

Product Number:	104570
Lot Number:	504500
Expiration:	14-Aug-2026
Platform(s):	HD-X and SR-X



Data below represents results generated on the

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	104546	504502	504502A
Detector Reagent	104548	505001	N/A
SBG Reagent	104550	505002	N/A
Sample Diluent	104552	504510	504510A
Calibrator Diluent Reagent	104745	504512	504512A
Control Diluent Reagent	104554	504511	504511A
Calibrator Concentrate	104566	405210	N/A
Control 1 Concentrate	104568	405211	N/A
Control 2 Concentrate	104569	405212	N/A
RGP 2.0 Reagent	104321	N/A ¹	N/A ¹
Activation Buffer Reagent	104324	N/A ¹	N/A ¹

¹ Reagents are not Kit Lot Specific

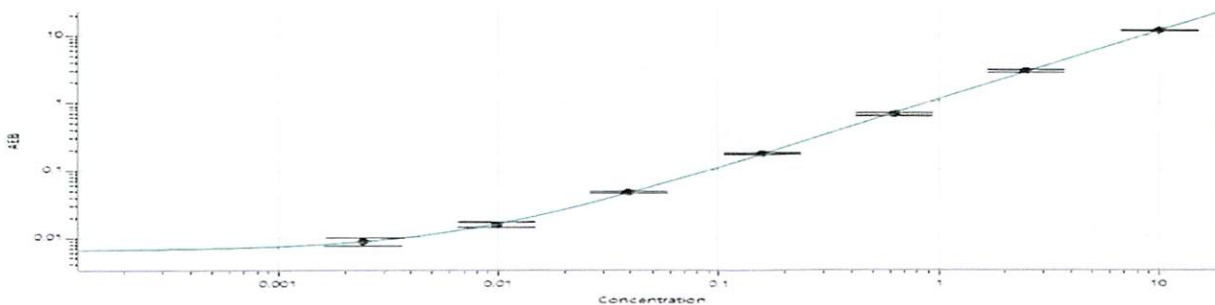
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	405211	0.811	0.517 - 0.959
Control 2 Concentrate	405212	7.92	5.22 - 9.70

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

Calibrator Concentrate Reconstitution Volume (uL)	970
Control 1 Concentrate Reconstitution Volume (uL)	970
Control 2 Concentrate Reconstitution Volume (uL)	970

Calibrator Levels (pg/mL)	
A	0
B	0.00244
C	0.00977
D	0.0391
E	0.156
F	0.625
G	2.50
H	10.0

Calibration Curve



Review/Approval

Dilish Patel
Name

Sr. Document QA
Title

ET 27 MAR 2025
Signature/Date