

Product Number:	102732
Lot Number:	503344
Expiration:	14-Feb-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	219501
Detector Reagent	102724	204801
SBG Reagent	102726	204802
Sample Diluent	102728	219511
Calibrator Diluent	102731	204514
Calibrator Concentrate	100658	106608
RGP Reagent	103159	N/A ¹

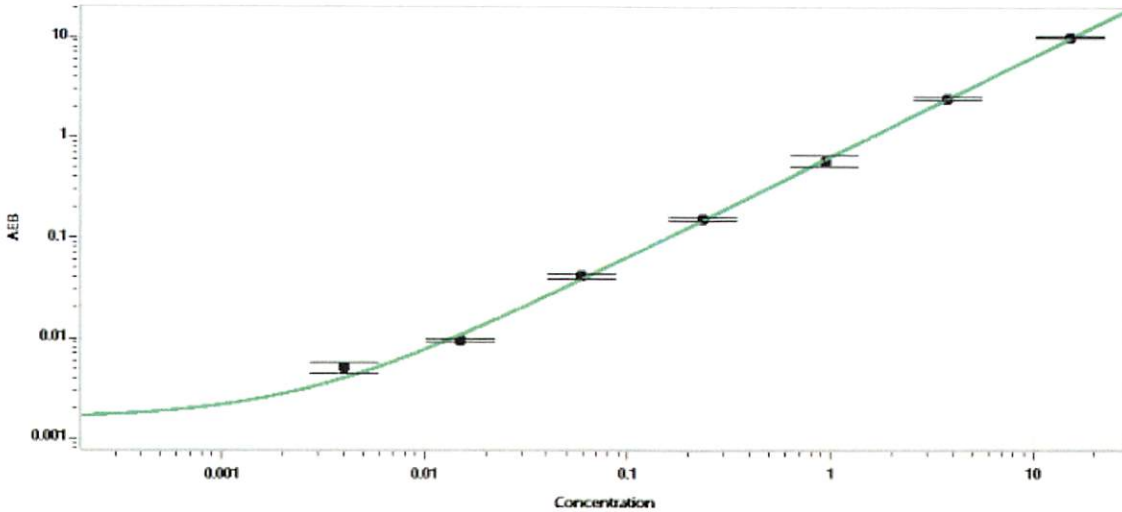
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	114108	0.139	0.068 - 0.168
Control 2	114108	3.45	1.92 - 4.04
Panel 2	112714	0.043	0.009 - 0.056
Panel 3	112714	2.15	0.767 - 2.36
Panel 4	112714	7.63	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)	444
---	-----

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

Title

Sonia Khillan / 14 OCT 2022

Signature/Date