

Product Number:	100794
Lot Number:	502127
Expiration:	13-Jun-2020
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	100781	916406
Detector Reagent	100782	916409
SBG Reagent	100783	916506
Sample Diluent	103202	928024
Calibrator Diluent	100793	930817
Calibrator Concentrate	100780	932311
SBG Diluent (SRX)†	103158	N/A1
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

† Specific to use on the SR-X platform - refer to the SR-X Kit Instructions (KI-0085).



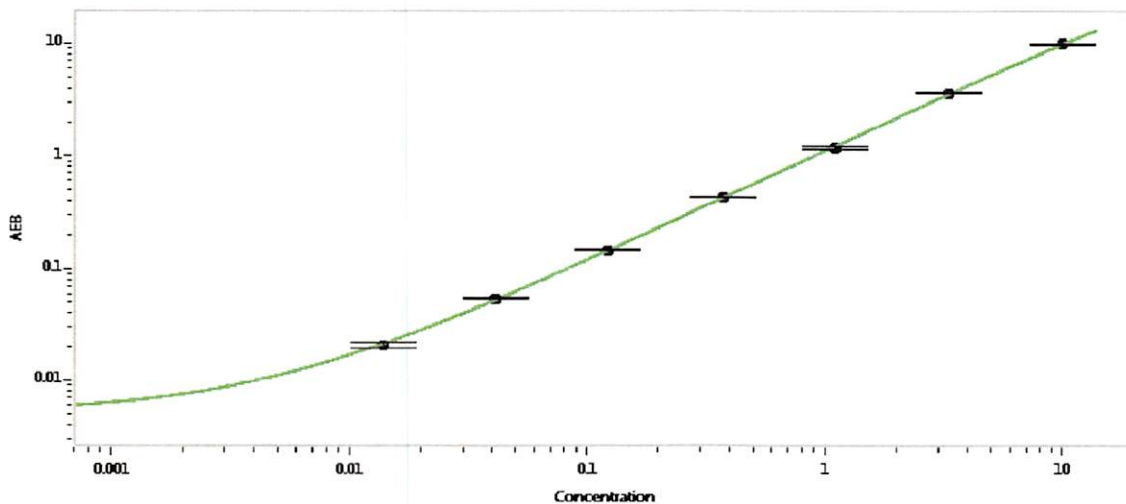
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-062316-02	9.61	7.17-10.8
Control 2	MF-062316-02	0.791	0.606-0.910
Panel 1	MF-110316-05	3.89	3.54-5.31
Panel 2	MF-110316-06	45.8	39.6-59.3
Panel 3	MF-110316-07	3.82	3.16-4.74

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

Calibrator Concentrate Value Assignment (pg/mL)	134
---	-----

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.014
C	0.041
D	0.123
E	0.370
F	1.11
G	3.33
H	10.0

Review/Approval

Patrizia Stadler Assoc. Director, QC
Name Title

P. Stadler 16 Dec 2019
Signature/Date