

Product Number:	102329
Lot Number:	502506
Expiration:	2-Jun-2021
Platform(s):	HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	102322	023107
Detector Reagent	102324	023301
SBG Reagent	102325	023302
Sample Diluent	102326	015424
Calibrator Diluent	102327	020424
Calibrator Concentrate	100849	006603
RGP Reagent	103159	N/A <sup>1</sup>

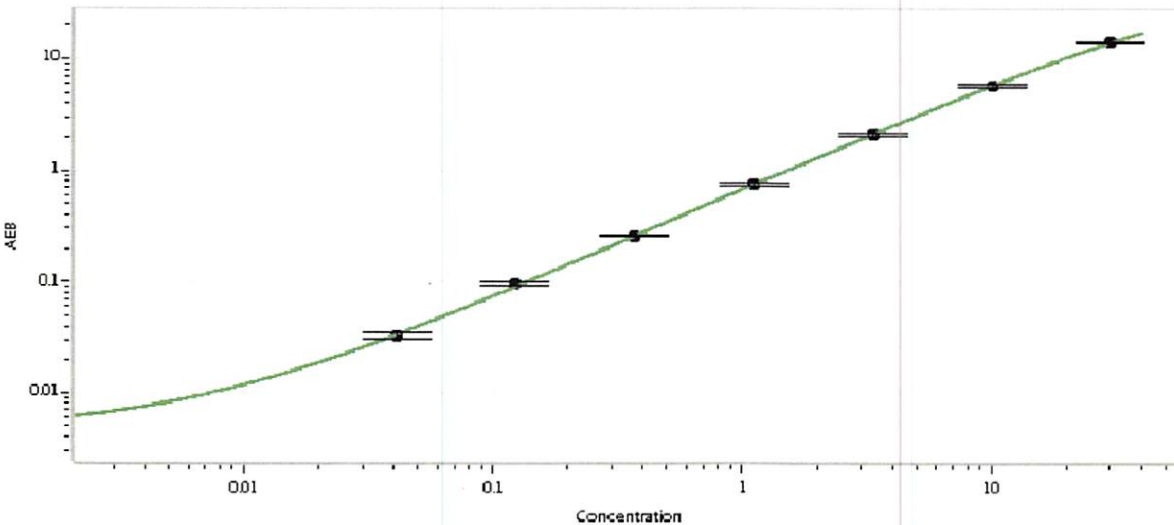
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-092716-01	0.239	0.139-0.285
Control 2	MF-092716-01	30.2	19.4-37.3
Panel 1	MF-092116-03	0.098	0.065-0.135
Panel 2	MF-092116-04	0.093	0.065-0.135
Panel 3	MF-092116-05	19.1	10.5-25.3

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

Calibrator Concentrate Value Assignment (pg/mL)	4586
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 03 Sep 2020

Name

Title

Signature/Date