

Product Number:	101132
Lot Number:	503719
Expiration:	29-Aug-2023
Platform(s):	HD-1/HD-X



Data below represents results generated on the Simoa® HD-1 / HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101119	227903
Detector Reagent	101121	228001
SBG Reagent	101131	228002
Sample Diluent	103319	227230
Calibrator Diluent	101125	227712
Calibrator Concentrate	101123	203904
RGP Reagent	103159	N/A ¹

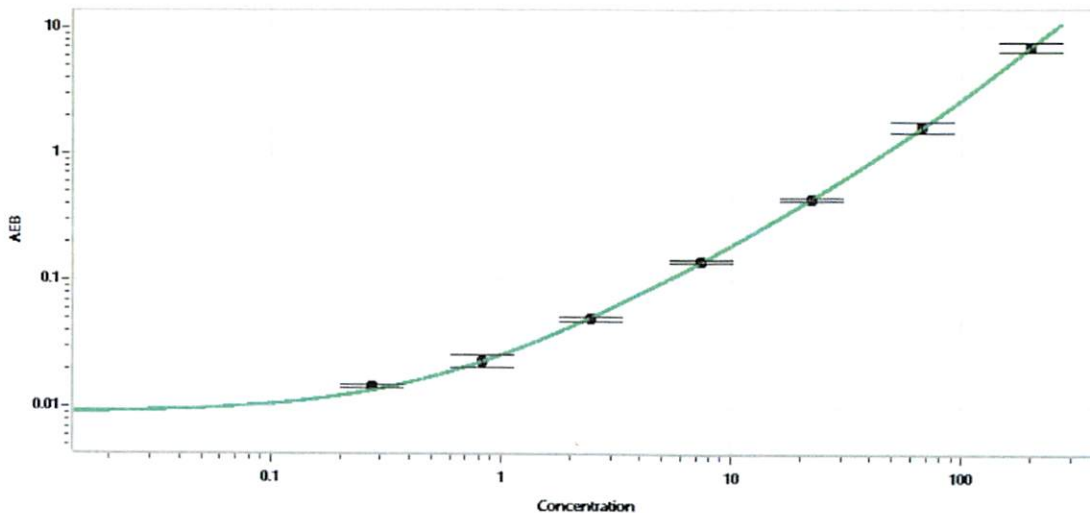
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	130713	8.99	6.43 - 10.1
Control 2	130713	298	151 - 332
Panel 1	106201	161	94.7 - 210
Panel 2	106201	99.2	59 - 123
Panel 3	106201	298	183 - 365
Panel 4	106201	295	171 - 332

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

Calibrator Concentrate Value Assignment (pg/mL)	5081
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Sonia Khillar
Name

QC
Title

Sonia Khillar 08MAR2023
Signature/Date