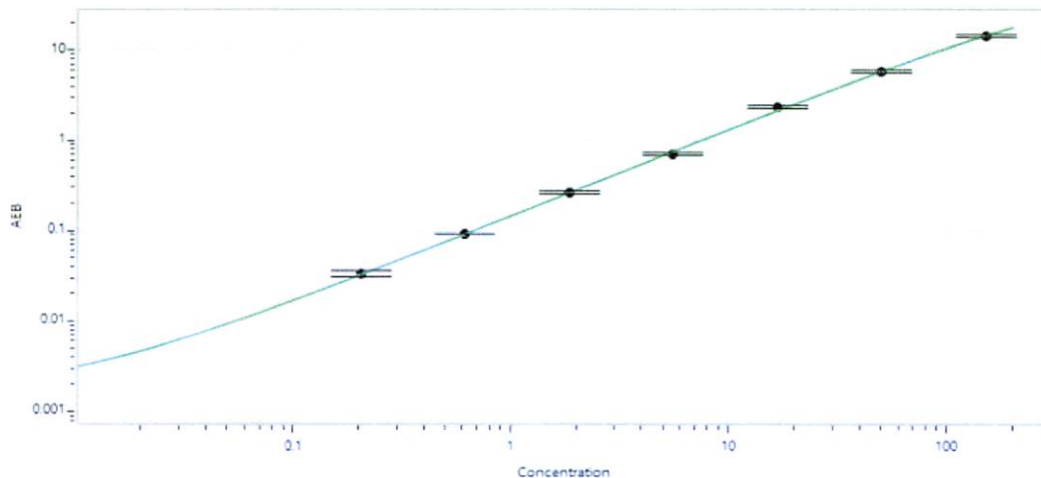
² RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	323002	4.13	3.50 - 5.24
Control 2	323002	121	81.4 - 128
Panel 1	205507	2.09	1.35 - 2.47
Panel 2	205507	33.5	22.4 - 46.5
Panel 3	205507	87.8	61.0 - 112

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

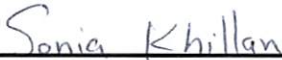

Calibrator Concentrate Value Assignment (pg/mL)	3252
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.206
C	0.617
D	1.85
E	5.56
F	16.7
G	50.0
H	150

Review/Approval

Name: Sonia Khillan Title: QC Signature/Date: Sonia Khillan 11SEP2023