

Product Number:	102732
Lot Number:	501605
Expiration:	19-Jan-2020
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	102722	902209
Detector Reagent	102724	901902
SBG Reagent	102726	901901
Sample Diluent	102728	901905
Calibrator Diluent	102731	901906
Calibrator Concentrate	100658	833101
RGP Reagent	103159	N/A <sup>1</sup>

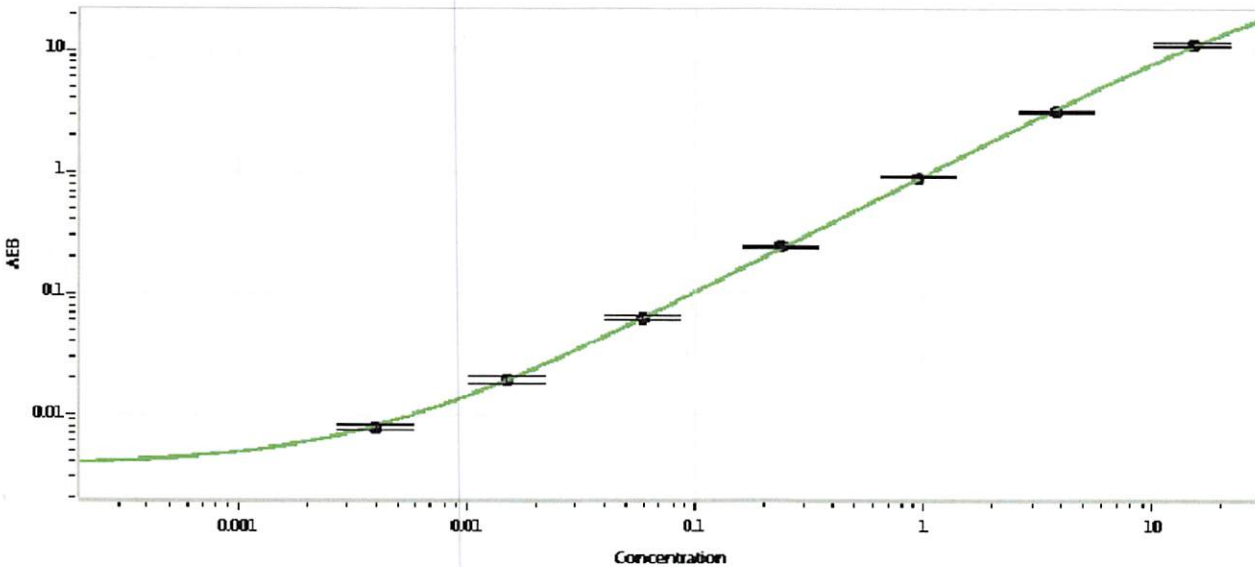
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-061217-01	0.173	0.134-0.289
Control 2	MF-061217-01	3.59	2.44-6.35
Panel 1	MF-061317-03	0.042	0.037-0.088
Panel 2	MF-061317-04	0.260	0.191-0.490
Panel 3	MF-061317-05	4.12	3.18-7.89

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

Calibrator Concentrate Value Assignment (pg/mL)	202
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.004
C	0.015
D	0.059
E	0.234
F	0.938
G	3.75
H	15.0

Review/Approval	Patrizia Stadler	Associate Director, QC	<i>P. Stadler</i>	23 Jan 2019
	Name	Title	Signature/Date	