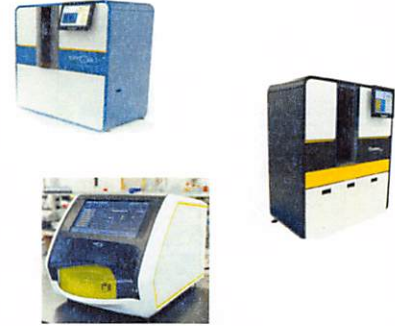


Product Number:	102215
Lot Number:	503428
Expiration:	23-Mar-2023
Platform(s):	HD-1/HD-x & SR-X



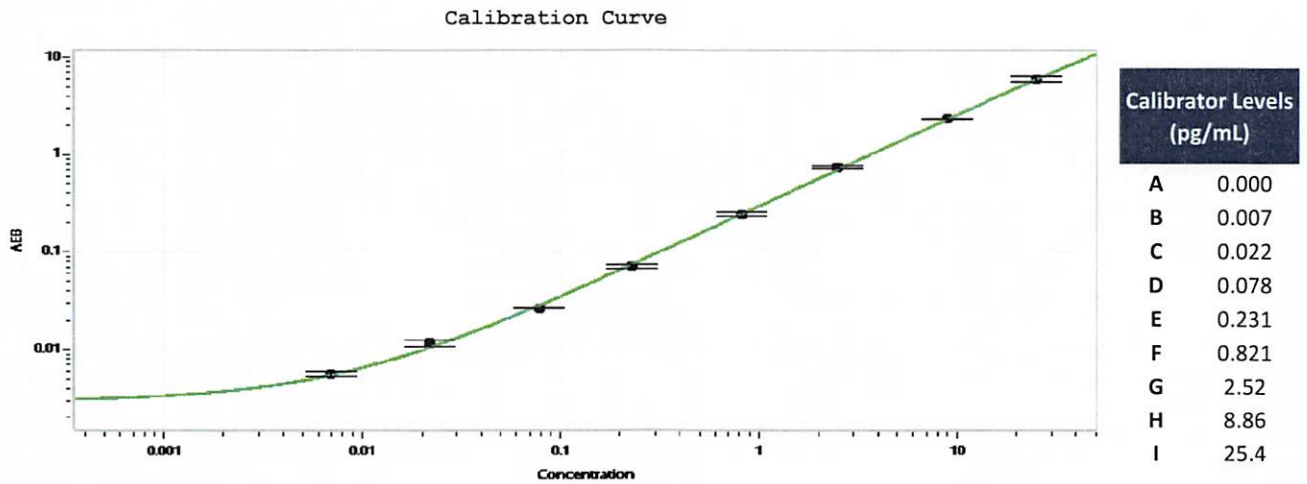
Component	Part Number	Lot Number
Bead Reagent	102211	210008
Detector Reagent	102212	210102
SBG Reagent	102213	210103
Assay Diluent	102214	210104
Calibrators	101310	210302
Control 1	100702	210303
Control 2	100703	210303
RGP Reagent	103159	N/A <sup>1</sup>

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

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	210303	0.075	0.062-0.092
Control 2	210303	9.24	6.69-10.6
Panel 1	019601	0.190	0.127-0.243
Panel 2	019601	2.23	1.56-2.34
Panel 3	019601	13.3	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 QC Manager Shivani Goel 09 May 2022

Name Title Signature/Date