

Product Number:	101622
Lot Number:	501828
Expiration:	30-Mar-2020
Platform(s):	HD-1 & SR-X

Component	Part Number	Lot Number
Bead Reagent	101624	912907
Detector Reagent	101623	913602
SBG Reagent	101625	913003
Sample Diluent	103203	908910
Calibrator Diluent	101627	912015
Calibrator Concentrate	100652	913004
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific



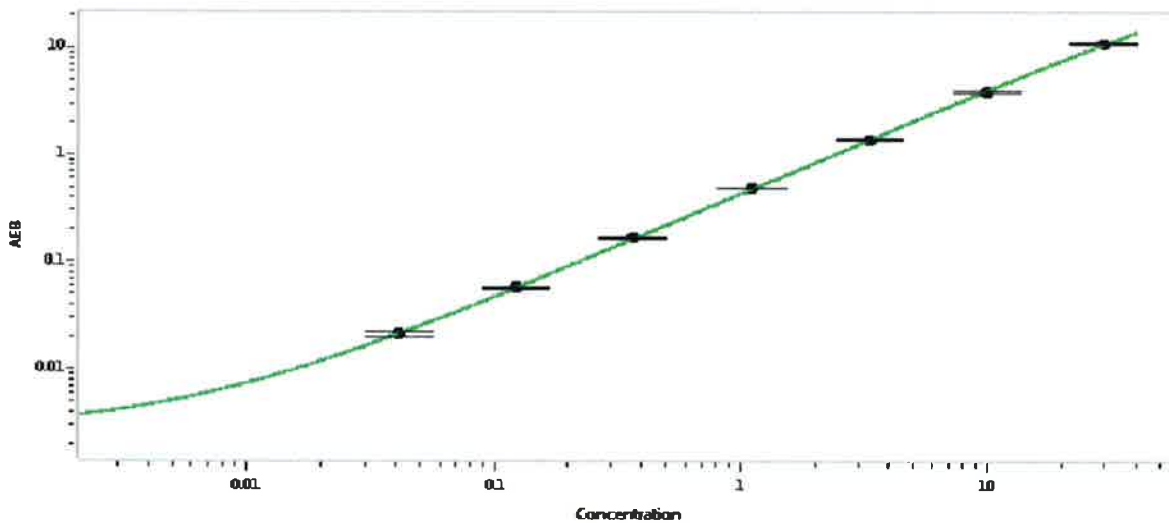
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	901702	1.14	0.923-1.38
Control 2	901702	19.5	14.9-22.4
Panel 1	MF-050918-01	55.1	41.6-67.5
Panel 2	MF-050918-02	7.00	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)	519
---	-----

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>	17 Jan 2019
	Name	Title	Signature/Date	