

Product Number:	102329
Lot Number:	501820
Expiration:	30-Mar-2020
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	102322	912407
Detector Reagent	102324	912406
SBG Reagent	102325	912608
Sample Diluent	102326	908909
Calibrator Diluent	102327	912014
Calibrator Concentrate	100849	80923
RGP Reagent	103159	N/A ¹

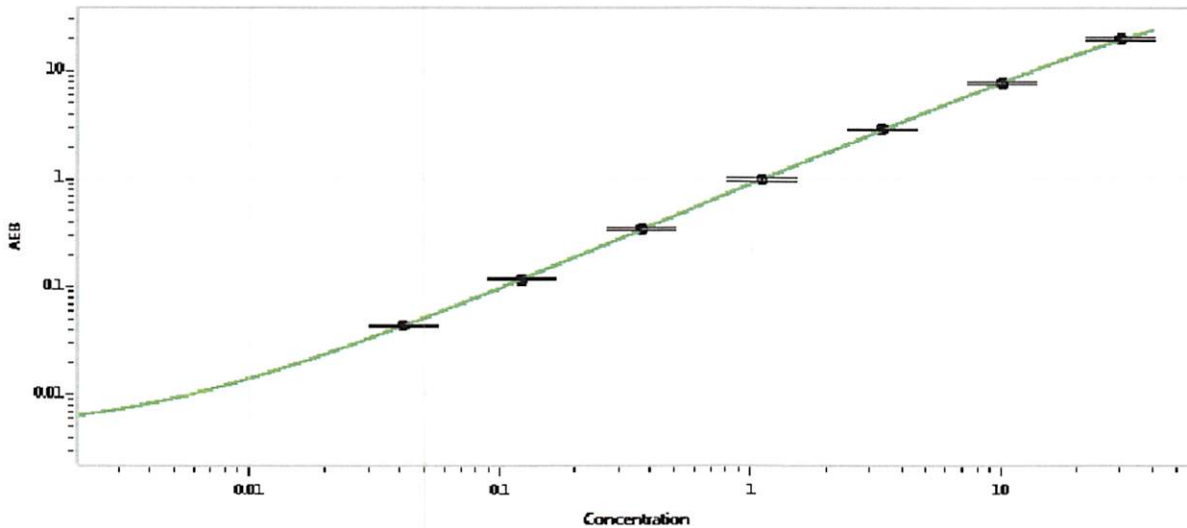
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-092716-01	0.181	0.108-0.300
Control 2	MF-092716-01	24.5	17.8-37.9
Panel 1	MF-092116-03	0.093	0.065-0.137
Panel 2	MF-092116-04	0.089	0.069-0.135
Panel 3	MF-092116-05	15.7	9.71-25.5

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

Calibrator Concentrate Value Assignment (pg/mL)	647 pg/mL
---	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval: Patrizia Stadler QC Assoc. Director
 Name: Patrizia Stadler Title: QC Assoc. Director Signature/Date: P. Stadler 09 MAY 19