

Product Number:	101552
Lot Number:	502218
Expiration:	21-Nov-2020
Platform(s):	HD-1/HD-x & SR-X



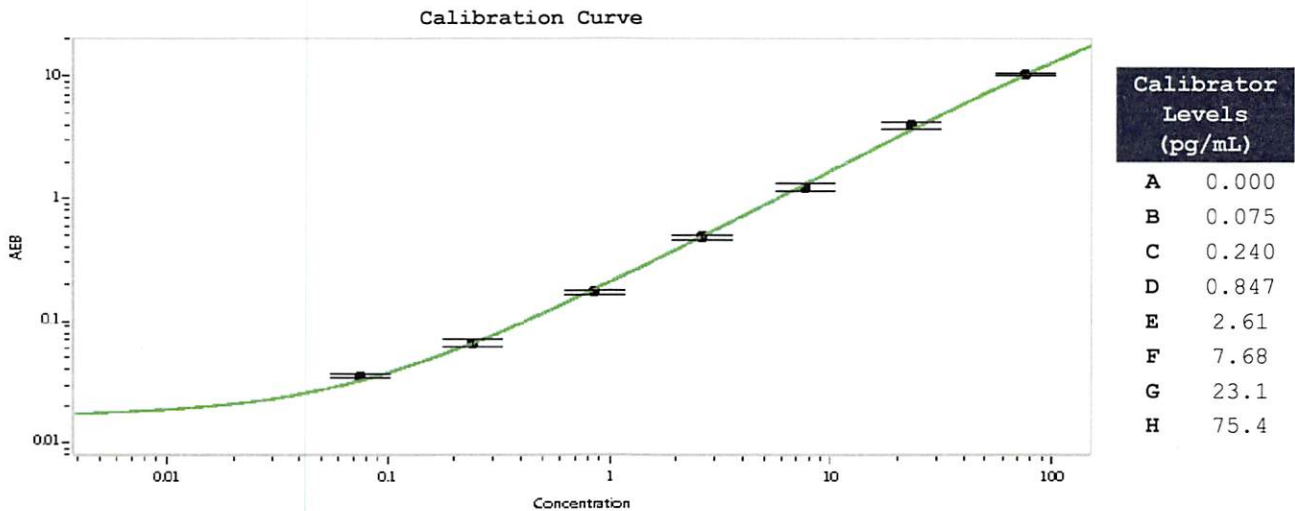
Data below represents results generated on the Simoa™ HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101554	934411
Detector Reagent	101553	933113
SBG Reagent	101555	933114
Sample Diluent	101556	933116
Calibrators	101559	003001
Controls	101560-101561	003002
SBG Diluent (SR-X) †	103158	N/A <sup>1</sup>
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	003002	0.954	0.754-1.16
Control 2	003002	43.5	33.2-49.8
Panel 1	916901	0.465	0.089-1.59
Panel 2	916902	1.03	0.538-2.13
Panel 3	916903	31.2	15.5-33.4
Panel 4	916904	96.3	63.0-102

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



Review/Approval	Patrizia Stadler Assoc. Director, QC Name Title	<i>P. Stadler</i> Signature/Date	10 Feb 2020 Date
-----------------	--	-------------------------------------	---------------------