

Product Number:	104570
Lot Number:	504348
Expiration:	12-Feb-2026
Platform(s):	HD-X



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

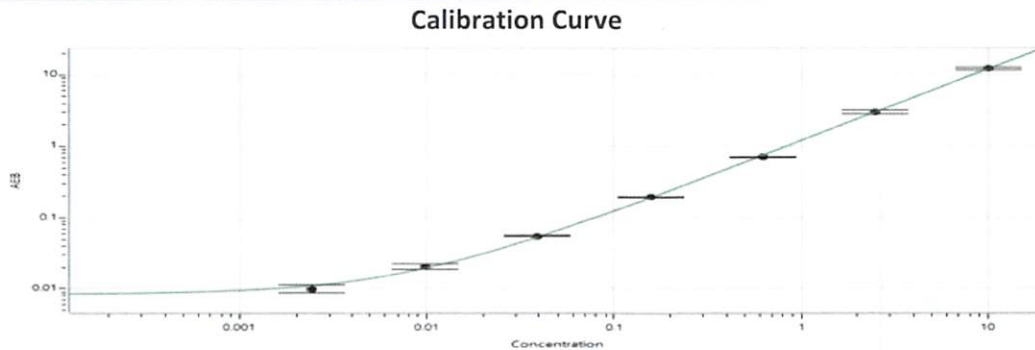
Component	Part Number	Lot Number
Bead Reagent	104546	422503
Detector Reagent	104548	422712
SBG Reagent	104550	422713
Sample Diluent	104552	422619
Calibrator Diluent Reagent	104745	422621
Control Diluent Reagent	104554	422620
Calibrator Concentrate	104566	406408
Control 1 Concentrate	104568	406409
Control 2 Concentrate	104569	406410
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific

Release Materials*			* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	
Control 1 Concentrate	406409	0.692	0.494 - 0.917
Control 2 Concentrate	406410	7.08	5.03 - 9.33

Calibrator Concentrate Reconstitution Volume (uL)	842
Control 1 Concentrate Reconstitution Volume (uL)	842
Control 2 Concentrate Reconstitution Volume (uL)	842

Calibrator Levels (pg/mL)	
A	0.000
B	0.00244
C	0.00977
D	0.0391
E	0.156
F	0.625
G	2.50
H	10.0



Review/Approval

TAM LEE
 DOC CONTROL II, QA
 Tam Lee
 20 SEP 2024

Name	Title	Signature/Date
------	-------	----------------