

Product Number:	101672
Lot Number:	503538
Expiration:	5-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	101674	219406
Detector Reagent	101673	219507
SBG Reagent	101675	219508
Sample Diluent	101676	219410
Calibrator Diluent	101677	218615
Calibrator Concentrate	101734	220910
RGP Reagent	103159	N/A ¹

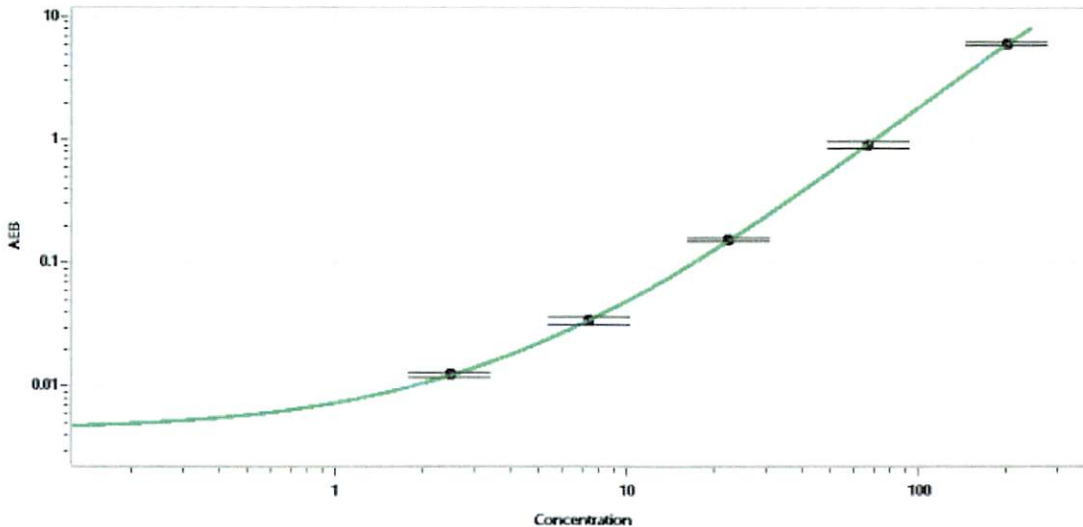
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	220910	19.5	15.0 - 25.0
Control 2	220910	388	300 - 500
Panel 1	024003	125	93.6 - 149
Panel 2	024003	275	198 - 326

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

Calibrator Concentrate Value Assignment (pg/mL)	3870
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.47
C	7.41
D	22.2
E	66.7
F	200

Review/Approval

Sonia Khilkin QC Manager Sonia Khilkin 13 MAR 2023
Name Title Signature/Date