

Product Number:	103400
Lot Number:	501893
Expiration:	16-Feb-2020
Platform(s):	HD-1/HD-X and SR-X

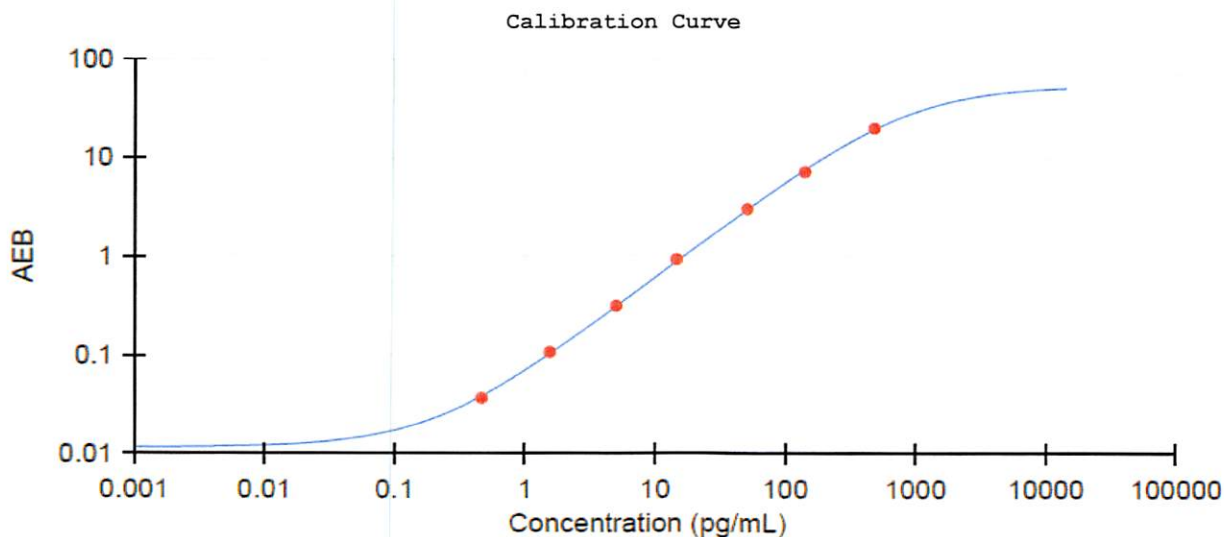


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102246	905611
Detector Reagent	102248	904703
SBG Reagent	102250	904702
Sample Diluent	103399	905601
Calibrators	102255	916301
Controls	102256-102257	916401
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific.

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	916401	4.18	3.30-5.02	
Control 2	916401	157	133-199	
Panel 1	823301	5.53	4.08-6.68	
Panel 2	823302	6.66	4.90-7.61	
Panel 3	823601	88.2	59.3-101	
Panel 4	823602	328	236-402	



Calibrator Levels (pg/mL)	
A	0.000
B	0.468
C	1.55
D	5.04
E	14.6
F	50.8
G	141
H	478

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>p.stadler</i>	08 JUL 2019
	Name	Title	Signature/Date	