

Product Number:	102732
Lot Number:	503411
Expiration:	14-Jul-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	102722	226502
Detector Reagent	102724	227301
SBG Reagent	102726	227302
Sample Diluent	102728	219512
Calibrator Diluent	102731	227313
Calibrator Concentrate	100658	204302
RGP Reagent	103159	N/A ¹

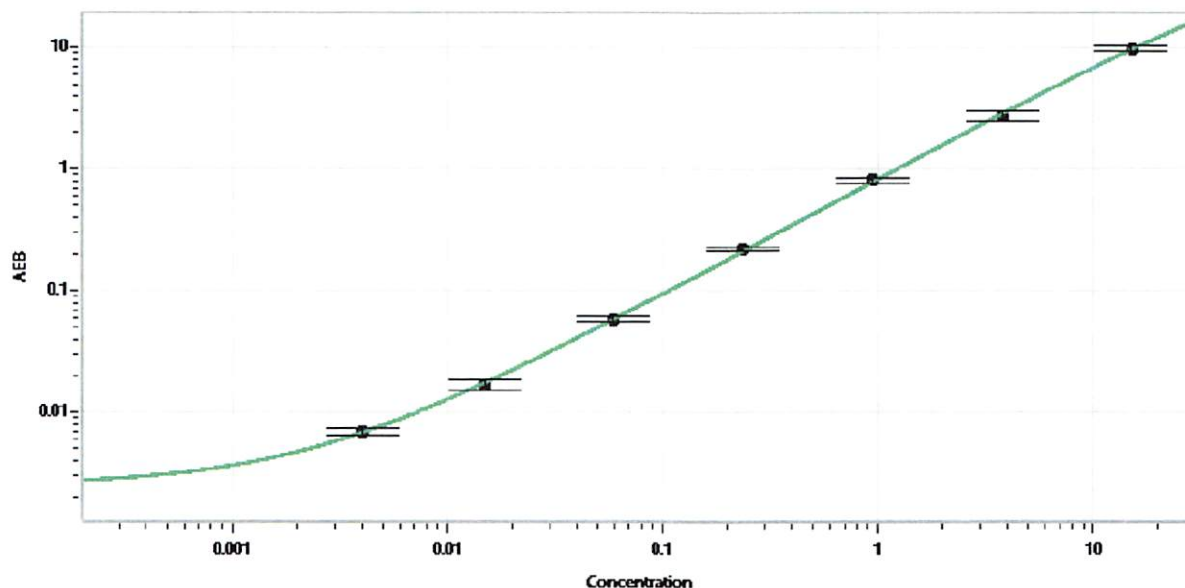
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	114108	0.083	0.068-0.168
Control 2	114108	2.78	1.92-4.04
Panel 2	112714	0.031	0.009-0.056
Panel 3	112714	1.67	0.767-2.36
Panel 4	112714	6.51	2.38-8.94

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

Calibrator Concentrate Value Assignment (pg/mL)	375
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.004
C	0.015
D	0.059
E	0.234
F	0.938
G	3.75
H	15.0

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia 12/16/2023
Signature/Date