

Product Number:	101552
Lot Number:	503320
Expiration:	14-Oct-2022
Platform(s):	HD-1/HD-X & SR-X

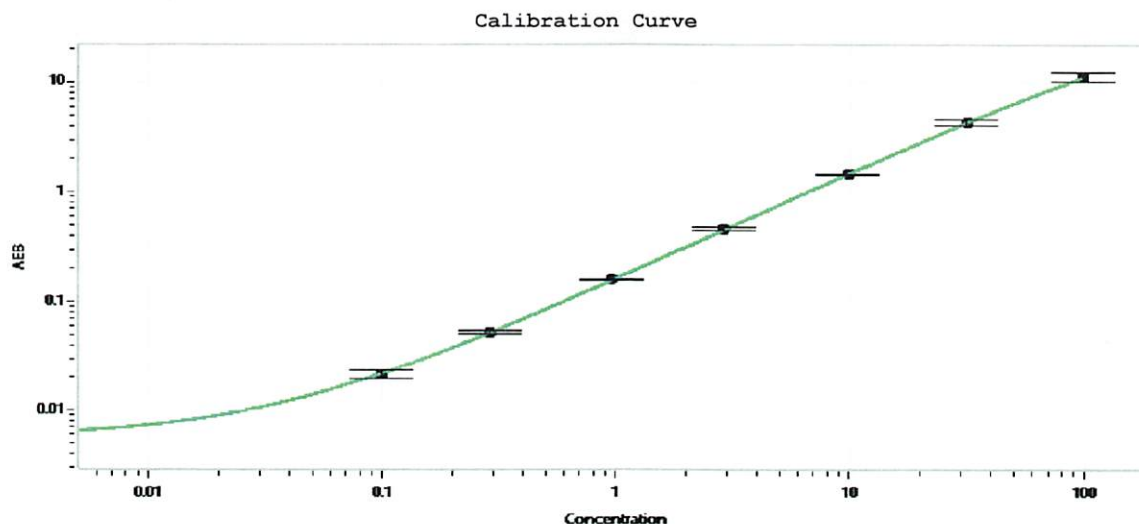


Component	Part Number	Lot Number
Bead Reagent	101554	128704
Detector Reagent	101553	135505
SBG Reagent	101555	128804
Sample Diluent	101556	128805
Calibrators	101559	136005
Control 1	101560	136006
Control 2	101561	136006
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Data below represents results generated on the Simoa® HD-X / HD-1 & SR-X Analyzer.

Release Materials*				
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	136006	0.993	0.805-1.21	* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
Control 2	136006	64.1	46.7-79.2	
Panel 1	117205	4.58	3.48 - 5.22	
Panel 2	117205	2.70	2.00 - 4.90	
Panel 3	117205	23.5	17.9 -26.8	
Panel 4	119612	38.6	22.6 - 41.6	



Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 10 Mar 2022
Signature/Date