

Product Number:	104428
Lot Number:	504024
Expiration:	13-Oct-2024
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	104409	330310
Detector Reagent	104411	329602
SBG Reagent	104413	329601
Sample Diluent	104415	329113
Calibrator Diluent	104418	328617
Control Diluent	104476	328619
Calibrator Concentrate	104483	330609
Control 1 Concentrate	104426	330611
Control 2 Concentrate	104427	330612
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific

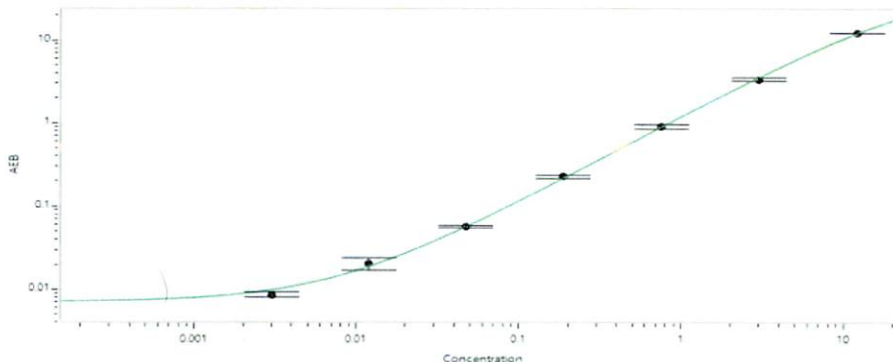
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	330611	0.208	0.103 - 0.227
Control 2 Concentrate	330612	5.14	3.08 - 6.89
Panel 1	328501	0.030	0.013 - 0.032
Panel 2	328501	0.677	0.379 - 0.704
Panel 3	328501	3.04	1.82 - 3.37
Panel 4	328501	6.07	3.81 - 7.08

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

Calibrator Concentrate Reconstitution Volume (uL)	858
Control 1 Concentrate Reconstitution Volume (uL)	858
Control 2 Concentrate Reconstitution Volume (uL)	858

Calibrator Levels (pg/mL)	
A	0
B	0.003
C	0.012
D	0.047
E	0.188
F	0.750
G	3.00
H	12.0

Calibration Curve



Review/Approval

ROBERT SOMOGYI

QC MANAGER

[Signature] 12/11/2024

Name

Title

Signature/Date