

Product Number:	101664
Lot Number:	503004
Expiration:	15-Jun-2022
Platform(s):	HD-1/HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101666	116606
Detector Reagent	101665	117003
SBG Reagent	101667	117004
Sample Diluent	101668	117005
Calibrator Diluent	101669	117001
Calibrator Concentrate	103217	202410
RGP Reagent	103159	N/A ¹

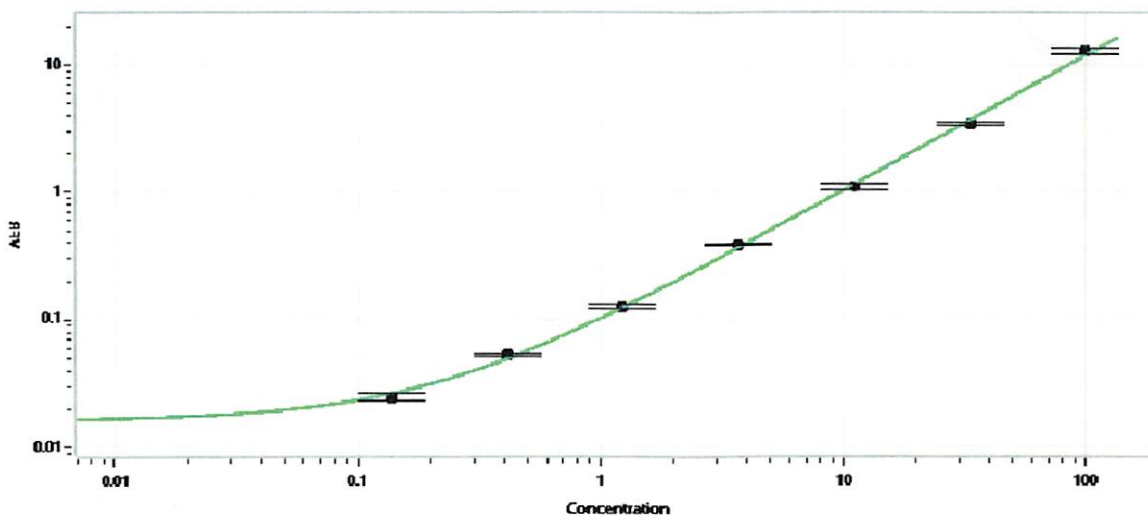
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	201802	0.876	0.704-1.30
Control 2	201802	65.5	47.0-70.9
Panel 1	202604	8.19	5.73-8.59
Panel 2	202604	45.6	33.2-49.8
Panel 3	202604	108	89.6-134
Panel 4	N/A	N/A	N/A

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

Calibrator Concentrate Value Assignment (pg/mL)	1417
---	------

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

Shivani Goel
Name

QC Manager
Title

Shivani Goel 11 Feb 2022
Signature/Date