

Product Number:	102732
Lot Number:	501514
Expiration:	20-May-2019
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	102722	833002
Detector Reagent	102724	833001
SBG Reagent	102726	833011
Sample Diluent	102728	832410
Calibrator Diluent	102731	826811
Calibrator Concentrate	100658	833101
RGP Reagent	103159	N/A ¹

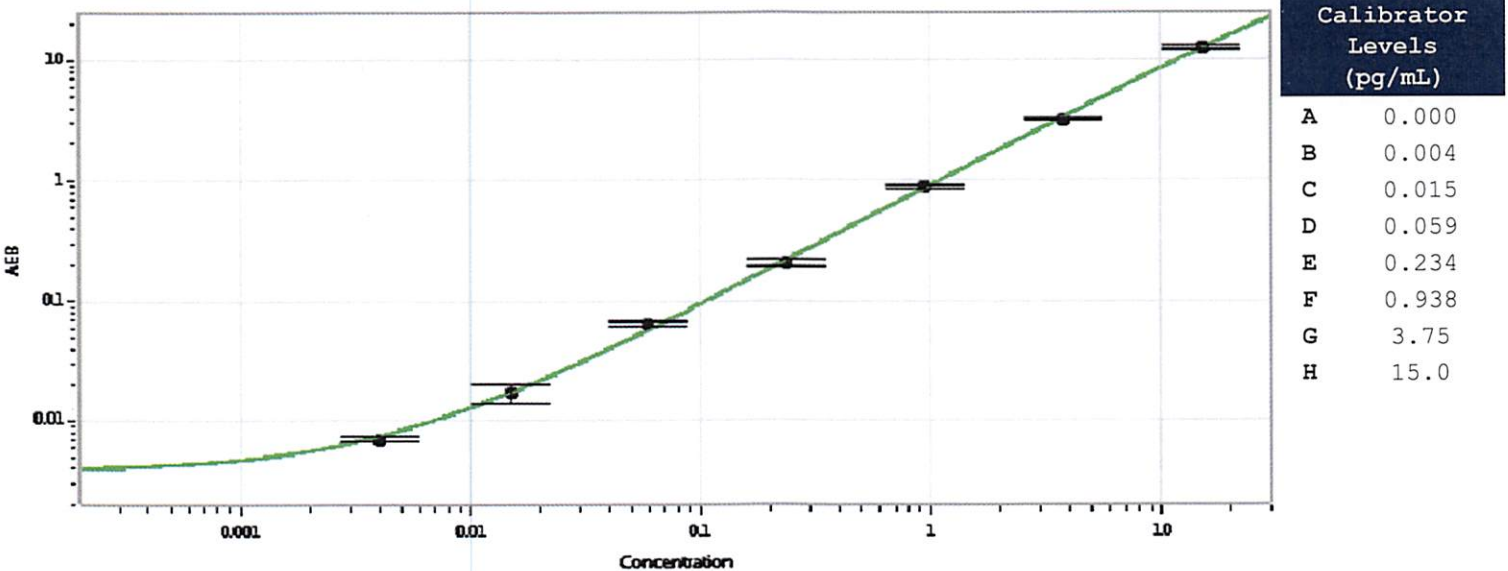
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-061217-01	0.191	0.134-0.289
Control 2	MF-061217-01	3.71	2.44-6.35
Panel 1	MF-061317-03	0.050	0.037-0.088
Panel 2	MF-061317-04	0.276	0.191-0.490
Panel 3	MF-061317-05	4.41	3.18-7.89

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

Calibrator Concentrate Value Assignment (pg/mL)	202
---	-----

Calibration Curve



Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date