

Product Number:	102336
Lot Number:	501819
Expiration:	1-Oct-2019
Platform(s):	HD-1



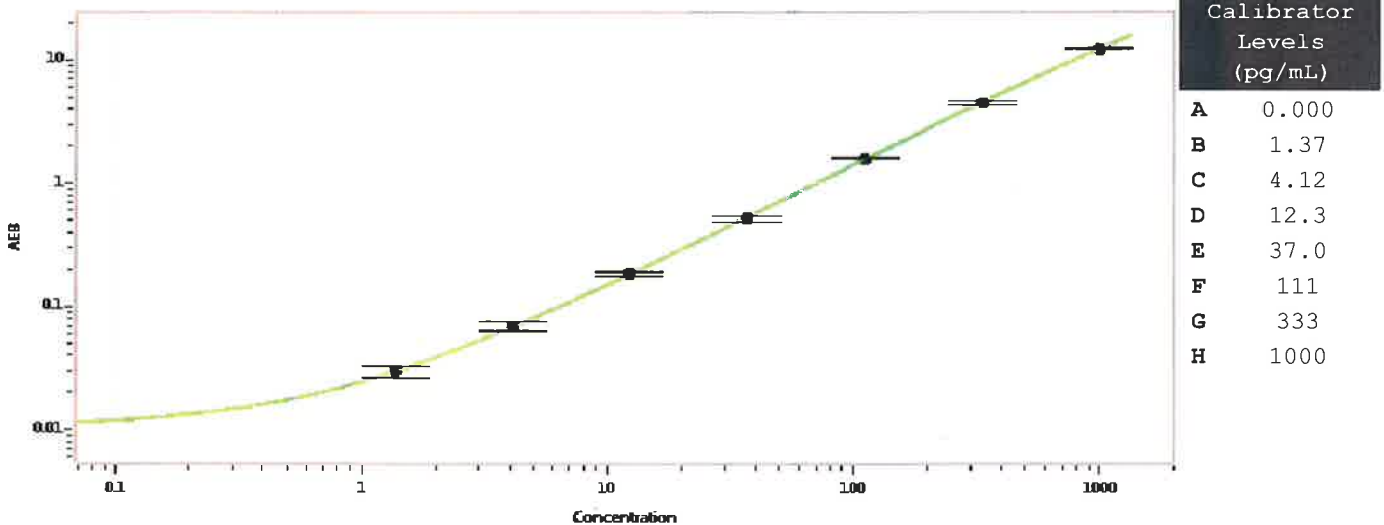
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102330	912002
Detector Stock	102331	912004
Calibrator Concentrate	102335	912103
Calibrator Diluent	102332	827417
Sample Diluent	102333	912003

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	20 ng/mL
-------------------------------------	----------

Calibration Curve



Review/Approval

Patrizia Stadler  
Name

QC Assoc. Director  
Title

*P. Stadler 13 May 2019*  
Signature/Date