

Product Number:	103277
Lot Number:	503211
Expiration:	7-Nov-2022
Platform(s):	HD-1/ HDX



Data below represents results generated on the Simoa® HD-1 / HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	103265	131104
Detector Reagent	103268	131105
SBG Reagent	103269	131213
Sample Diluent	103270	131206
Calibrator Diluent	103266	131214
Calibrator Concentrate	103272	032905
RGP Reagent	103159	N/A ¹

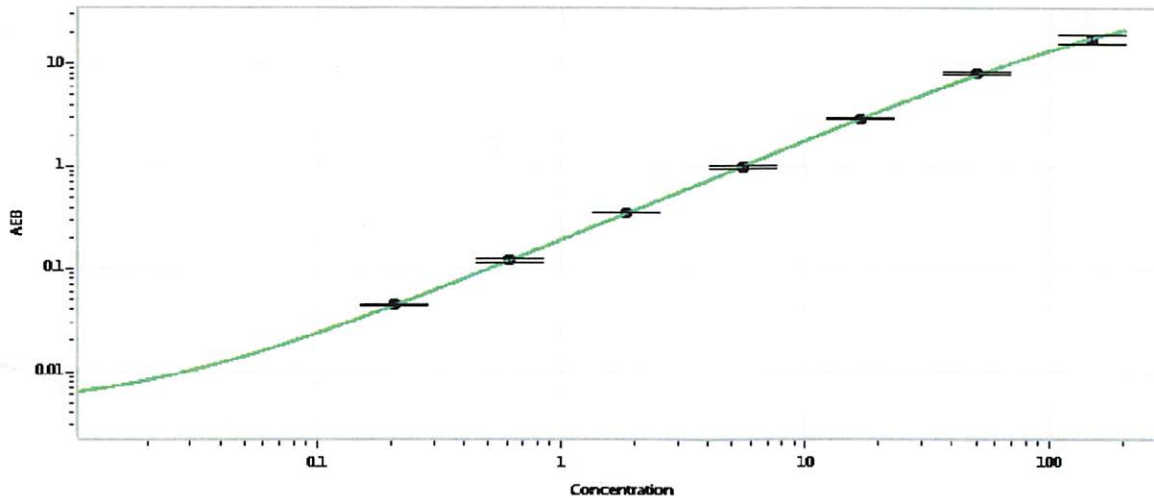
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	203709	3.17	2.76-4.27
Control 2	203709	81.8	67.6-116
Panel 1	205507	1.64	1.30-2.52
Panel 2	205507	32.8	22.8-47.6
Panel 3	205507	78.3	63.4-114

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

Calibrator Concentrate Value Assignment (pg/mL)	5358
---	------

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

Shivani Goel
Name

QC Manager
Title

Shivani Goel 08 April 2022
Signature/Date