

Product Number:	101478
Lot Number:	503369
Expiration:	27-Feb-2023
Platform(s):	HD-1/HD-X & SR-X



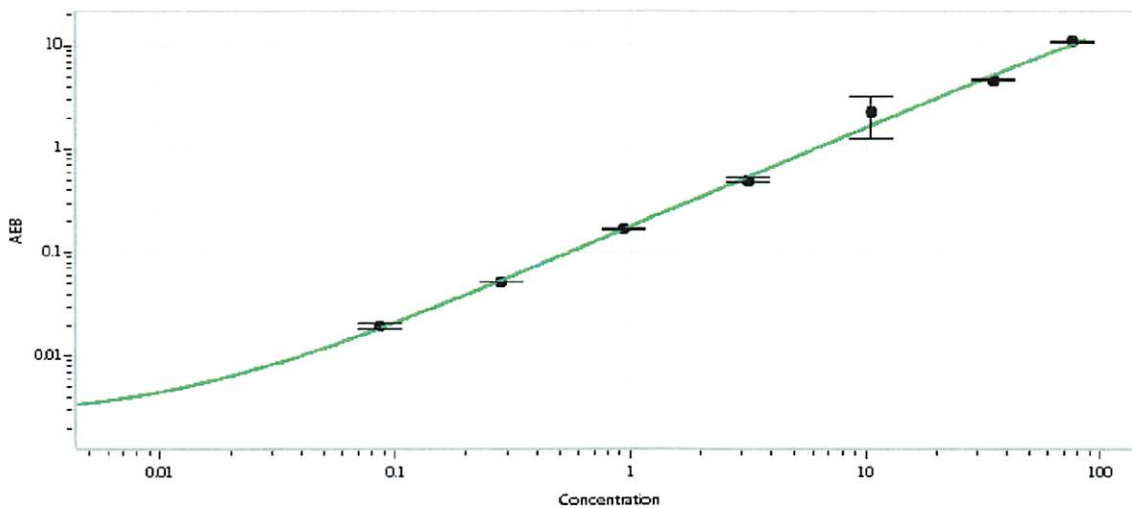
Data below represents results generated on the Simoa® HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101480	209605
Detector Reagent	101481	209807
SBG Reagent	101482	209808
Sample Diluent	101483	209810
Calibrators	100758-100765	210006
Control 1	100768	210007
Control 2	100769	210007
RGP Reagent	103159	N/A1

¹ RGP is not Kit Lot Specific

Release Materials*				Ranges
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	210007	2.98	2.31-3.56	
Control 2	210007	68.5	51.3-76.9	
Panel 2	122912	4.772	3.63-5.85	
Panel 3	122912	57.27	41.8-69.7	

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.087
C	0.281
D	0.947
E	3.18
F	10.6
G	35.1
H	76.1

Review/Approval: Shivani Goel Name, QC Manager Title, Shivani Goel 06 May 2022 Signature/Date