

Product Number:	101622
Lot Number:	503656
Expiration:	28-Apr-2023
Platform(s):	HD-1 / HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101624	211809
Detector Reagent	101623	220703
SBG Reagent	101625	220901
Sample Diluent	103732	220903
Calibrator Diluent	101627	218618
Calibrator Concentrate	100652	126802
RGP Reagent	103159	N/A <sup>1</sup>

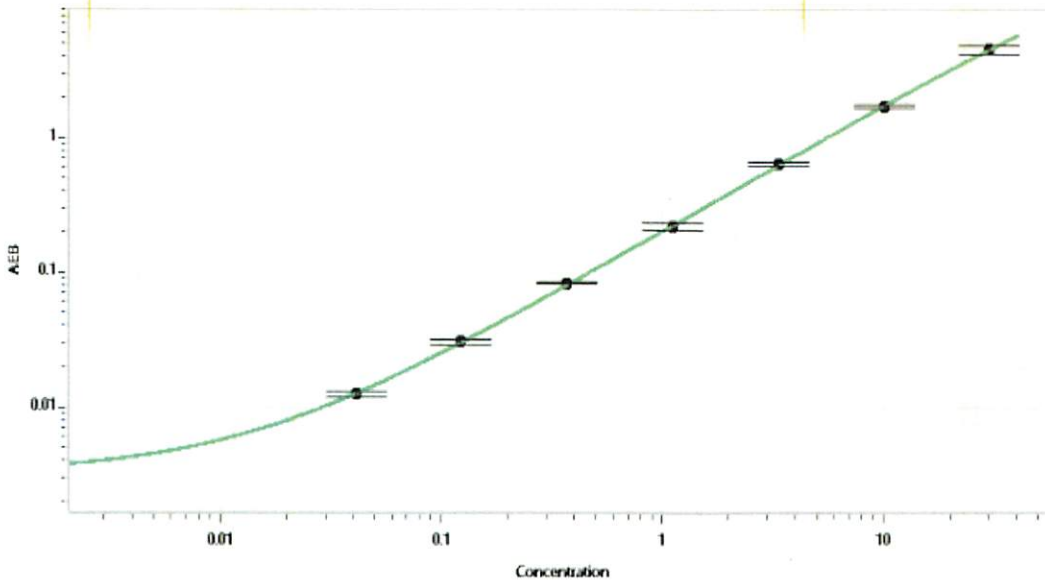
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	126802	2.27	1.50 - 2.50
Control 2	126802	20.0	15.0 - 25.0
Panel 1	029301	45.3	39.8 - 59.7
Panel 2	029401	5.14	4.35 - 6.53
Panel 3	029301	1.39	1.22 - 1.82

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

Calibrator Concentrate Value Assignment (pg/mL)	541
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval: Sonia Khillan BC Manager Sonia Khillan / 15SEP2022  
 Name Title Signature/Date