

Product Number:	103186
Lot Number:	502845
Expiration:	30-Mar-2022
Platform(s):	HD-1/ HD-X

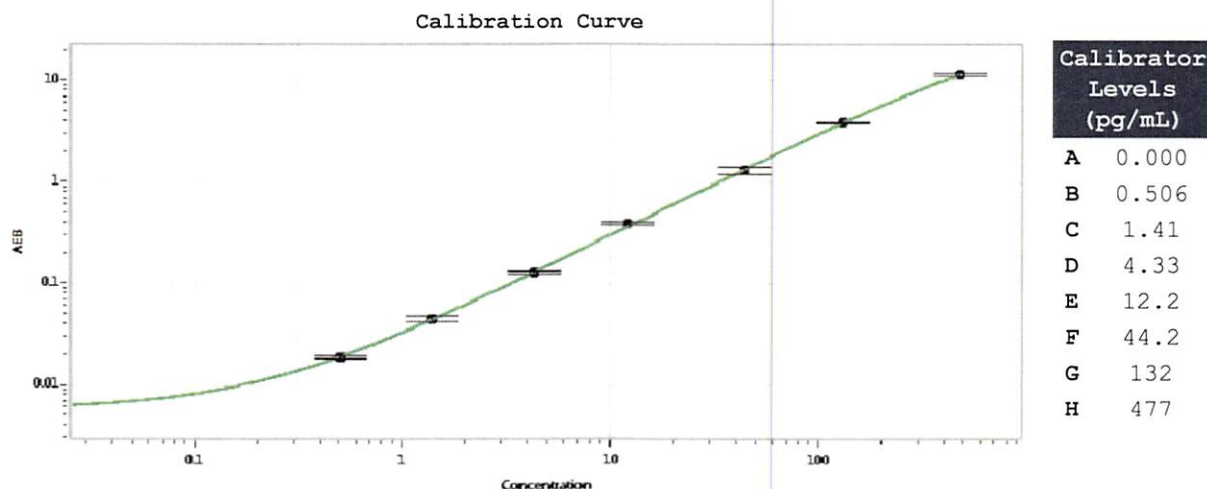
Component	Part Number	Lot Number
Bead Reagent	102246	109016
Detector Reagent	102248	108913
SBG Reagent	102250	108909
Sample Diluent	102252	108914
Calibrators	102255	109806
Control 1	102256	109807
Control 2	102257	109807
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



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

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	109807	4.90	3.44-5.87	
Control 2	109807	252	174-298	
Panel 1	008501	14.1	8.59-20.5	
Panel 2	008501	12.0	7.39-15.1	
Panel 3	008501	69.3	43.8-91.8	
Panel 4	009401	203	144-253	



Review/Approval

Junpil Park

QC Supervisor

Name

Title

Signature/Date

[Signature] 14 APR 2021