

Product Number:	102215
Lot Number:	504457
Expiration:	24-Mar-2026
Platform(s):	SR-X and HD-1/HD-X



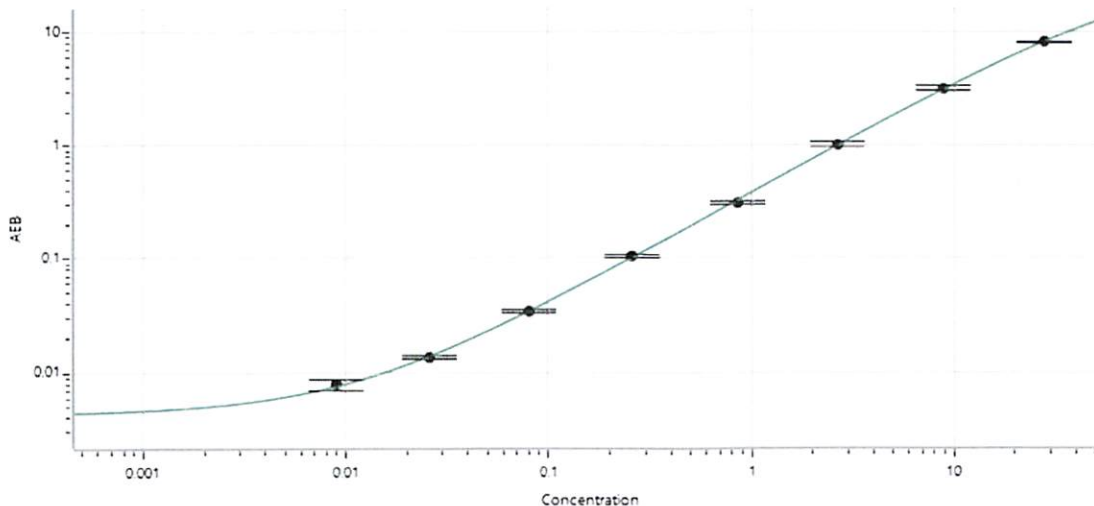
Data below represents results generated on the Simoa® HD-X.

Component	Part Number	Lot Number
Bead Reagent	104297	433012
Detector Reagent	104338	433006
SBG Reagent	104340	433008
Sample Diluent	104342	433015
Calibrators	104343	433004
Control 1	104344	433005
Control 2	104345	433005
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

Release Materials*			* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	
Control 1	433005	0.083	0.063 - 0.098
Control 2	433005	9.09	7.55 - 11.3

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.009
C	0.026
D	0.080
E	0.259
F	0.852
G	2.64
H	8.89
I	27.9

Review/Approval: Dipesh Patel Name, Doc Control II QA Title, [Signature] Signature/Date, 11 DEC 2024 Signature/Date