

Product Number:	102318
Lot Number:	501252
Expiration:	29-Sep-2018
Platform(s):	HD-1



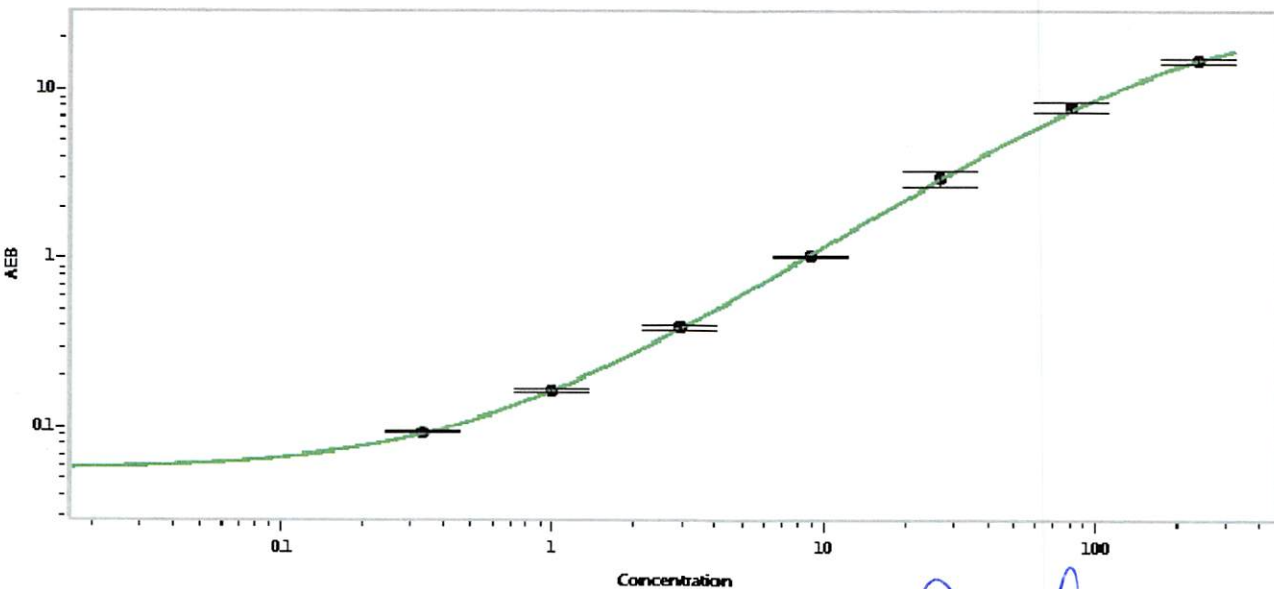
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102309	MF-042318-03
Detector Stock	102310	MF-051618-01
Calibrator Concentrate	102314	MF-092916-02
Calibrator Diluent	102311	MF-041718-03
Sample Diluent	102312	MF-041718-02

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Concentrate	101957
RGP Reagent	103159

Calibrator Concentrate Target Value	4,800 pg/mL
--	-------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.329
C	0.988
D	2.96
E	8.89
F	26.7
G	80.0
H	240

James Drummey
Senior Director, Manufacturing Operations
On behalf of:

[Handwritten Signature]
May 18, 2018

Review/Approval	Sivasankari Murugan	QC Supervisor	
	Name	Title	Signature/Date