

Product Number:	101154
Lot Number:	503331
Expiration:	5-Dec-2022
Platform(s):	HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	101141	202707
Detector Reagent	101143	202808
SBG Reagent	101153	202809
Sample Diluent	103320	133919
Calibrator Diluent	101147	202207
Calibrator Concentrate	101145	105003
RGP Reagent	103159	N/A ¹

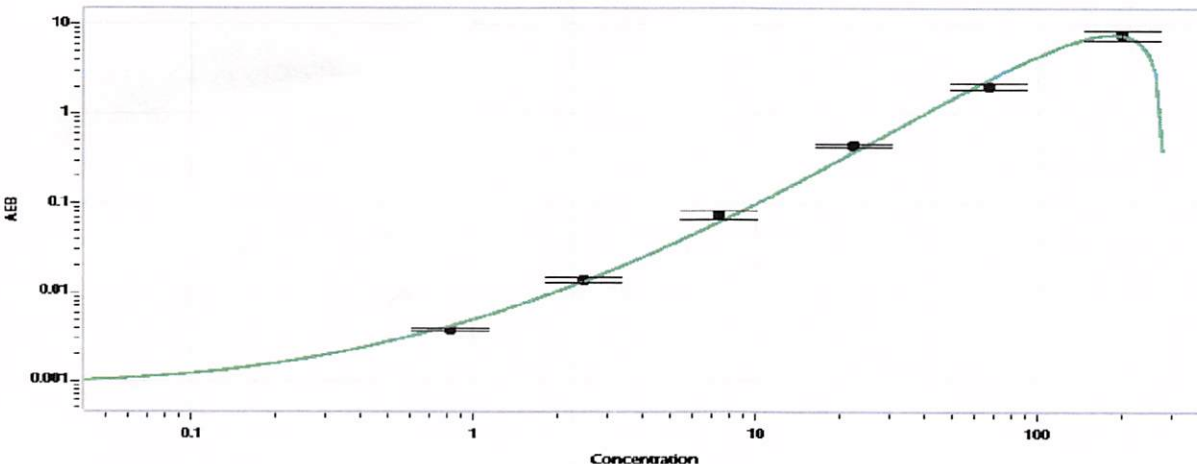
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	918402	37.1	32.3-50.2
Control 2	918402	312	258-456
Panel 1	113008	74.7	63.7-102
Panel 2	113008	41.3	35.4-53.2
Panel 3	115408	247	214-322
Panel 4	113008	182	156-237

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

Calibrator Concentrate Value Assignment (pg/mL)	3841
---	------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.823
C	2.47
D	7.41
E	22.2
F	66.7
G	200

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Soni PAB 27JUL2022
Signature/Date