

Product Number:	102713
Lot Number:	501608
Expiration:	10-Apr-2019
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	102706	812002
Detector Stock	102707	810602
Calibrator Concentrate	102710	MF-041618-05
Calibrator Diluent	102708	MF-041018-15
Sample Diluent	102709	MF-041018-16

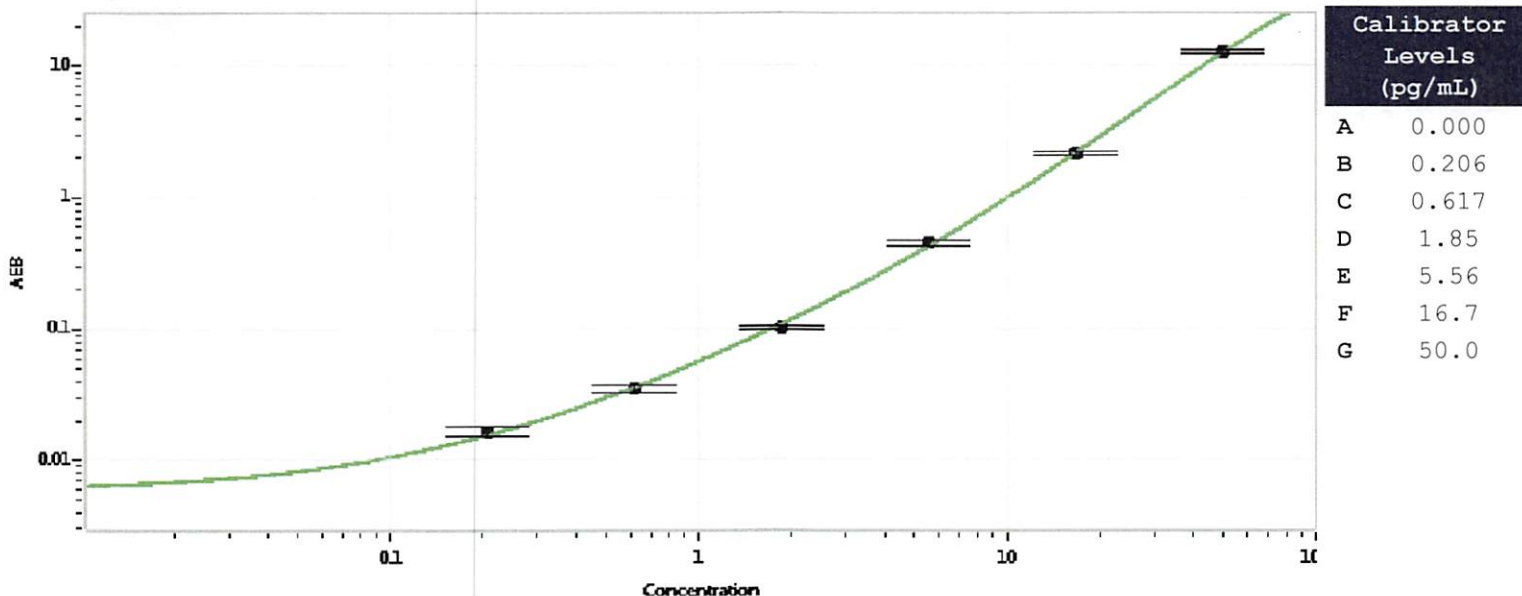


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Generic Components - Included With Kit But Not Lot Number Specific		
Discovery Bead Diluent		101954
Discovery Detector Diluent		101955
Discovery SBG Diluent		101956
Discovery SBG Stock		103297
RGP Reagent		103159

Calibrator Concentrate Target Value	2 ng/mL
-------------------------------------	---------

Calibration Curve



Review/Approval

Patrizia Stadler QC Assoc. Director

P. Stadler 28 Jan 2019

Name

Title

Signature/Date