

<b>Product Number:</b>	101580
<b>Lot Number:</b>	504515
<b>Expiration:</b>	5-Aug-2026
<b>Platform(s):</b>	SR-X and HD-1/HD-X



*Data below represents results generated on the Simoa<sup>®</sup> HD-X Analyzer.*

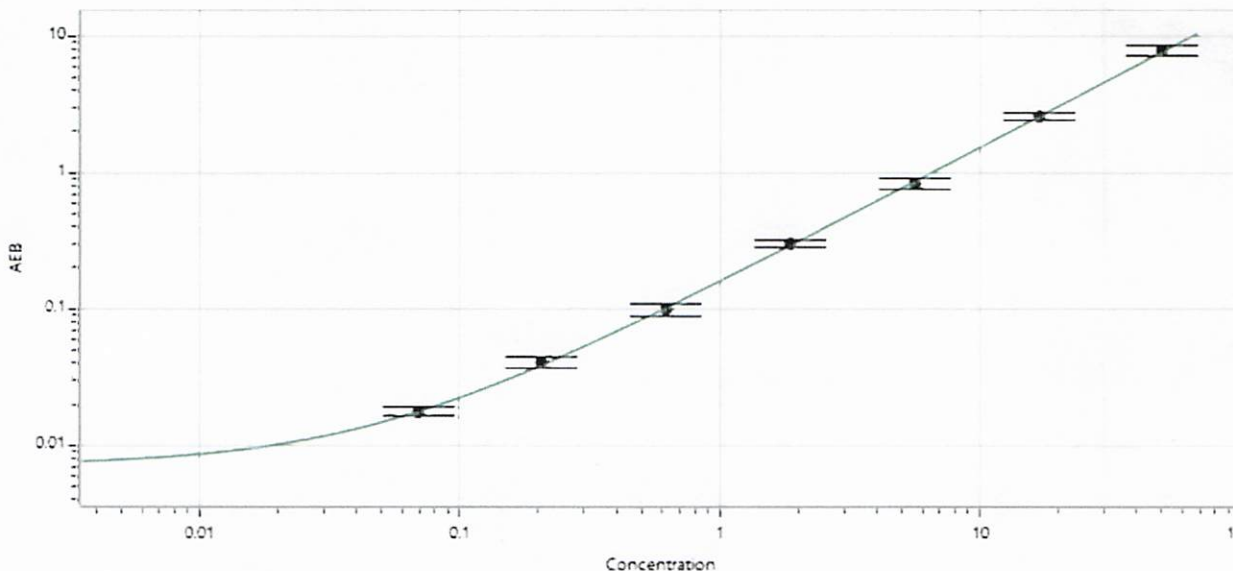
Component	Part Number	Lot Number
Bead Reagent	104236	507207
Detector Reagent	104327	507209
SBG Reagent	104329	507302
Sample Diluent	104331	503611
Calibrator Diluent	101585	503602
Calibrator Concentrate	104332	507708
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
<b>Control 1</b>	206502	2.58	2.00 - 3.72
<b>Control 2</b>	206502	43.3	33.0 - 61.4

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

**Calibrator Concentrate Value Assignment (pg/mL)** 1558



**Calibrator Levels (pg/mL)**

A	0.000
B	0.069
C	0.206
D	0.617
E	1.85
F	5.56
G	16.7
H	50.0

Review/Approval

Dipesh Patel Sr. Doc. Control QA [Signature] 26MAR2025 [Date]  
Name Title Signature/Date