

Product Number:	101605
Lot Number:	503310
Expiration:	19-Dec-2022
Platform(s):	HD-1 / HD-X



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

Component	Part Number	Lot Number
Bead Reagent	101607	202307
Detector Reagent	101606	203711
SBG Reagent	101608	202506
Sample Diluent	103200	202507
Calibrator Diluent	101610	135311
Calibrator Concentrate	100895	202206
RGP Reagent	103159	N/A <sup>1</sup>

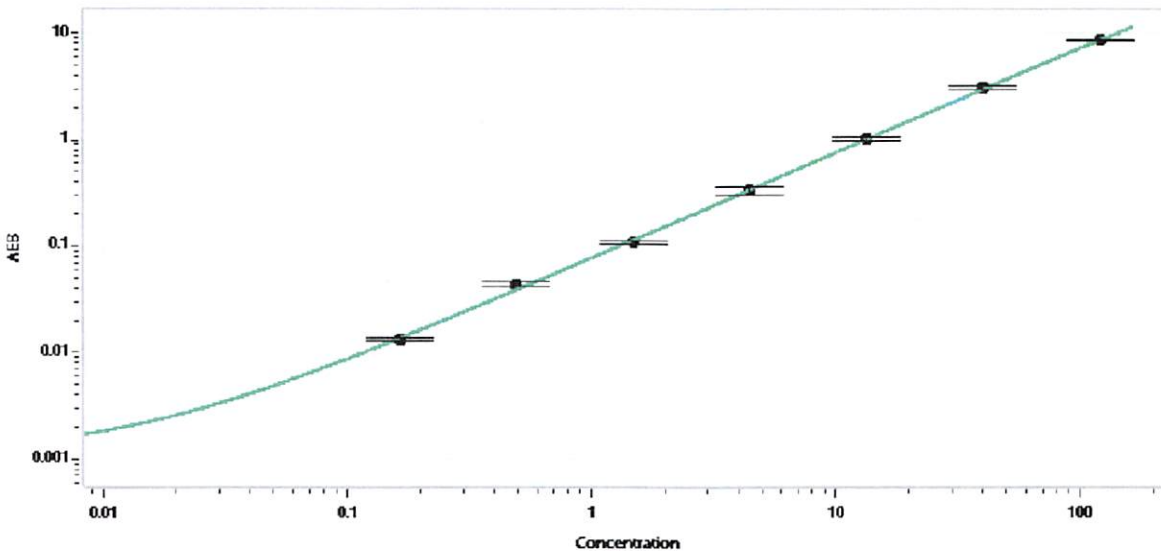
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	202606	1.47	0.907-1.60
Control 2	202606	84.7	70.2-124
Panel 2	124703	1.38	0.985-1.95
Panel 3	124703	8.82	6.49-13.3
Panel 4	124703	46.8	38.5-78.1

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

Calibrator Concentrate Value Assignment (pg/mL)	3130
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.165
C	0.494
D	1.48
E	4.44
F	13.3
G	40.0
H	120

Review/Approval: Sonia Khillan Name, QC Manager Title, Sonia Khillan / 31 AUG 2022 Signature/Date