

Product Number:	100988
Lot Number:	503283
Expiration:	2-Oct-2022
Platform(s):	HD-1/HD-X



Component	Part Number	Lot Number
Bead Reagent	100974	204505
Detector Reagent	100976	127512
SBG Reagent	100987	204509
Sample Diluent	103734	204511
Calibrator Diluent	100980	133925
Calibrator Concentrate	100978	201206
RGP Reagent	103159	N/A <sup>1</sup>

Data below represents results generated on the Simoa® HD-1 / HD-X Analyzer.

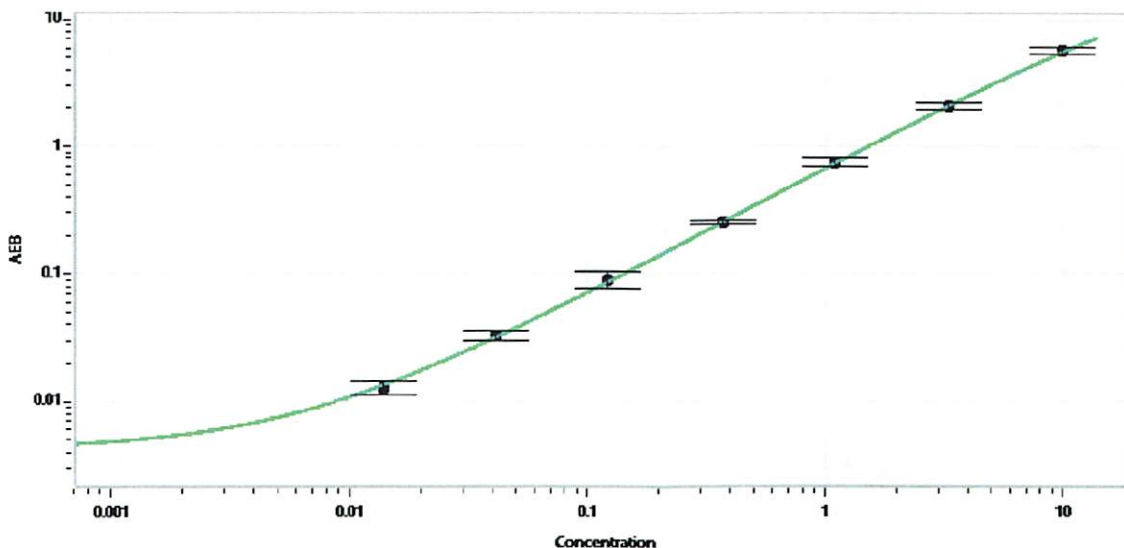
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108314	0.159	0.076-0.256
Control 2	108314	13.0	10.1-15.8
Panel 1	030001	0.171	0.103-0.238
Panel 3	030001	25.0	19.5-39.0
Panel 4	030001	0.141	0.109-0.216

\* 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 Name, QC Manager Title, Shivani Goel 11 Mar 2022 Signature/Date