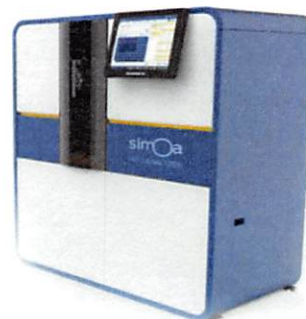


<b>Product Number:</b>	101588
<b>Lot Number:</b>	501274
<b>Expiration:</b>	29-Nov-2018
<b>Platform(s):</b>	HD-1



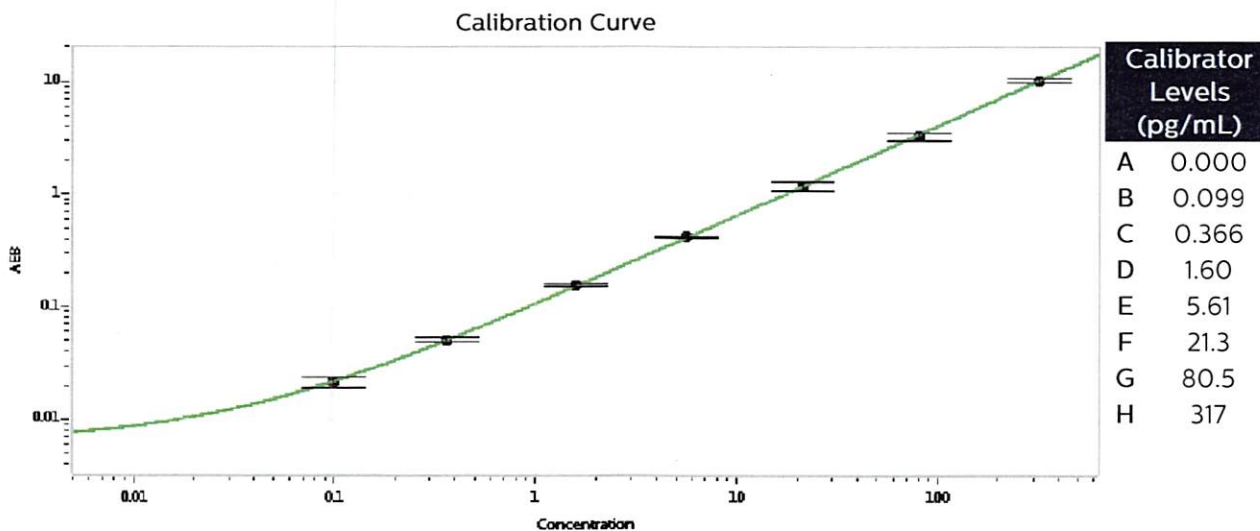
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101590	81513
Detector Reagent	101589	81502
SBG Reagent	101591	81514
Sample Diluent	101592	81493
Calibrators	101110	80330
Controls	100551-100552	80331
RGP Reagent	103159	N/A <sup>1</sup>

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

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80331	0.202	0.114-0.297
Control 2	80331	152	120-199
Panel 1	MF-021716-05	2.08	1.66-2.50
Panel 2	MF-021716-06	4.87	4.18-6.27
Panel 3	MF-021716-07	117	95.7-144

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



Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date  
08 Jun 2018