

Product Number:	104073
Lot Number:	503989
Expiration:	2-Nov-2024
Platform(s):	HD-X



Component	Part Number	Lot Number
Bead Reagent	104067	330619
Detector Reagent	104068	331011
SBG Reagent	104069	331003
Sample Diluent	104071	331010
Calibrators	102255	331001
Control 1	102256	331002
Control 2	102257	331002
RGP Reagent	103159	N/A <sup>1</sup>

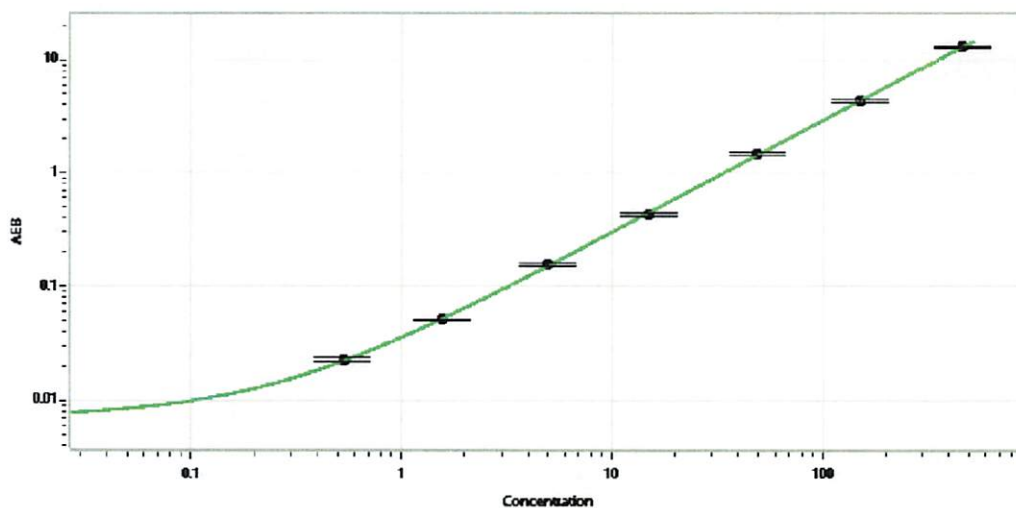
Data below represents results generated on the Simoa HD-X® Analyzer.

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

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	331002	14.7	11.5 - 17.3
Control 2	331002	612	472 - 708
Panel 1	224402	10.4	7.86 - 11.8
Panel 2	224402	7.44	5.60 - 8.59
Panel 3	224402	72.4	56.1 - 84.1
Panel 4	301701	560	462 - 719

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.531
C	1.57
D	4.94
E	14.9
F	49.2
G	150
H	463

Review/Approval

*Sonia Khillan*  
Name

QC  
Title

*Sonia Khillan 04/DEC/2023*  
Signature/Date