

Product Number:	101132
Lot Number:	503330
Expiration:	26-Nov-2022
Platform(s):	HD-1/HD-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101119	203812
Detector Reagent	101121	203902
SBG Reagent	101131	203903
Sample Diluent	103319	133922
Calibrator Diluent	101125	202208
Calibrator Concentrate	101123	203904
RGP Reagent	103159	N/A ¹

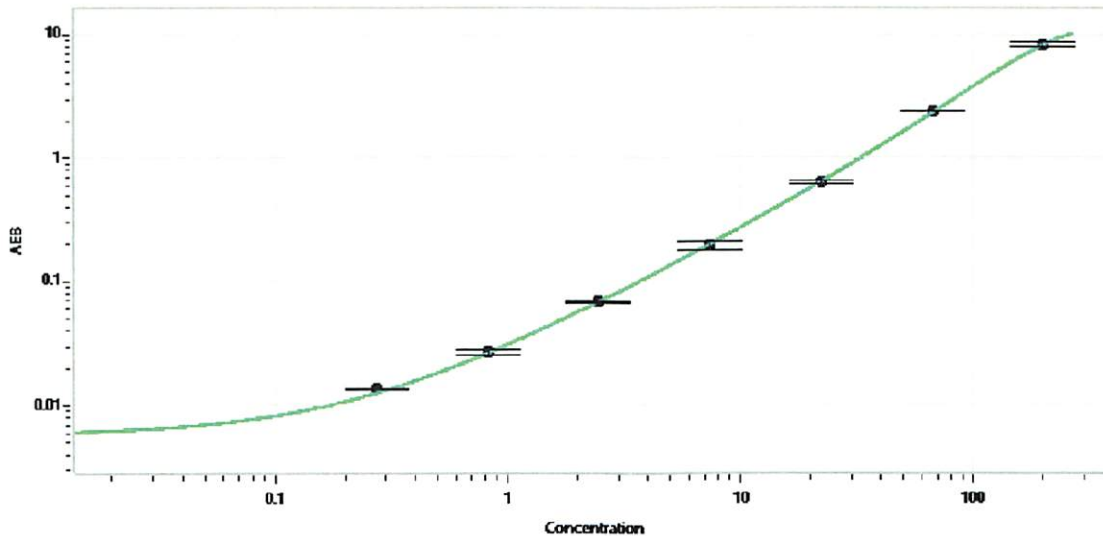
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	130713	7.17	6.39-9.89
Control 2	130713	241	163-298
Panel 1	106201	124	88.5-213
Panel 2	106201	71.8	56.2-123
Panel 3	106201	227	185-352
Panel 4	106201	226	189-296

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

Calibrator Concentrate Value Assignment (pg/mL)	5081
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval: Shivani Goel QC Shivani Goel 01Mar2022
Name Title Signature/Date