

Product Number:	100988
Lot Number:	503169
Expiration:	24-Apr-2022
Platform(s):	HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	100974	113704
Detector Reagent	100976	127512
SBG Reagent	100987	113711
Sample Diluent	103734	113507
Calibrator Diluent	100980	111409
Calibrator Concentrate	100978	201206
RGP Reagent	103159	N/A <sup>1</sup>

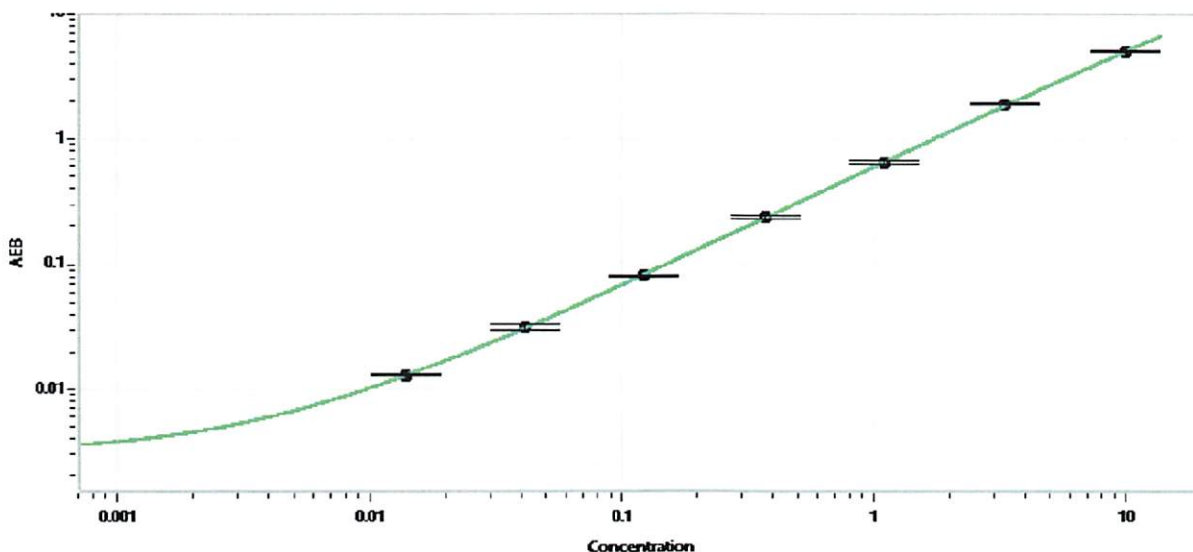
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108314	0.166	0.013-0.293
Control 2	108314	13.4	2.53-20.9
Panel 1	030001	0.140	0.034-0.278
Panel 3	030001	30.6	5.90-49.1
Panel 4	030001	0.163	0.046-0.262

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

Calibrator Concentrate Value Assignment (pg/mL)	136
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.014
C	0.041
D	0.123
E	0.370
F	1.11
G	3.33
H	10.0

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 02 Feb 2022

Name

Title

Signature/Date