

Product Number: 102732
Lot Number: 501605
Expiration: 17-Jan-2020
Platform(s): HD-1 & SR-X



Component	Part Number	Lot Number
Bead Reagent	102722	902209
Detector Reagent	102724	901902
SBG Reagent	102726	901901
Sample Diluent	102728	901905
Calibrator Diluent	102731	901906
Calibrator Concentrate	100658	833101
RGP Reagent	103159	N/A ¹

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

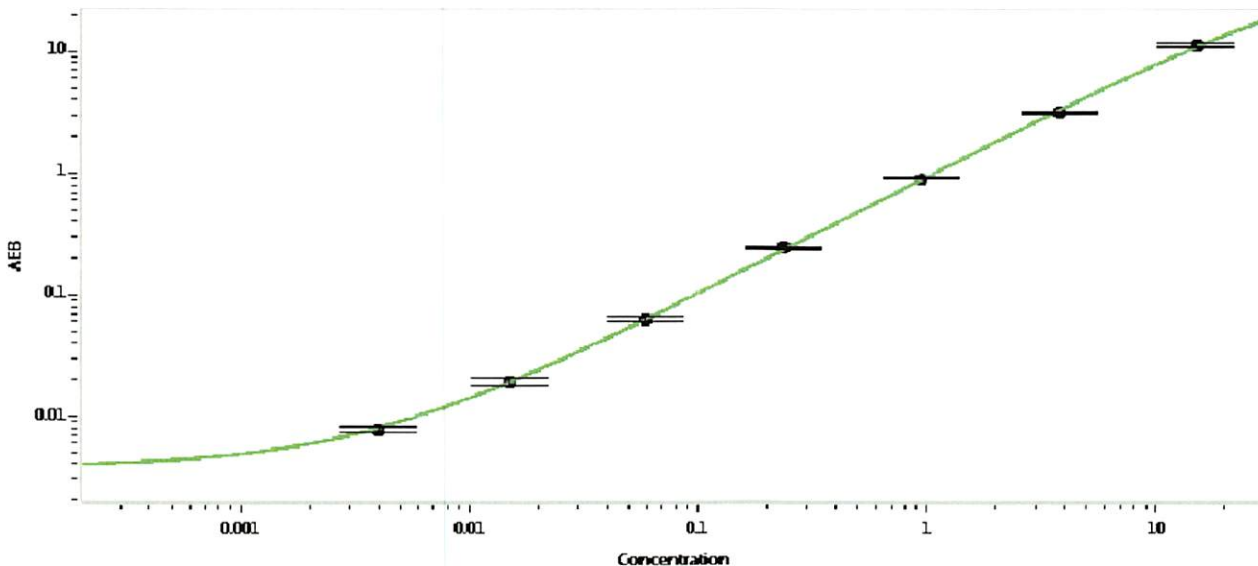
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-061217-01	0.173	0.134-0.289
Control 2	MF-061217-01	3.59	2.44-6.35
Panel 1	MF-061317-03	0.042	0.037-0.088
Panel 2	MF-061317-04	0.260	0.191-0.490
Panel 3	MF-061317-05	4.12	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



Calibrator Levels (pg/mL)	
A	0.000
B	0.004
C	0.015
D	0.059
E	0.234
F	0.938
G	3.75
H	15.0

Review/Approval: Patrizia Stadler Associate Director, QC
 Name: Patrizia Stadler Title: QC Signature/Date: p. Stadler 23 Jun 2019