

Product Number:	101622
Lot Number:	502358
Expiration:	07Oct2020
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	928906
Detector Reagent	101623	928809
SBG Reagent	101625	928802
Sample Diluent	103203	928021
Calibrator Diluent	101627	928013
Calibrator Concentrate	100652	006408
RGP Reagent	103159	N/A ¹

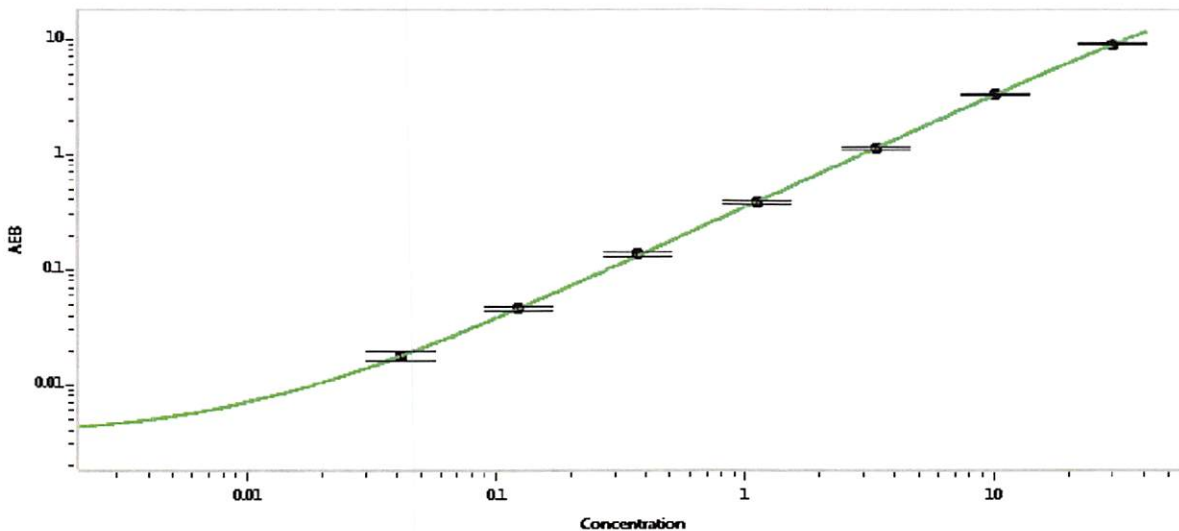
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	901702	1.05	0.944-1.42
Control 2	901702	17.2	15.4-23.1
Panel 1	MF-050918-01	49.9	41.6-67.5
Panel 2	MF-050918-02	6.99	5.03-8.84
Panel 3	MF-050918-03	1.09	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)	675
---	-----

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

Shivani Goel

QC Manager

Shivani Goel 11 May 2020

Name

Title

Signature/Date