

Product Number:	104292
Lot Number:	503947
Expiration:	21-Jun-2024
Platform(s):	HD-1 / HD-X



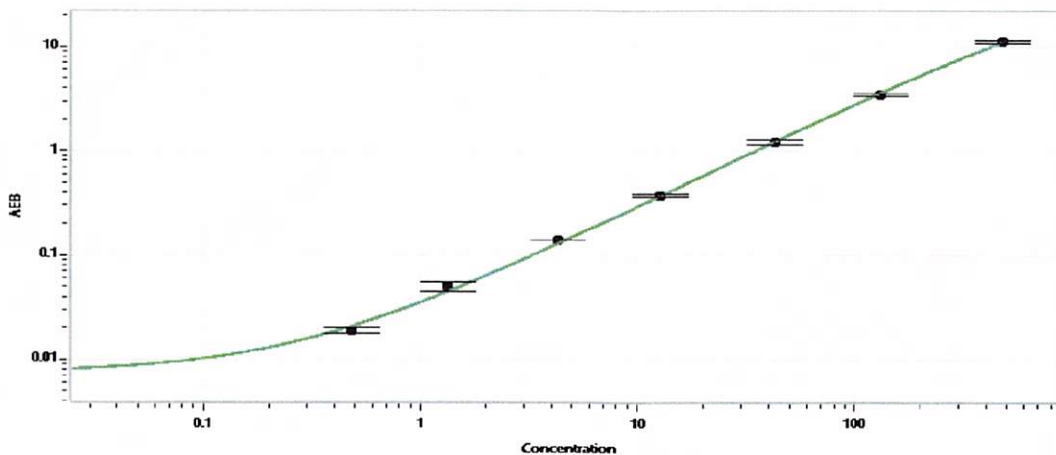
Component	Part Number	Lot Number	Control Number
Bead Reagent	104067	319905	319905A
Detector Reagent	104068	320706	N/A
SBG Reagent	104069	317209	317209A
Sample Diluent	104071	317210	317210A
Calibrators	102255	317804	317804A
Control 1	102256	317805	317805A
Control 2	102257	317805	317805A
RGP Reagent	103159	N/A ¹	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	317805A	11.5	9.88 - 14.8
Control 2	317805A	510	417 - 625
Panel 1	224402	9.88	7.86- 11.8
Panel 2	224402	6.94	5.60-8.59
Panel 3	224402	70.6	56.1-84.1
Panel 4	301701	549	462 - 719

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

Calibration Curve



Review/Approval Sonia Khillan QC Manager Sonia Khillan - 16 AUG 2023
 Name Title Signature/Date