

Product Number: 102556
 Lot Number: 501817
 Expiration: 29-Mar-2020
 Platform(s): HD-1



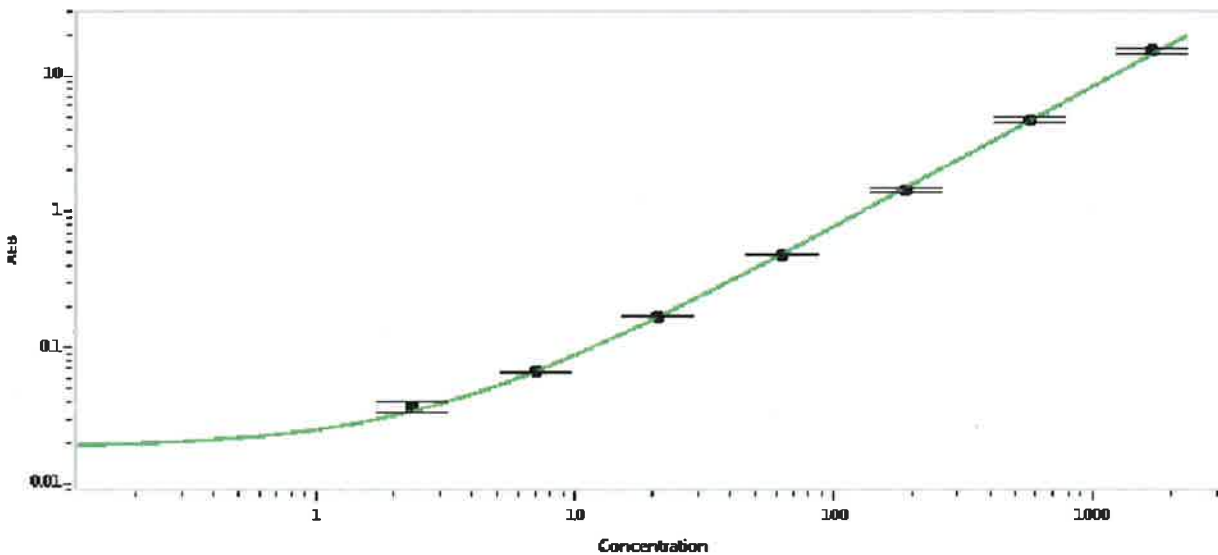
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102549	911909
Detector Stock	102550	911912
Calibrator Concentrate	102553	900707
Calibrator Diluent	102551	910308
Sample Diluent	102552	908905

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: 40 ng/mL

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.33
C	7.00
D	21.0
E	63.0
F	189
G	567
H	1700

Review/Approval: Patrizia Stadler QC Assoc. Director
 Name: Patrizia Stadler Title: QC Assoc. Director
 Signature/Date: Patrizia Stadler 14 MAY 2019