

Product Number:	103186
Lot Number:	502212
Expiration:	17-Jan-2021
Platform(s):	HD-1/HD-x SRx

Component	Part Number	Lot Number
Bead Reagent	102246	002705
Detector Reagent	102248	002701
SBG Reagent	102250	002711
Sample Diluent	102252	002704
Calibrators	102255	935001
Controls	102256 - 102257	935701
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

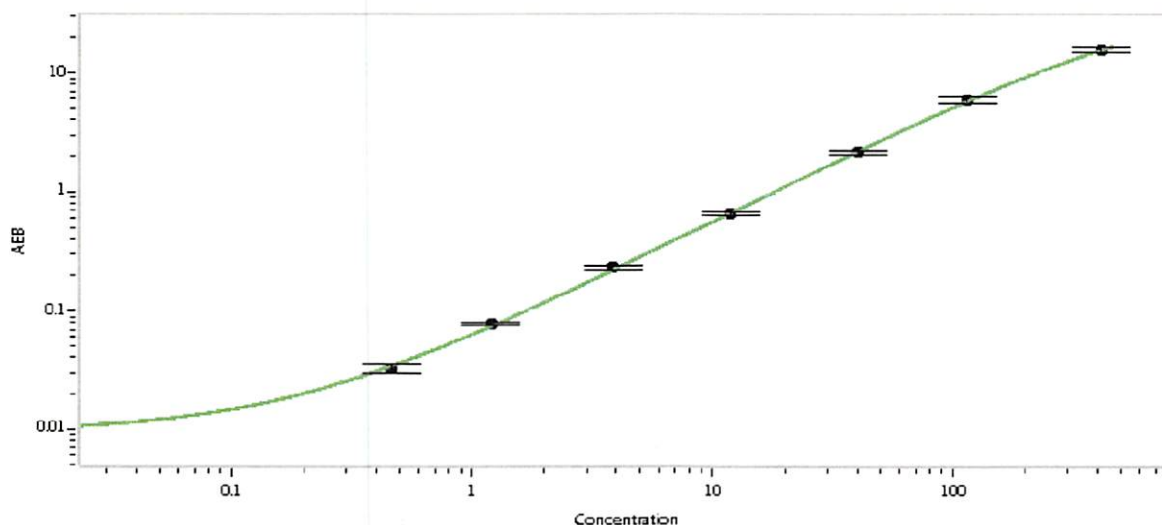


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	935701	4.79	3.07-5.31
Control 2	935701	156	132-198
Panel 1	922601	21.0	14.1-21.1
Panel 2	922602	7.24	4.94-7.41
Panel 3	922603	41.7	28.4-42.6
Panel 4	922604	171	129-204

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

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.461
C	1.20
D	3.87
E	11.8
F	40.2
G	115
H	415

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	09 Feb 2020
	Name	Title	Signature	Date