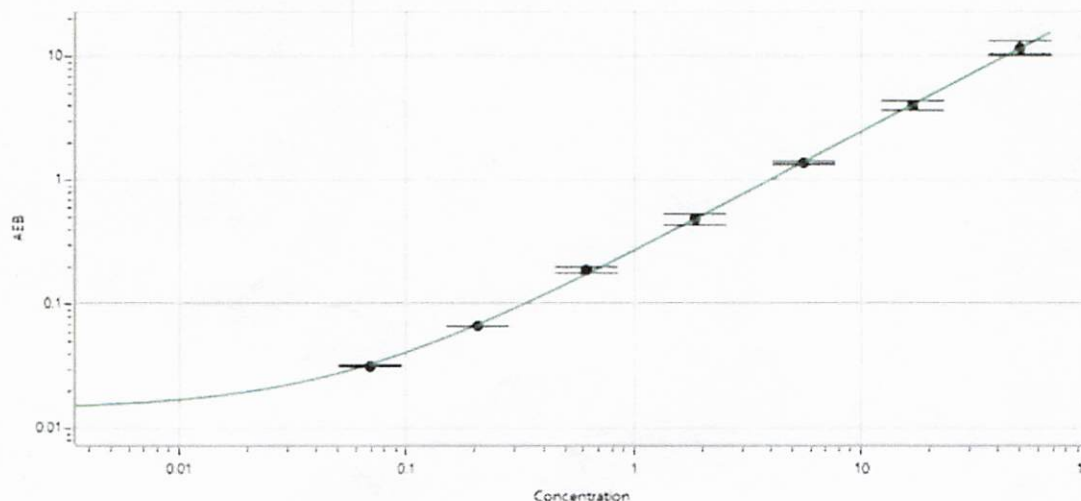


Component	Part Number	Lot Number
Bead Reagent	104236	514003
Detector Reagent	104327	514705
SBG Reagent	104329	516001
Sample Diluent	104331	503613
Calibrator Diluent	101585	510114
Calibrator Concentrate	104332	507708
RGP Reagent	103159	N/A ¹

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Spec. Range (pg/mL)	
Control 1	206502	3.09	2.00 - 3.72	
Control 2	206502	50.7	33.0 - 61.4	

1558



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

Dillesh P. H. Sr. Doc. Control - QA at / 13 Jun 2025
Name Title Signature/Date