

Product Number:	102860
Lot Number:	503714
Expiration:	2-May-2023
Platform(s):	HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	102852	230605
Detector Reagent	102854	231208
SBG Reagent	102855	231209
Sample Diluent	102856	231211
Calibrator Diluent	102859	231212
Calibrator Concentrate	102857	308007
RGP Reagent	103159	N/A <sup>1</sup>

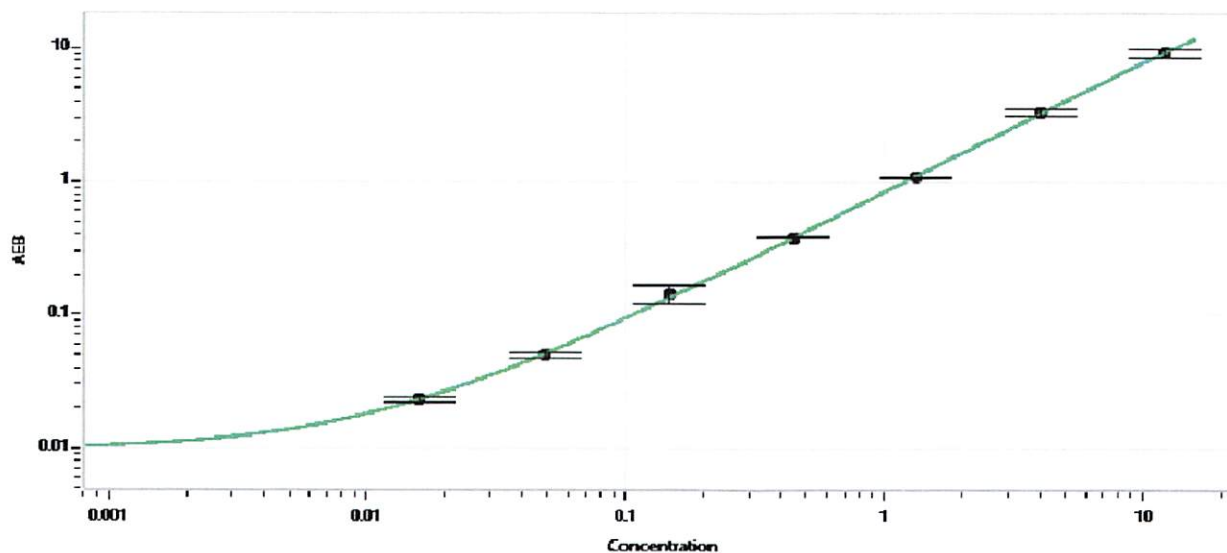
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	121506	0.233	0.229 - 0.344
Control 2	121506	4.30	3.93 - 5.90
Panel 1	231306	0.283	0.227 - 0.340
Panel 2	234108	2.48	2.47 - 3.71
Panel 3	234108	9.06	8.61 - 13.2

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

Calibrator Concentrate Value Assignment (pg/mL)	2329
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.016
C	0.049
D	0.148
E	0.444
F	1.33
G	4.00
H	12.0

Review/Approval

*Matt Koulouris*  
Name

QC Supervisor  
Title

*[Signature]* 05-Apr-2023  
Signature/Date