

Product Number:	102583
Lot Number:	502205
Expiration:	21-Nov-2020
Platform(s):	HD-1/HD-X



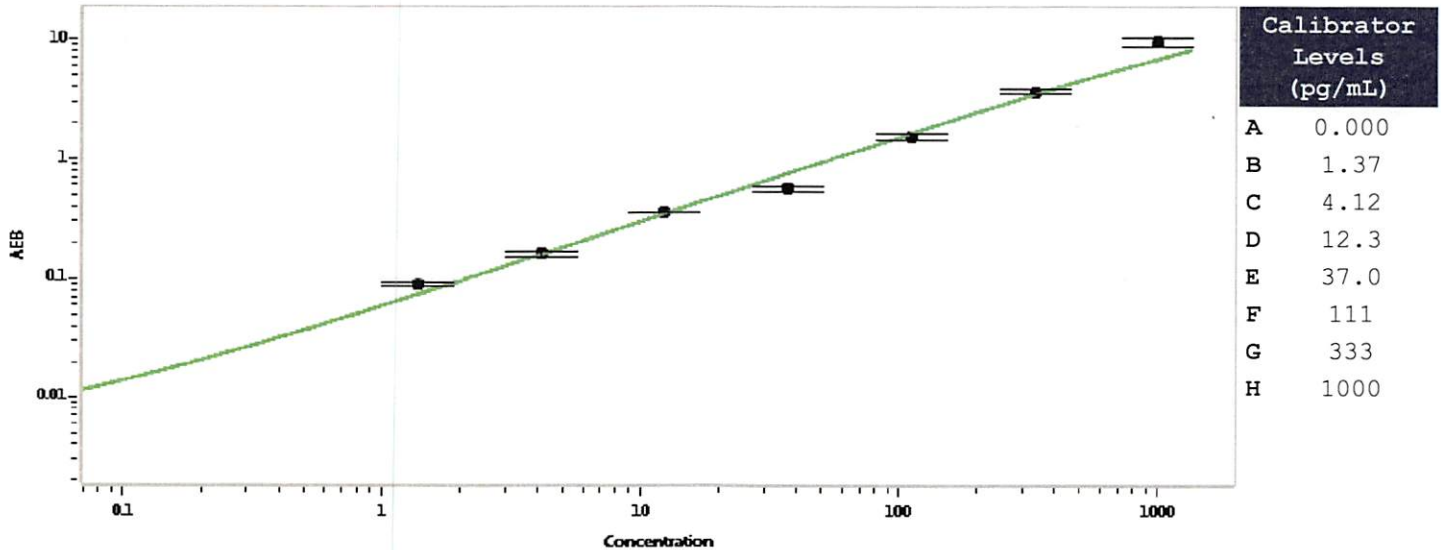
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102576	002209
Detector Stock	102577	002205
Calibrator Concentrate	102580	912804
Calibrator Diluent	102578	932925
Sample Diluent	102579	933813

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	10 ng/mL
-------------------------------------	----------

Calibration Curve



Review/Approval

Patrizia Stadler

QC Asso. Director

*P. Stadler* 30 Jan 2020

Name

Title

Signature/Date