

Product Number:	101154
Lot Number:	501065
Expiration:	29-May-2018
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	101141	73521
Detector Reagent	101143	73522
SBG Reagent	101153	73523
Sample Diluent	101151	73500
Calibrator Diluent	101147	73334
Calibrator Concentrate	101145	73524
RGP Reagent	103159	N/A ¹

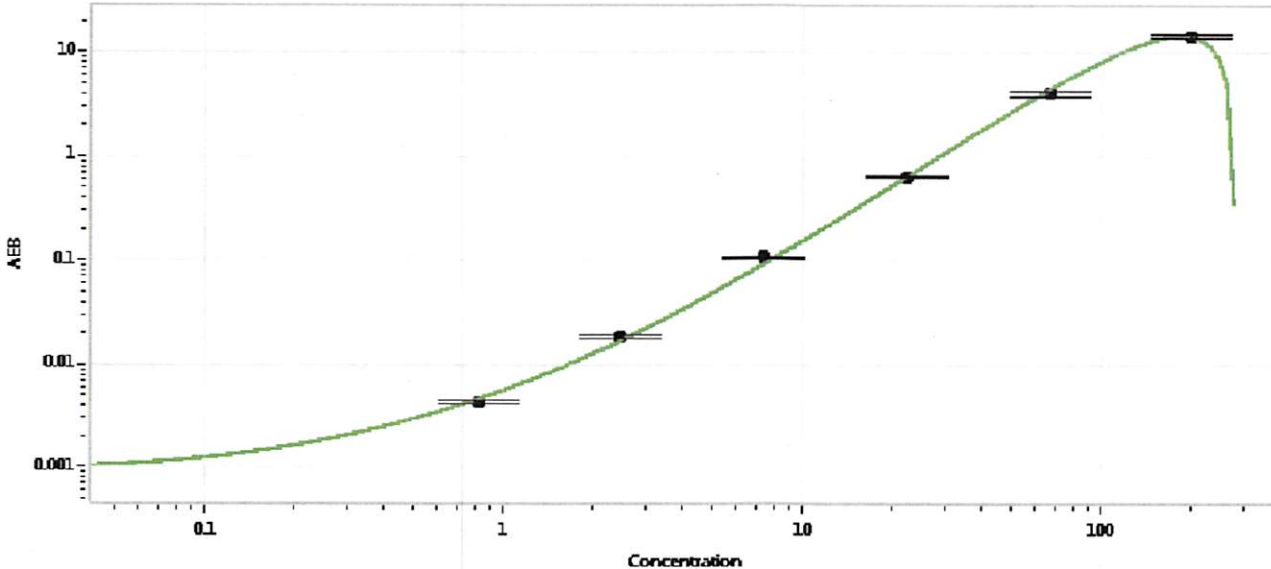
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	41952	58.6	36.7-80.4
Control 2	41952	339	208-456
Panel 1	1816	88.6	42.6-115
Panel 2	1712	91.9	47.3-104
Panel 3	2004	132	78.3-165

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

Calibrator Concentrate Value Assignment (pg/mL)	14,199
---	--------

Calibration Curve



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

Review/Approval: Sivasankari Murugan Name, QC Supervisor Title, [Signature] Signature/Date, 31 Jan 2018 Date