

Product Number:	102215
Lot Number:	503449
Expiration:	23-Mar-2023
Platform(s):	HD-1/HD-x & SR-X



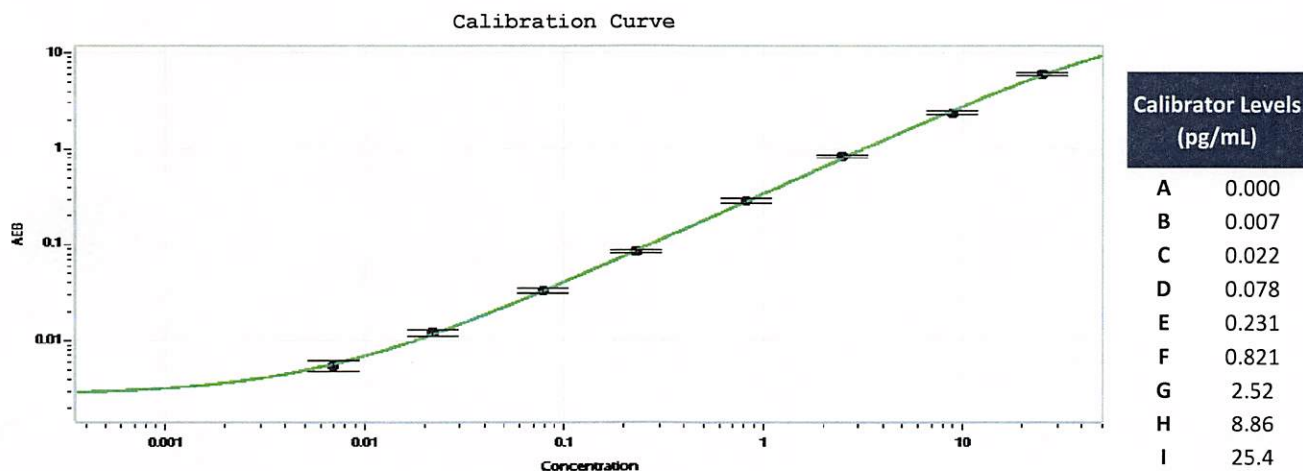
Component	Part Number	Lot Number
Bead Reagent	102211	210701
Detector Reagent	102212	210802
SBG Reagent	102213	210803
Assay Diluent	102214	210804
Calibrators	101310	210302
Control 1	100702	210303
Control 2	100703	210303
RGP Reagent	103159	N/A ¹

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

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	210303	0.079	0.062-0.092
Control 2	210303	8.81	6.69-10.6
Panel 1	019601	0.194	0.127-0.243
Panel 2	019601	2.30	1.56-2.34
Panel 3	019601	13.1	9.44-14.2

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



Review/Approval _____

Shivani Goel
Name

QC Manager
Title

Shivani Goel 09 May 2022
Signature/Date