

Product Number:	101635
Lot Number:	503443
Expiration:	12-Jul-2023
Platform(s):	HD-1/ HD-X



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

Component	Part Number	Lot Number
Bead Reagent	101637	219301
Detector Reagent	101636	219606
SBG Reagent	101638	219302
Sample Diluent	101639	219303
Calibrator Diluent	101640	219307
Calibrator Concentrate	100661	219404
RGP Reagent	103159	N/A <sup>1</sup>

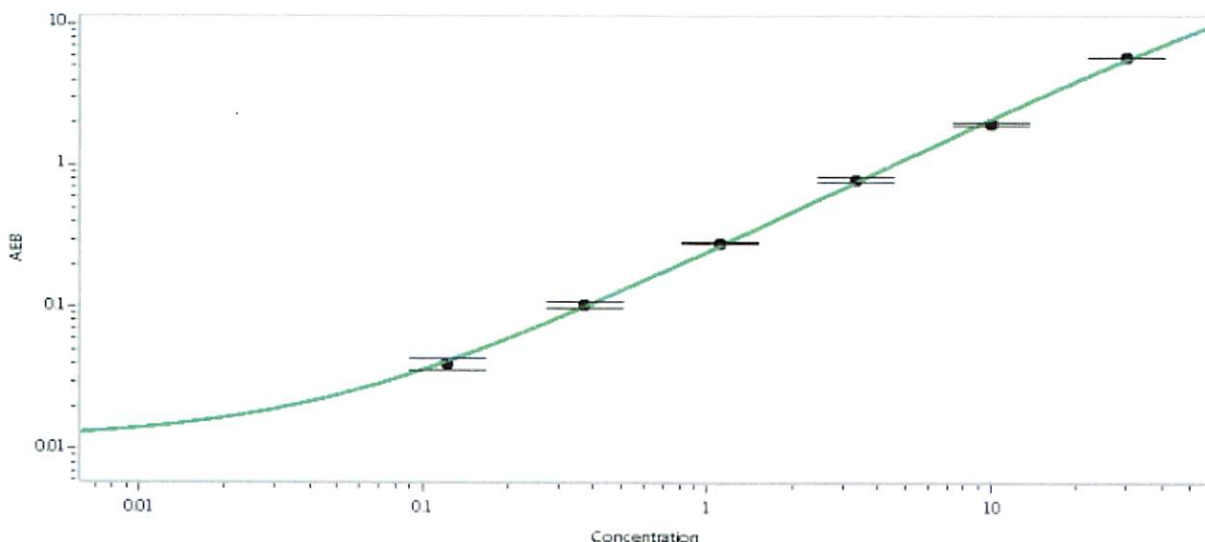
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	129010	0.560	0.389 - 0.605
Control 2	129010	54.2	38.9 - 59.7
Panel 1	035301	0.309	0.206 - 0.492
Panel 2	035301	3.72	2.65 - 4.99
Panel 3	035301	48.1	28.3 - 55.8

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

Calibrator Concentrate Value Assignment (pg/mL)	2055
---	------

Calibration Curve



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

Review/Approval: Sonia Khillan Name: Sonia Khillan Title: QC Signature/Date: Sonia Rolk 04 APR 2023