

Product Number:	101622
Lot Number:	502058
Expiration:	27-Feb-2020
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	101624	923803
Detector Reagent	101623	923110
SBG Reagent	101625	923204
Sample Diluent	103203	912023
Calibrator Diluent	101627	905824
Calibrator Concentrate	100652	906001
RGP Reagent	103159	N/A ¹

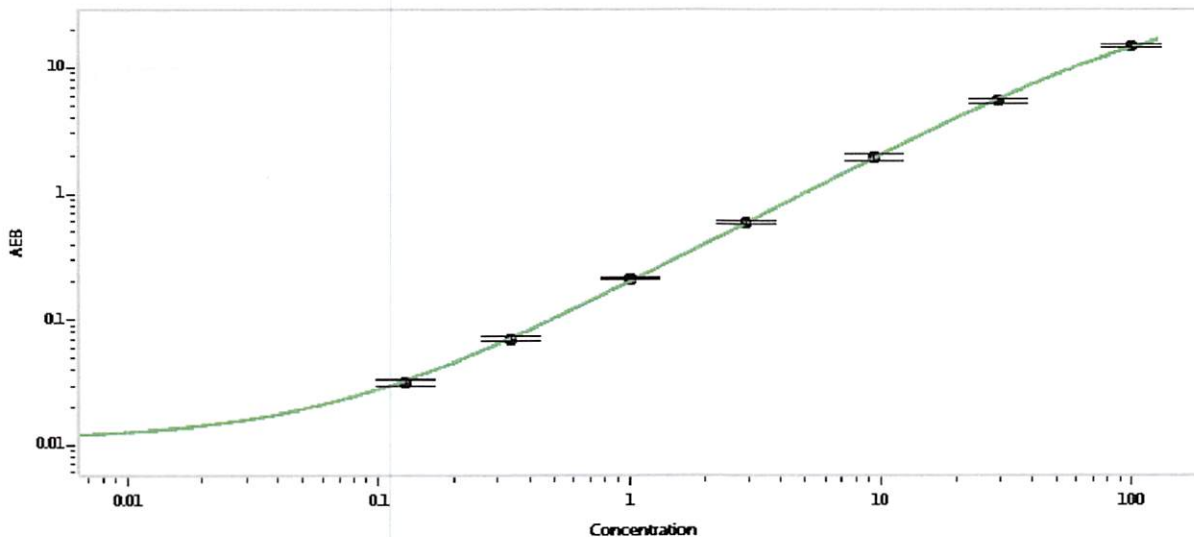
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	901702	1.28	0.934-1.40
Control 2	901702	20.9	15.1-22.7
Panel 1	MF-050918-01	60.7	41.6-67.5
Panel 2	MF-050918-02	8.28	5.03-8.84
Panel 3	MF-050918-03	1.15	0.644-1.24

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

Calibrator Concentrate Value Assignment (pg/mL)	864
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	10 Oct 2019
	Name	Title	Signature/Date	