

Product Number:	103714
Lot Number:	503356
Expiration:	15-Feb-2023
Platform(s):	HD-1 / HD-X & SR-X



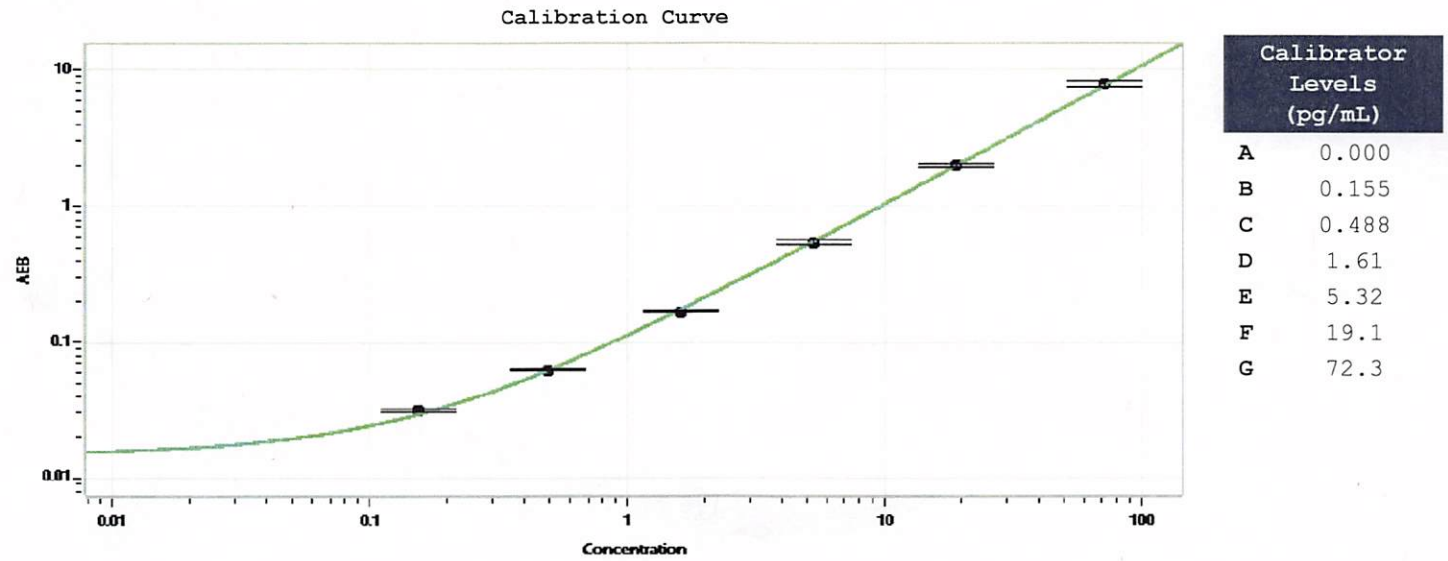
Data below represents results generated on the Simoa® HD-X

Component	Part Number	Lot Number
Bead Reagent	103698	206104
Detector Reagent	103700	206105
SBG Reagent	103702	206106
Sample Diluent	103704	206107
Reference Calibrator	103711	207204
Control 1	103712	207205
Control 2	103713	207205
RGP Reagent	103159	N/A <sup>1</sup>

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

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	207205	2.99	2.32-3.47
Control 2	207205	100	79.3-119
Panel 1	013903	0.770	0.258-1.11
Panel 2	013903	1.78	0.841-2.44
Panel 3	013903	2.54	1.79-4.31

\* 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 23 Mar 2022 Signature/Date