

Product Number:	100794
Lot Number:	501203
Expiration:	5-Oct-2018
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	80952
Detector Reagent	100782	80950
SBG Reagent	100783	80951
Sample Diluent	103202	80864
Calibrator Diluent	100793	80793
Calibrator Concentrate	100780	80254
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific.

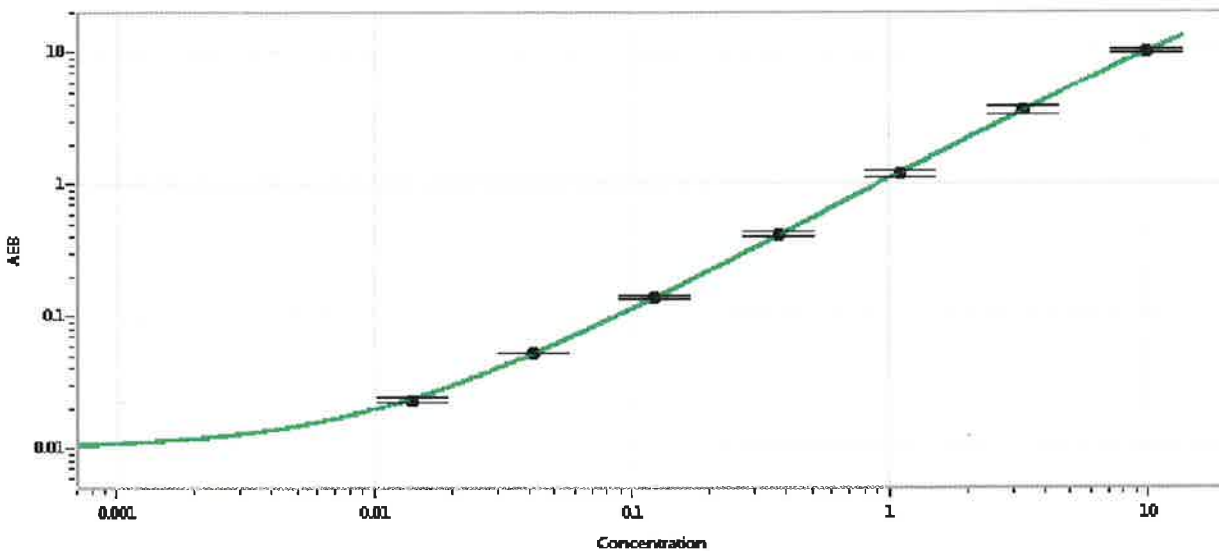
† Specific to use on the SRX platform – refer to the SRX Kit Instructions (KI-0085).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-062316-02	9.88	4.95-10.9
Control 2	MF-062316-02	0.800	0.493-0.877
Panel 1	MF-110316-05	3.58	2.54-6.00
Panel 2	MF-110316-06	35.4	32.6-54.9
Panel 3	MF-110316-07	2.66	1.90-5.31

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

Calibrator Concentrate Value Assignment (pg/mL)	746
---	-----

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

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 13 APR 2018
Signature/Date