

<b>Product Number:</b>	105187
<b>Lot Number:</b>	504422
<b>Expiration:</b>	21-Apr-2026
<b>Platform(s):</b>	HD-X



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

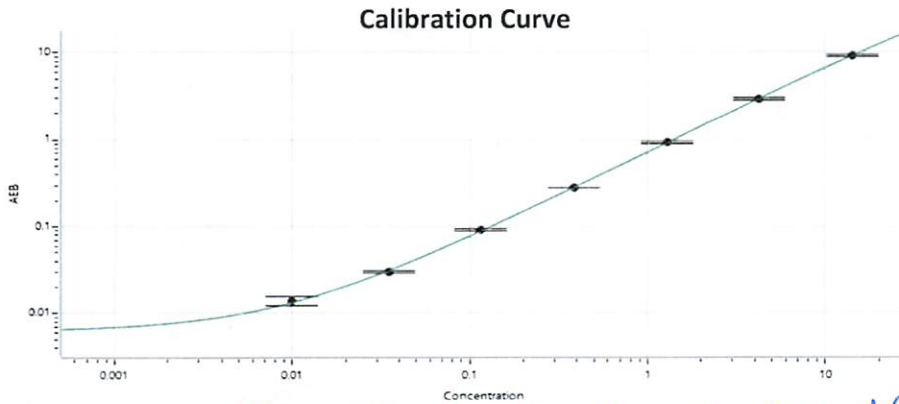
Component	Part Number	Lot Number
Bead Reagent	105139	429601
Detector Reagent	105141	429802
SBG Reagent	105143	429803
Sample Diluent	105144	429513
Calibrator Diluent	105172	429514
Control Diluent	105173	429515
Calibrator Concentrate	105183	412323
Control 1 Concentrate	105185	412324
Control 2 Concentrate	105186	412325
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	412324	0.700	0.432 - 0.893
Control 2 Concentrate	412325	16.7	12.3 - 22.9

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

<b>Calibrator Concentrate Reconstitution Volume (uL)</b>	875
<b>Control 1 Concentrate Reconstitution Volume (uL)</b>	875
<b>Control 2 Concentrate Reconstitution Volume (uL)</b>	875



Calibrator Levels (pg/mL)	
A	0.000
B	0.010
C	0.035
D	0.115
E	0.385
F	1.28
G	4.26
H	14.2

Review/Approval: TAM LEE DOC CONTROL II, QA Man Lee 22 NOV 2024  
Name Title Signature/Date