

Product Number:	104364
Lot Number:	503994
Expiration:	22-Oct-2025
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	104352	323506
Detector Reagent	104353	323406
SBG Reagent	104354	323704
Sample Diluent	104355	326511
Calibrator Diluent	104358	324103
Control Diluent	104359	324104
Calibrator Concentrate	104361	317309
Control 1 Concentrate	104362	317310
Control 2 Concentrate	104363	317311
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

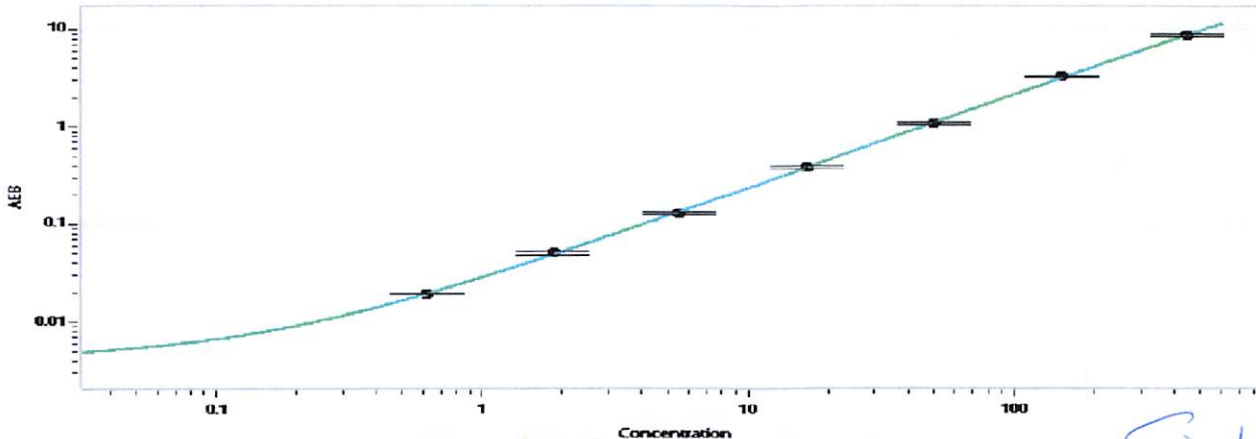
¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	317310	4.30	2.97 - 6.15
Control 2	317311	380	264 - 479
Panel 1	224402	9.46	7.86 - 11.8
Panel 2	224402	6.70	5.60 - 8.59
Panel 3	224402	74.9	56.1 - 84.1
Panel 4	301701	598	462 - 719

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

Calibrator and Controls Concentrate Reconstitution Volume (uL)	1236
--	------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.617
C	1.85
D	5.56
E	16.7
F	50.0
G	150
H	450

Review/Approval

Sonia Khilka
Name

QC Manager
Title

Sonia Khilka 09 OCT 2023
Signature/Date