

Product Number:	101212
Lot Number:	502327
Expiration:	19-Dec-2020
Platform (s):	HD-1/HDX

Component	Part Number	Lot Number
Bead Reagent	101198	010506
Detector Reagent	101200	014202
SBG Reagent	101211	010302
Sample Diluent	101209	935314
Calibrator Diluent	101204	010304
Calibrator Concentrate	101202	010701
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



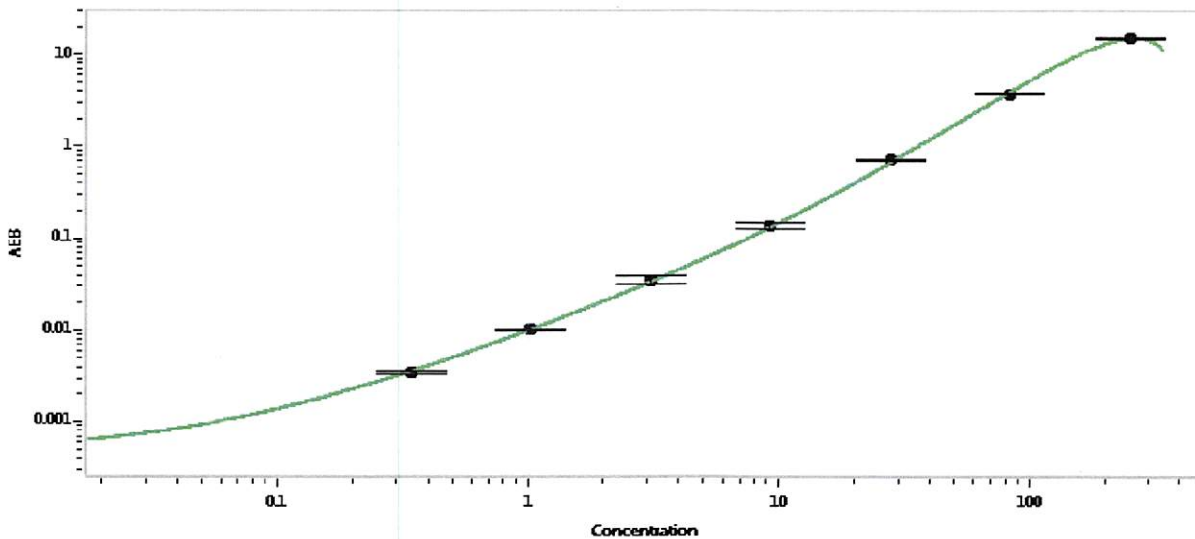
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	828802	81.7	55.1-82.6
Control 2	828802	527	354-532
Panel 1	MF-052318-05	200	141-211
Panel 2	MF-052318-06	52.3	35.9-53.8
Panel 3	MF-052318-07	97.5	72.7-109

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

Calibrator Concentrate Value Assignment (pg/mL)	3768
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.343
C	1.03
D	3.09
E	9.26
F	27.8
G	83.3
H	250

Review/Approval

Patrizia Stadler QC Assoc Director

P. Stadler

01 Jun 2020

Name

Title

Signature/Date