

Product Number:	103186
Lot Number:	501706
Expiration:	16-Feb-2020
Platform(s):	HD-1 & SR-X

Component	Part Number	Lot Number
Bead Reagent	102246	906503
Detector Reagent	102248	904703
SBG Reagent	102250	904702
Sample Diluent	102252	904701
Reference Calibrators	102255	906401
Controls	102256 - 102257	906402
RGP Reagent	103159	N/A ¹

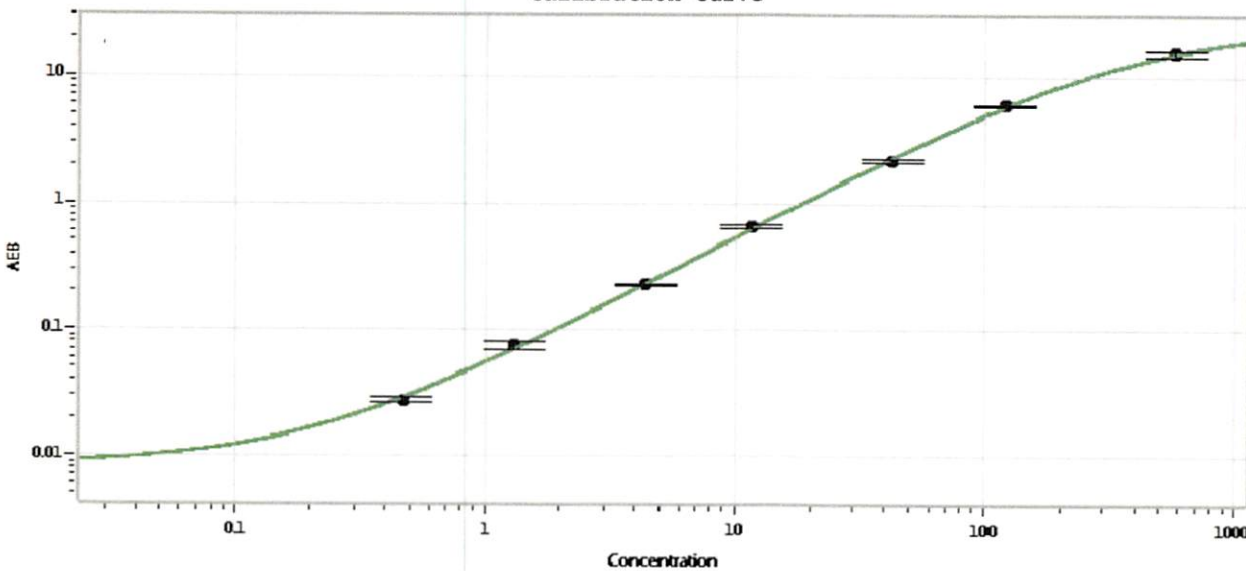
¹ RGP is not Kit Lot Specific



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

	Release Materials*			* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	906402	3.70	3.06-4.94	
Control 2	906402	134	129-193	
Panel 1-A	823301	4.85	4.08-6.68	
Panel 1-B	823302	5.50	4.90-7.61	
Panel 2	823601	65.1	59.3-101	
Panel 3	823602	254	236-402	

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.466
C	1.42
D	4.79
E	14.0
F	47.6
G	142
H	472

Review/Approval

Patrizia Stadler

QC Assoc. Director

P. Stadler

18 OCT 2019

Name

Title

Signature/Date