

Product Number:	101154
Lot Number:	501809
Expiration:	18-Dec-2019
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	101141	912108
Detector Reagent	101143	909107
SBG Reagent	101153	909106
Sample Diluent	103320	907312
Calibrator Diluent	101147	908816
Calibrator Concentrate	101145	73524
RGP Reagent	103159	N/A ¹

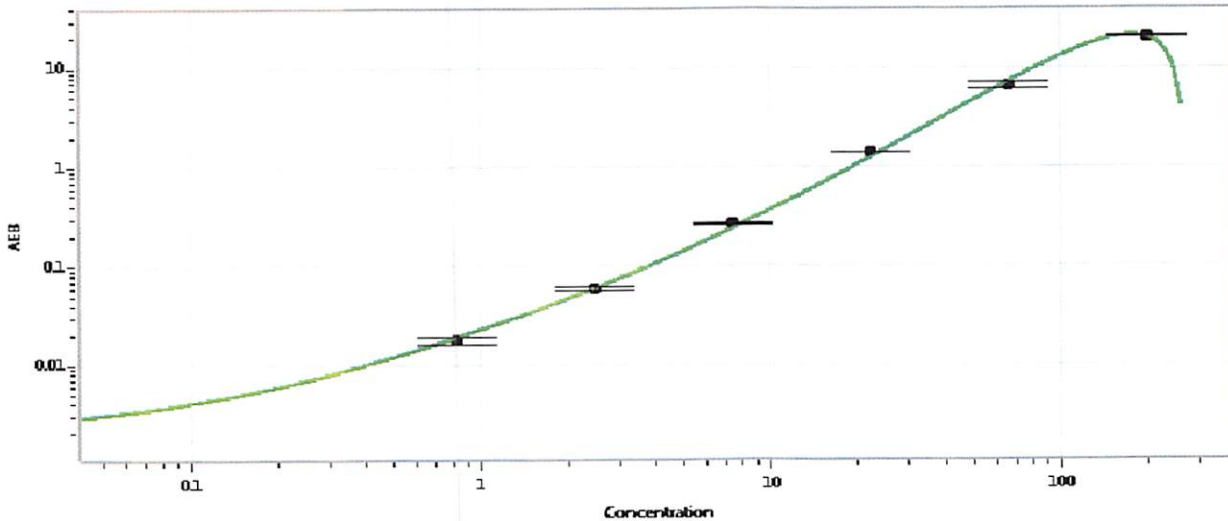
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	41952	68.1	51.8-77.6
Control 2	41952	362	269-404
Panel 1	909901	92.8	68.6-103
Panel 2	909902	91.7	66.7-100
Panel 3	909903	263	186-278
Panel 4	909904	221	180-270

* 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

Patrizia Stadler QC Assoc. Director
Name Title

P. Stoolle 07MAY2019
Signature/Date