

Product Number: 103337
Lot Number: 501655
Expiration: 24-Oct-2019
Platform(s): HD-1 & SR-X



Component	Item Number	Lot Number
Bead Reagent	103334	833909
Detector Reagent	101091	833904
SBG Reagent	101101	833902
Sample Diluent	103336	933903
Calibrator Diluent	101095	829715
Calibrator Concentrate	101093	828403
RGP Reagent	103159	N/A ¹

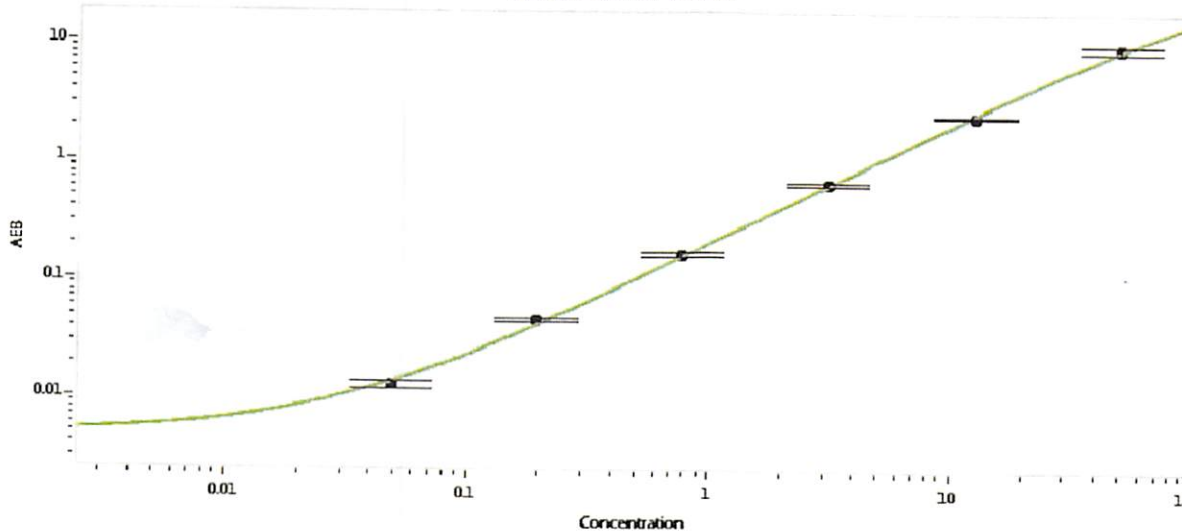
Data below represents results generated on the Simoa™ HD-1 Analyzer.

¹ 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	MF-092215-03	9.06	6.93-10.4	
Control 2	MF-092215-03	87.1	68.9-103	
Panel 1	MF-033018-03	3.35	2.56-3.84	
Panel 2	MF-033018-04	5.96	4.86-7.29	
Panel 3	MF-033018-05	62.1	46.9-70.4	

Calibrator Concentrate Value Assignment (pg/mL) 1,291

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.049
C	0.195
D	0.781
E	3.13
F	12.5
G	50.0

Review/Approval

Patrizia Stadler QC Assoc. Director

P. Stadler 08Mar2019

Name

Title

Signature/Date