

Product Number:	103562
Lot Number:	502879
Expiration:	21-Mar-2022
Platform(s):	HD-1/HD-x & SR-X



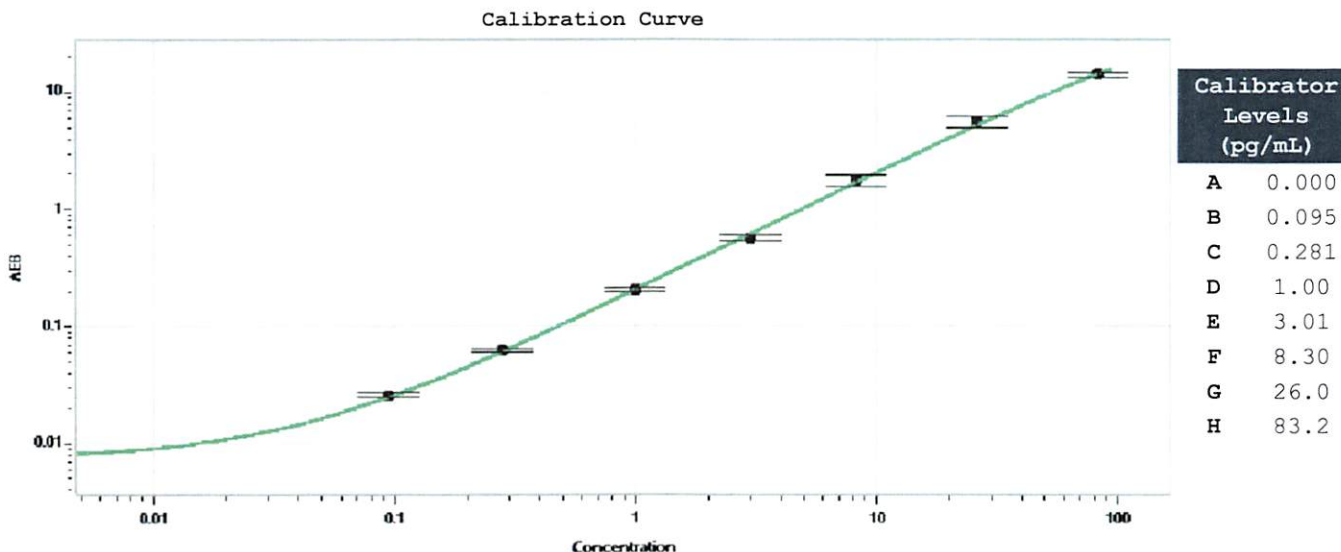
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101554	110905
Detector Reagent	101553	110909
SBG Reagent	101555	110910
Sample Diluent	101556	110911
Calibrators	101559	112004
Control 1	101560	112005
Control 2	101561	112005
SBG Diluent (SR-X) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	112005	1.15	0.643-1.26
Control 2	112005	58.3	43.5-76.4
Panel 2	916902	1.24	0.635-1.32
Panel 3	916903	27.3	13.0-36.2
Panel 4	916904	78.5	37.0-118

* 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 12 May 2021
Signature/Date