

Product Number:	102329
Lot Number:	503384
Expiration:	5-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	102322	206302
Detector Reagent	102324	206402
SBG Reagent	102325	206401
Sample Diluent	102326	133938
Calibrator Diluent	102327	204613
Calibrator Concentrate	100849	126310
RGP Reagent	103159	N/A ¹

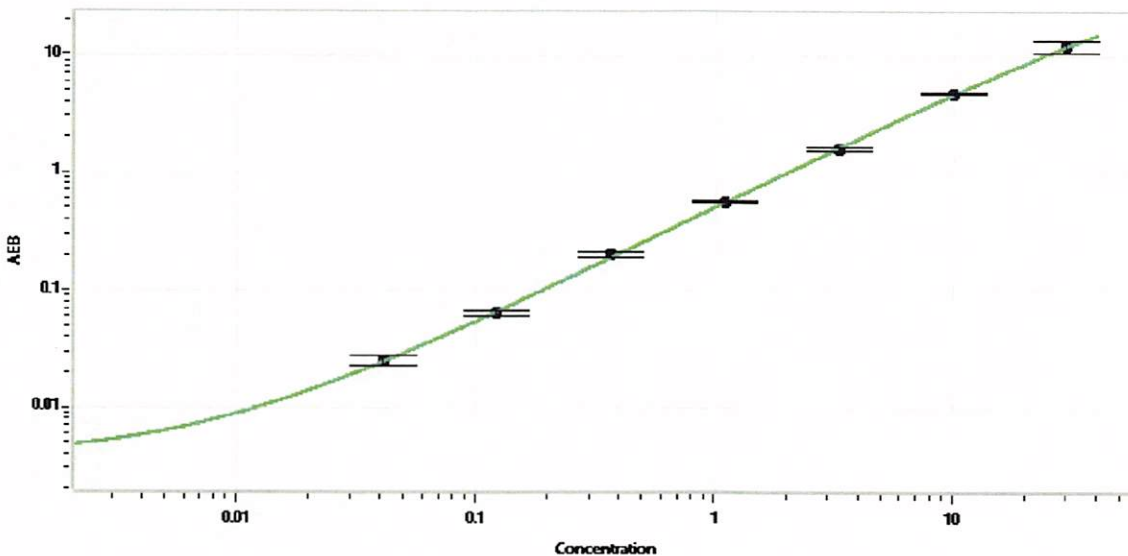
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	126310	0.203	0.150-0.250
Control 2	126310	21.5	18.8-31.3
Panel 1	027607	0.142	0.050-0.176
Panel 2	027607	0.153	0.071-0.213
Panel 3	027607	11.4	7.45-18.2

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

Calibrator Concentrate Value Assignment (pg/mL)	519
---	-----

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

Shivani Goel

QC Manager

Shivani Goel 06Apr2022

Name

Title

Signature/Date