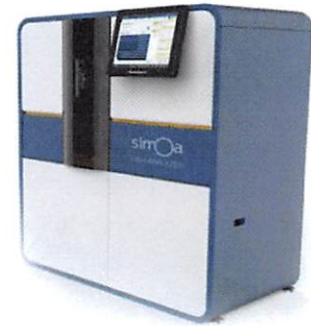


Product Number:	103186
Lot Number:	501484
Expiration:	29-Aug-2019
Platform(s):	HD-1 & SR-X



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

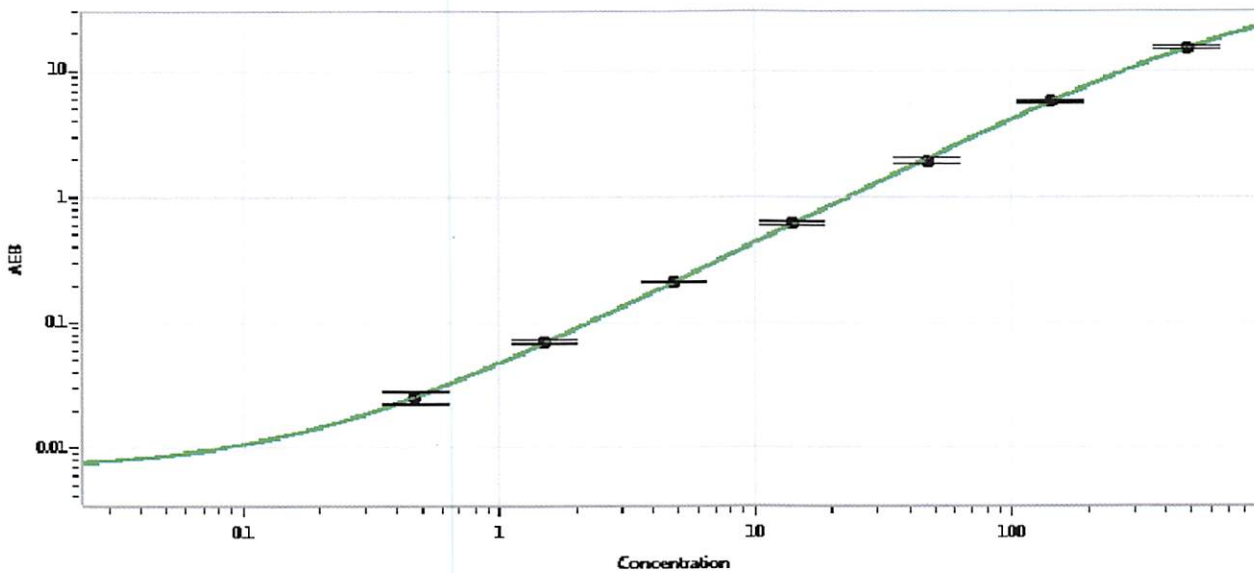
Component	Part Number	Lot Number
Bead Reagent	102246	824903
Detector Reagent	102248	824203
SBG Reagent	102250	824101
Sample Diluent	102252	826205
Reference Calibrators	102255	830302
Controls	102256 - 102257	830201
RGP Reagent	103159	N/A <sup>1</sup>

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

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	830201	4.32	3.20-5.08
Control 2	830201	159	122-183
Panel 1-A	823301	5.41	3.99-6.68
Panel 1-B	823302	6.56	4.74-7.59
Panel 2	823601	90.3	56.9-100
Panel 3	823602	337	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	0.468
C	1.49
D	4.81
E	13.8
F	46.8
G	142
H	482

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]* 08NOV2018  
Signature/Date