

Product Number:	101478
Lot Number:	501811
Expiration:	13-Feb-2020
Platform(s):	HD-1 & SR-X



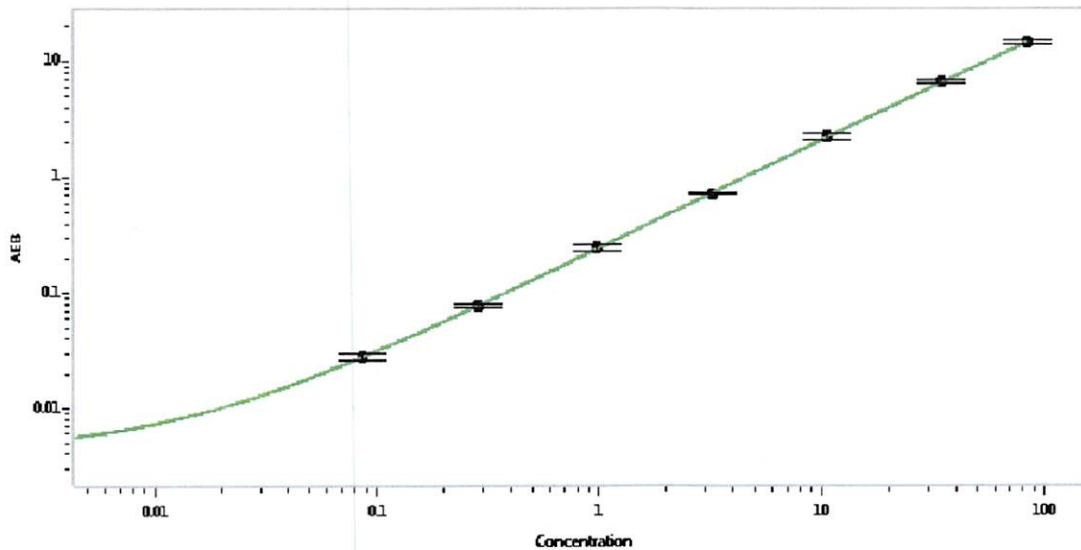
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101480	912205
Detector Reagent	101481	912207
SBG Reagent	101482	912208
Sample Diluent	101483	904417
Calibrators	100758-100765	911302/912101
Controls	100768-100769	912102
RGP Reagent	103159	N/A <sup>1</sup>

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

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	912102	2.73	2.29-3.60	
Control 2	912102	69.5	54.3-81.5	
Panel 1	MF-051017-04	0.575	0.508-0.855	
Panel 2	MF-042117-06	4.36	3.89-6.07	
Panel 3	MF-042117-07	58.7	45.5-73.7	

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.086
C	0.288
D	0.980
E	3.23
F	10.6
G	34.6
H	82.9

QC  
 Review/Approval Patrizia Stadler Assoc. Director P. Stodler 10 MAY 2019  
 Name Title Signature/Date