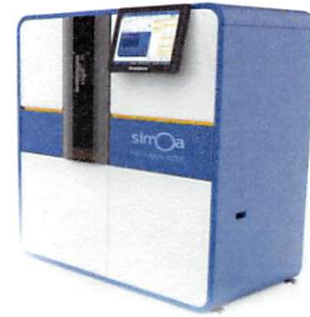


Product Number:	101588
Lot Number:	501228
Expiration:	21-Sep-2018
Platform(s):	HD-1



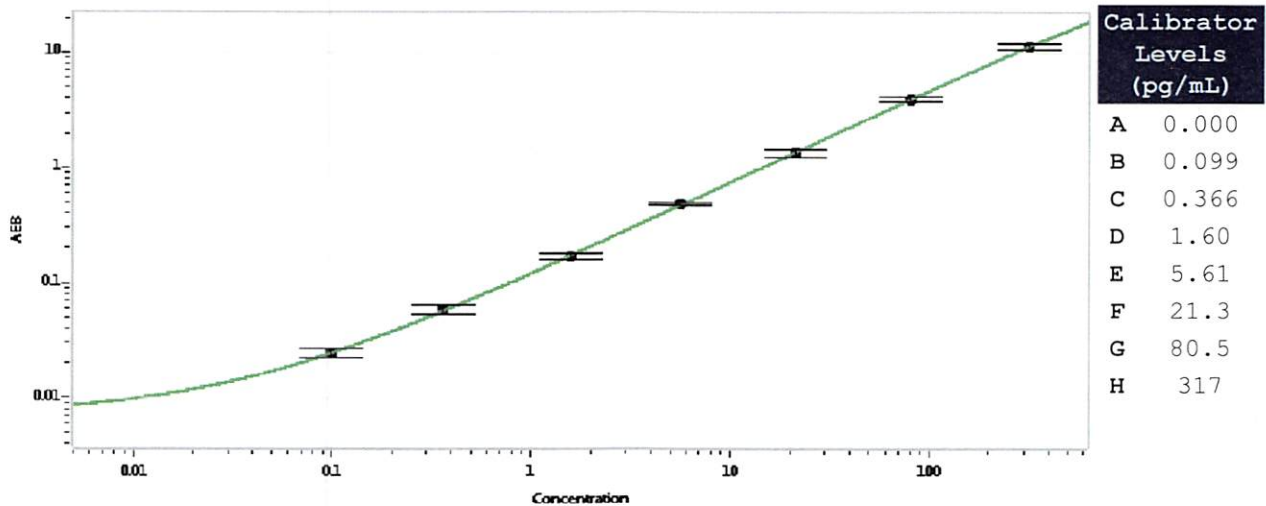
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101590	81137
Detector Reagent	101589	80804
SBG Reagent	101591	81150
Sample Diluent	101592	80801
Calibrators	101110	80330
Controls	100551-100552	80331
RGP Reagent	103159	N/A ¹

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	80331	0.208	0.117-0.295	
Control 2	80331	156	121-200	
Panel 1	MF-021716-05	2.23	1.66-2.50	
Panel 2	MF-021716-06	4.82	4.18-6.27	
Panel 3	MF-021716-07	112	95.7-144	

Calibration Curve



Review/Approval Sivasankari Murugan QC Supervisor
Name Title

S. Murugan 27/09/2018
Signature/Date