

Product Number:	103377
Lot Number:	502146
Expiration:	29-Oct-2020
Platform(s):	HD-1/HD-X and SR-X



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

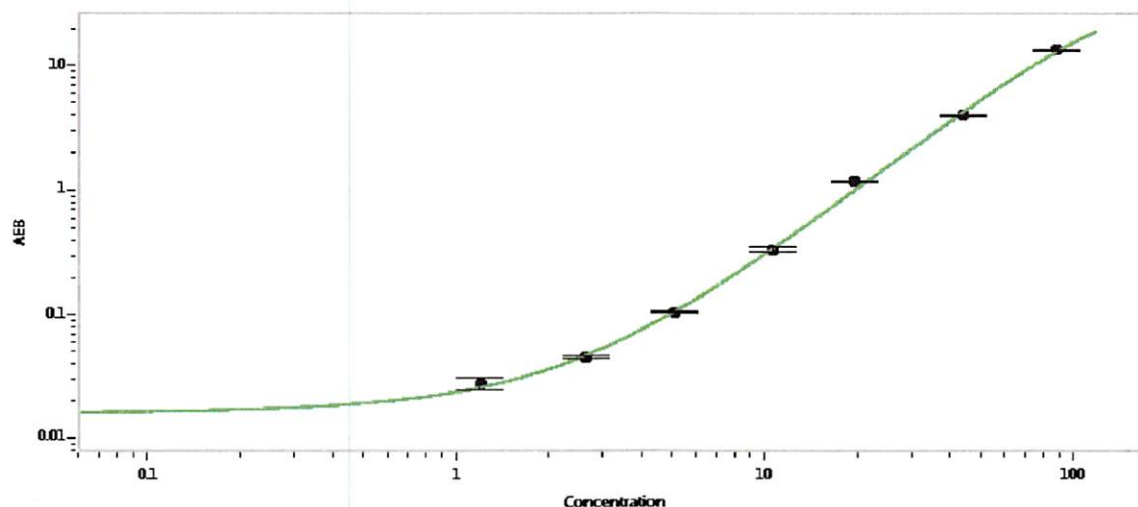
Component	Part Number	Lot Number
Bead Reagent	103364	934019
Detector Reagent	103366	933107
SBG Reagent	103368	932917
Sample Diluent	103370	933001
Calibrators	103374	901102
Controls	103375/103376	901103/901801
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific.

† Specific to use on the SRX platform - refer to the SRX Kit Instructions (KI-0045).

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	901103	20.6	16.1-24.2	
Control 2	901801	155	125-188	
Panel 1	900701	33.4	29.4-44.1	
Panel 2	900702	42.1	35.0-52.6	
Panel 3	900703	25.0	20.6-30.9	

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.20
C	2.64
D	5.08
E	10.6
F	19.7
G	44.2
H	88.0

Review/Approval

Patrizia Stadler Assoc. Director, QC
Name Title

P. Stadler 21 Dec 2019
Signature/Date