

Product Number:	103186
Lot Number:	501450
Expiration:	29-Aug-2019
Platform(s):	HD-1 & SR-X



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

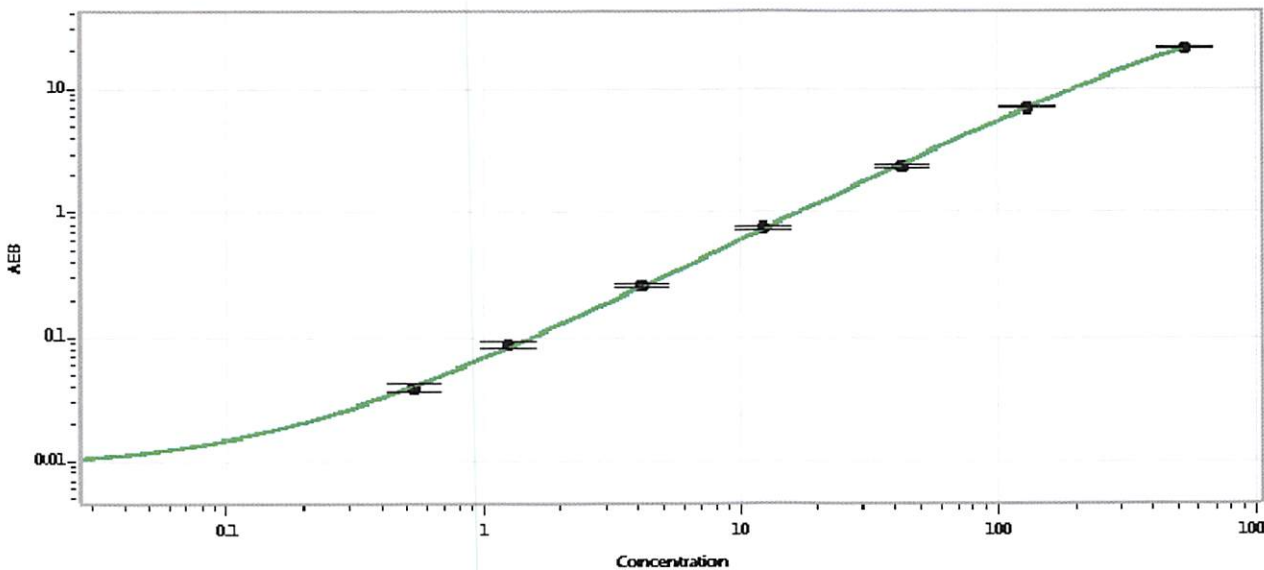
Component	Part Number	Lot Number
Bead Reagent	102246	824903
Detector Reagent	102248	824203
SBG Reagent	102250	824101
Sample Diluent	102252	826205
Reference Calibrators	102255	827701   828307
Controls	102256 - 102257	827702
RGP Reagent	103159	N/A <sup>1</sup>

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

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	827702	3.68	2.83-4.65
Control 2	827702	149	120-180
Panel 1-A	823301	5.59	3.99-6.68
Panel 1-B	823302	6.37	4.74-7.59
Panel 2	823601	91.8	56.9-100
Panel 3	823602	336	226-402

\* 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	0.538
C	1.25
D	4.13
E	12.2
F	42.2
G	131
H	530

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date  
16 Oct 2018