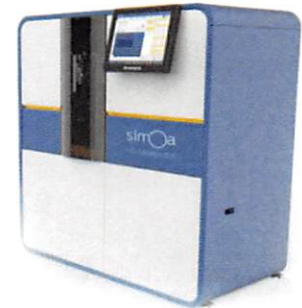


Product Number:	100794
Lot Number:	501689
Expiration:	25-Jan-2020
Platform(s):	HD-1 and SR-X



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

Component	Part Number	Lot Number
Bead Reagent	100781	905704
Detector Reagent	100782	905609
SBG Reagent	100783	905610
Sample Diluent	103202	902928
Calibrator Diluent	100793	904414
Calibrator Concentrate	100780	80254
SBG Diluent (SRX) †	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> Reagents are not Kit Lot Specific.

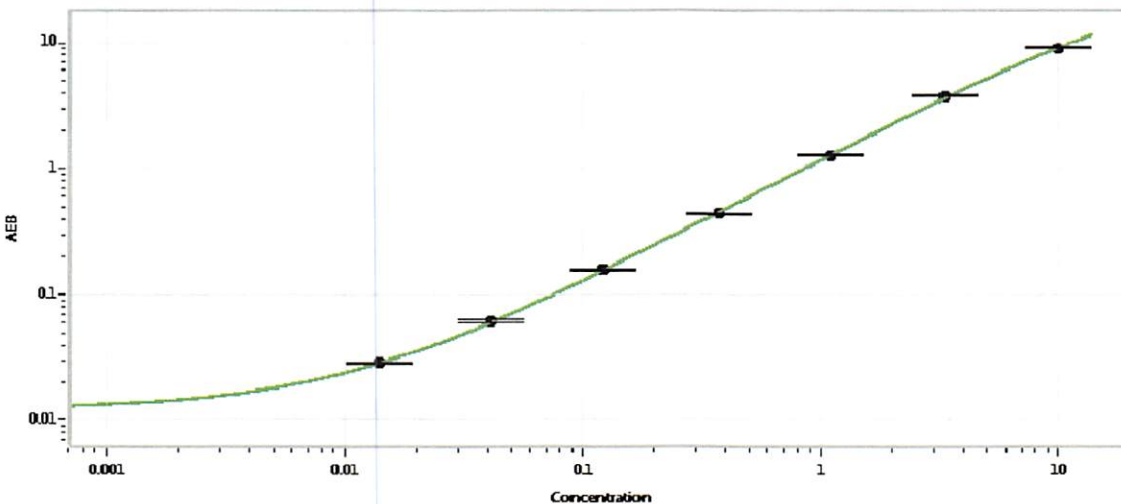
† Specific to use on the SR-X platform - refer to the SR-X Kit Instructions (KI-0085).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-062316-02	8.61	7.11-10.7
Control 2	MF-062316-02	0.764	0.595-0.893
Panel 1	MF-110316-05	3.78	3.61-5.41
Panel 2	MF-110316-06	44.2	40.0-60.0
Panel 3	MF-110316-07	3.49	3.14-4.71

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

Calibrator Concentrate Value Assignment (pg/mL)	829
---	-----

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

*P. Stadler* 04 March 2019

Name

Title

Signature/Date