

Product Number:	101580
Lot Number:	502507
Expiration:	2-Jun-2021
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	101582	023303
Detector Reagent	101581	023401
SBG Reagent	101583	023402
Sample Diluent	103201	015425
Calibrator Diluent	101585	023403
Calibrator Concentrate	100655	932607
RGP Reagent	103159	N/A ¹

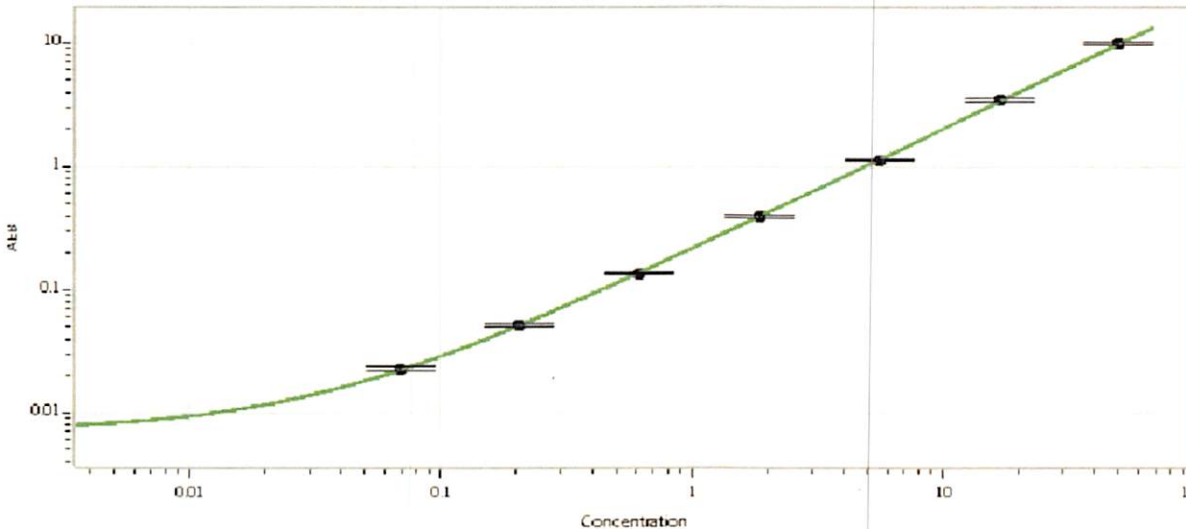
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-042817-01	2.65	2.08-3.12
Control 2	MF-051617-02	48.0	35.1-52.7
Panel 1	MF-051017-03	29.0	19.9-32.1
Panel 2	MF-032217-11	1.70	1.33-2.23
Panel 3	MF-033018-06	2.65	1.68-3.50

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

Calibrator Concentrate Value Assignment (pg/mL)	1311
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.069
C	0.206
D	0.617
E	1.85
F	5.56
G	16.7
H	50.0

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 03 Sep 2020
	Name	Title	Signature/Date