

Product Number:	101664
Lot Number:	503535
Expiration:	20-Apr-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	101666	218915
Detector Reagent	101665	219202
SBG Reagent	101667	219203
Sample Diluent	101668	219204
Calibrator Diluent	101669	219208
Calibrator Concentrate	103217	219209
RGP Reagent	103159	N/A ¹

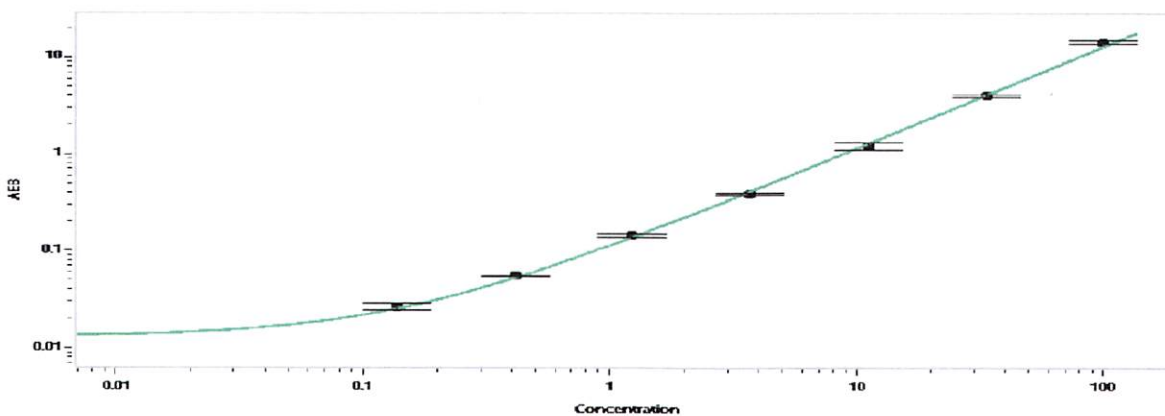
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	201802	1.10	0.681-1.45
Control 2	201802	64.1	45.8-78.9
Panel 1	202604	8.17	5.87-9.46
Panel 2	202604	47.6	31.3-58.5
Panel 3	202604	123	81.7-158

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

Calibrator Concentrate Value Assignment (pg/mL)	1323
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval

Sonia Khillan @C Manager *Sonia Khillan* / 29 JUL 2022
Name Title Signature/Date