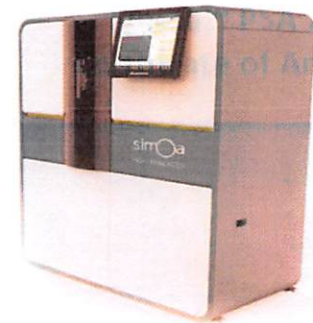


Product Number:	101478
Lot Number:	501821
Expiration:	19-Oct-2019
Platform(s):	HD-1 & SR-X



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

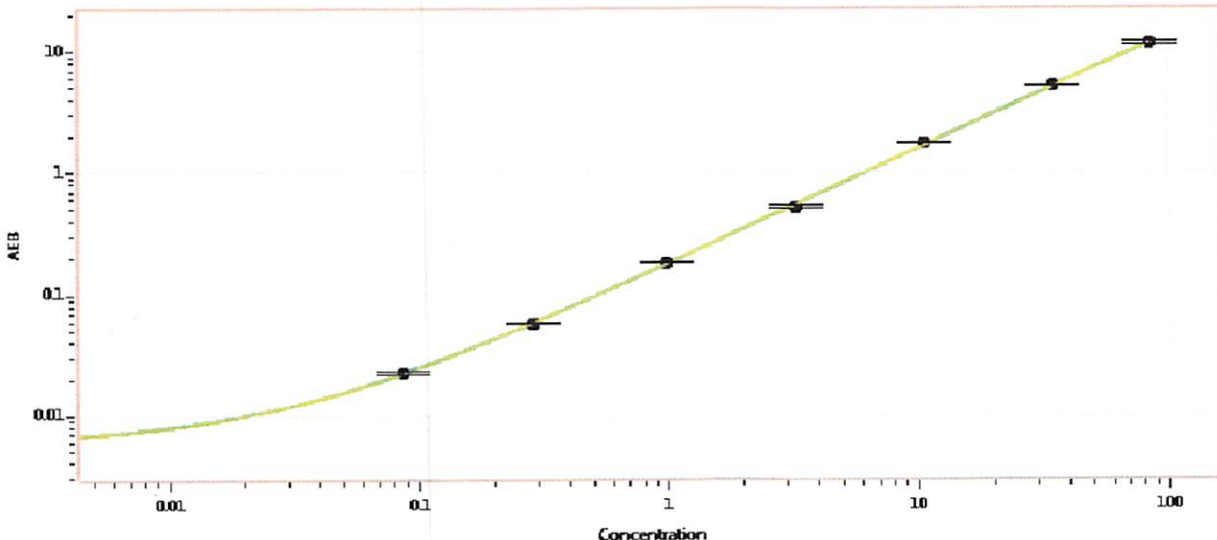
Component	Part Number	Lot Number
Bead Reagent	101480	831217
Detector Reagent	101481	829812
SBG Reagent	101482	829211
Sample Diluent	101483	829609
Calibrators	100758-100765	911301/912101
Controls	100768 & 100769	912102
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	912102	3.02	2.29-3.60
Control 2	912102	69.4	54.3-81.5
Panel 1	MF-051017-04	0.609	0.508-0.855
Panel 2	MF-042117-06	4.32	3.89-6.07
Panel 3	MF-042117-07	52.6	45.5-73.7

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

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

Review/Approval

Patrizia Stadler
Name

QC Assoc. Director
Title

P. Stadler 10/17/19
Signature/Date