

Product Number:	103180
Lot Number:	502295
Expiration:	26-Feb-2021
Platform(s):	HD-1/HD-x

Component	Part Number	Lot Number
Bead Reagent	102246	006304
Detector Reagent	102248	006206
SBG Reagent	102250	006205
Sample Diluent	102252	006310
Calibrators	102255	924701
Controls	102256 - 102257	924803
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



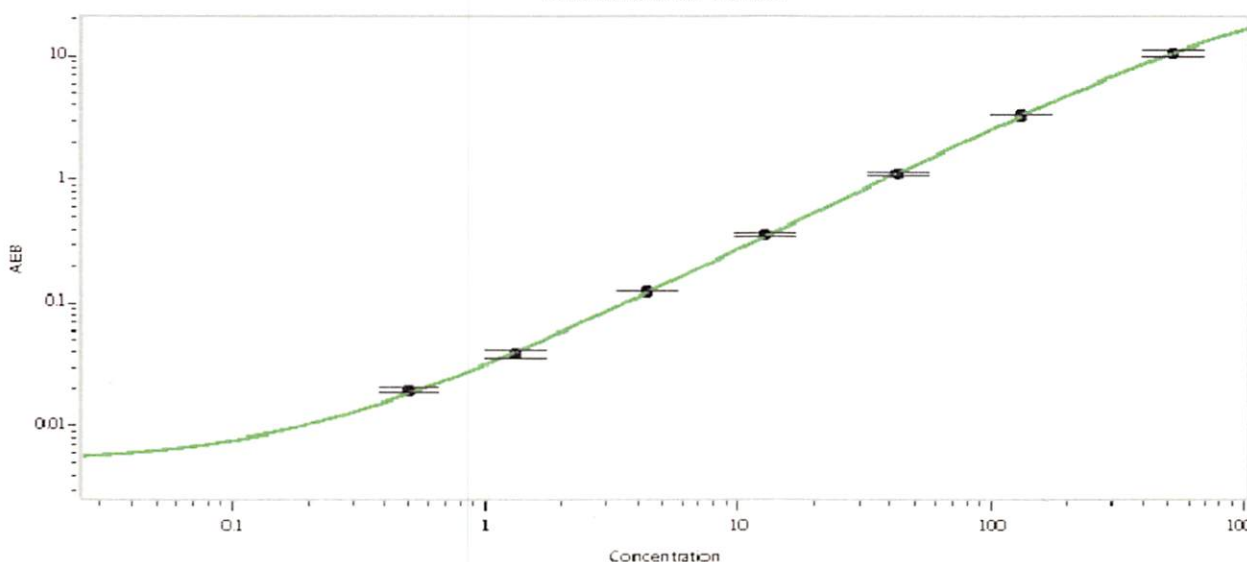
Data below represents results generated on the Simoa™ HD-X Analyzer.

Release Materials*

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	924803	3.46	2.50-4.49
Control 2	924803	143	116-173
Panel 1	922601	15.7	14.1-21.1
Panel 2	922602	5.48	4.94-7.41
Panel 3	922603	36.1	28.4-42.6
Panel 4	922604	156	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	0.503
C	1.33
D	4.32
E	12.8
F	42.3
G	130
H	521

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler

23 Mar 2020

Name

Title

Signature/Date