

<b>Product Number:</b>	105236
<b>Lot Number:</b>	504426
<b>Expiration:</b>	21-Apr-2026
<b>Platform(s):</b>	HD-X



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

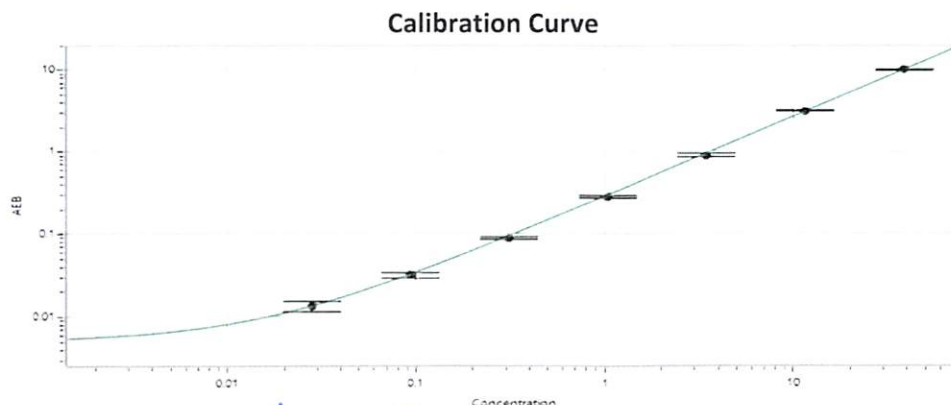
Component	Part Number	Lot Number
Bead Reagent	105215	429801
Detector Reagent	105217	429703
SBG Reagent	105219	429704
Sample Diluent	105220	429516
Calibrator Diluent	105222	429520
Control Diluent	105223	429519
Calibrator Concentrate	105232	411422
Control 1 Concentrate	105234	411424
Control 2 Concentrate	105235	411425
RGP 2.0 Reagent	104321	N/A <sup>1</sup>
Activation Buffer Reagent	104324	N/A <sup>1</sup>

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

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	411424	1.94	1.31 - 2.44
Control 2 Concentrate	411425	86.5	64.9 - 121

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

Calibrator Concentrate Reconstitution Volume (uL)	960
Control 1 Concentrate Reconstitution Volume (uL)	960
Control 2 Concentrate Reconstitution Volume (uL)	960



Calibrator Levels (pg/mL)	
A	0.000
B	0.028
C	0.094
D	0.312
E	1.04
F	3.46
G	11.5
H	38.4

Review/Approval Dhruvi Patel DOC Control I, QA SDP 05 Dec 2024  
 Name Title Signature/Date