

Product Number:	104542
Lot Number:	504054
Expiration:	18-Oct-2024
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	104515	331715
Detector Reagent	104517	331907
SBG Reagent	104519	331905
Sample Diluent	104521	332109
Calibrator Diluent	104524	329115
Control Diluent	104525	329116
Calibrator Concentrate	104538	333907
Control 1 Concentrate	104540	333909
Control 2 Concentrate	104541	333910
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific

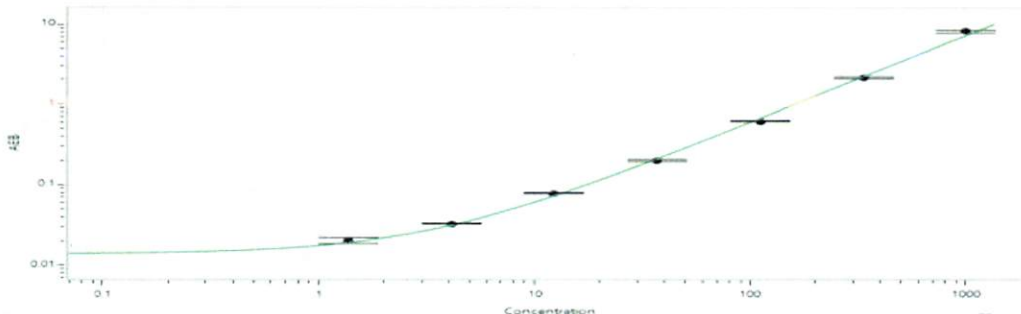
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	333909	309	184 - 486
Control 2 Concentrate	333910	4068	2156 - 5857
Panel 1	329315	382	242 - 450
Panel 2	329315	4565	3166 - 7069
Panel 3	N/A	N/A	N/A
Panel 4	N/A	N/A	N/A

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

Calibrator Concentrate Reconstitution Volume (uL)	2000
Control 1 Concentrate Reconstitution Volume (uL)	2000
Control 2 Concentrate Reconstitution Volume (uL)	2000

Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.12
D	12.3
E	37.0
F	111
G	333
H	1000

Calibration Curve



Review/Approval

ROBERT SOMOGYI
Name

QC MANAGER
Title

[Signature] 12 JAN 2024
Signature/Date