

Product Number:	100860
Lot Number:	501151
Expiration:	22-Aug-2018
Platform(s):	HD-1



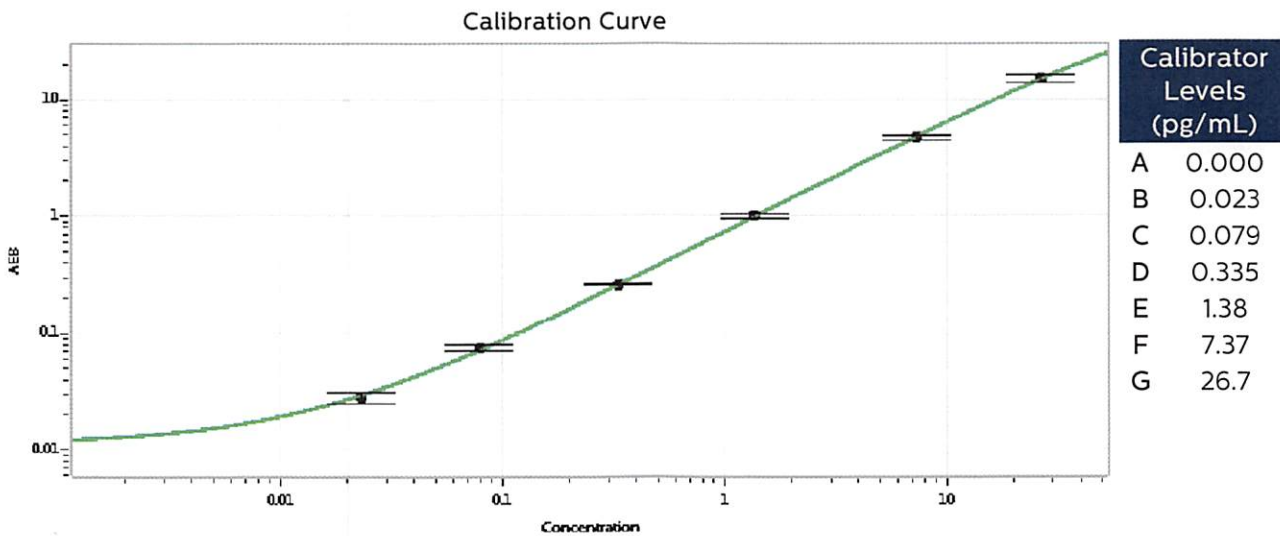
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	100865	80542
Detector Reagent	100862	80532
SBG Reagent	100868	80541
Sample Diluent	100870	80530
Calibrators	101111	72500/72790
Controls	100875 & 100876	71230
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	71230	14.7	10.0-15.4
Control 2	71230	0.207	0.167-0.250
Panel 1	MF-011916-05	1.95	1.32-2.35
Panel 2	MF-011916-06	15.0	6.74-15.7

\* 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 28 FEB 2018