

Product Number:	101212
Lot Number:	501411
Expiration:	28-Feb-2019
Platform(s):	HD-1



Component	Part Number	Lot Number
Bead Reagent	101198	824306
Detector Reagent	101200	824007
SBG Reagent	101211	824008
Sample Diluent	101209	824109
Calibrator Diluent	101204	824110
Calibrator Concentrate	101202	72780
RGP Reagent	103159	N/A ¹

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

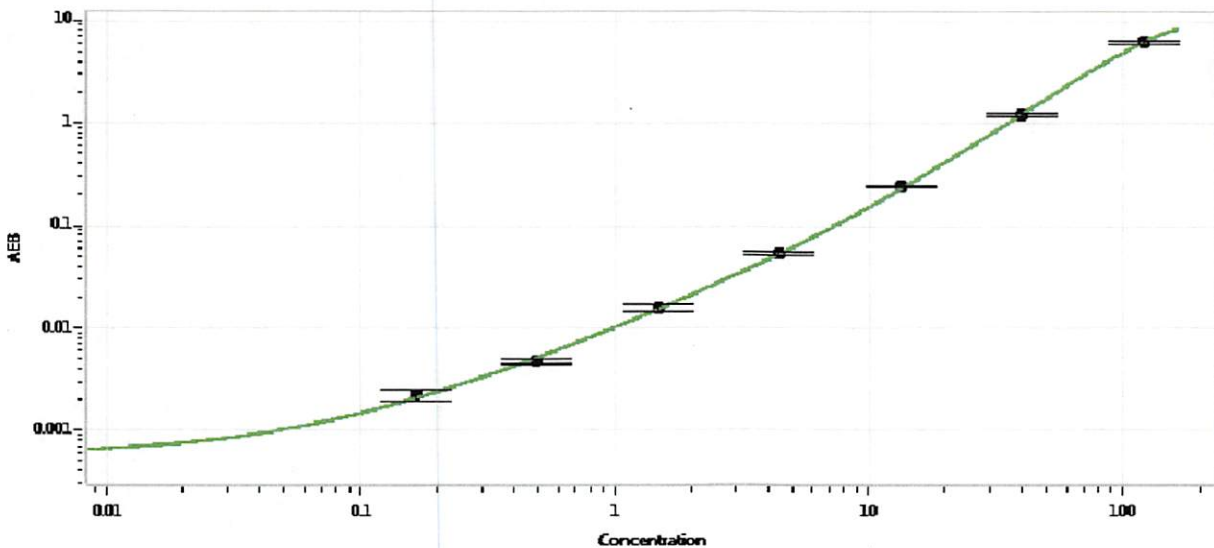
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 2	828802	68.7	57.4-86.2
Control 3	828802	531	455-683
Panel 1	MF-052318-05	190	144-216
Panel 2	MF-052318-06	46.3	353-52.9
Panel 3	MF-052318-07	97.9	73.9-111

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

Calibrator Concentrate Value Assignment (pg/mL)	2,037
---	-------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.165
C	0.494
D	1.48
E	4.44
F	13.3
G	40.0
H	120

Review/Approval

Sivasankari Murugar QC Supervisor
Name Title

[Signature] 19 OCT 2018
Signature/Date