

Product Number:	102184
Lot Number:	502259
Expiration:	21-Nov-2020
Platform(s):	HD-1/ HD-X



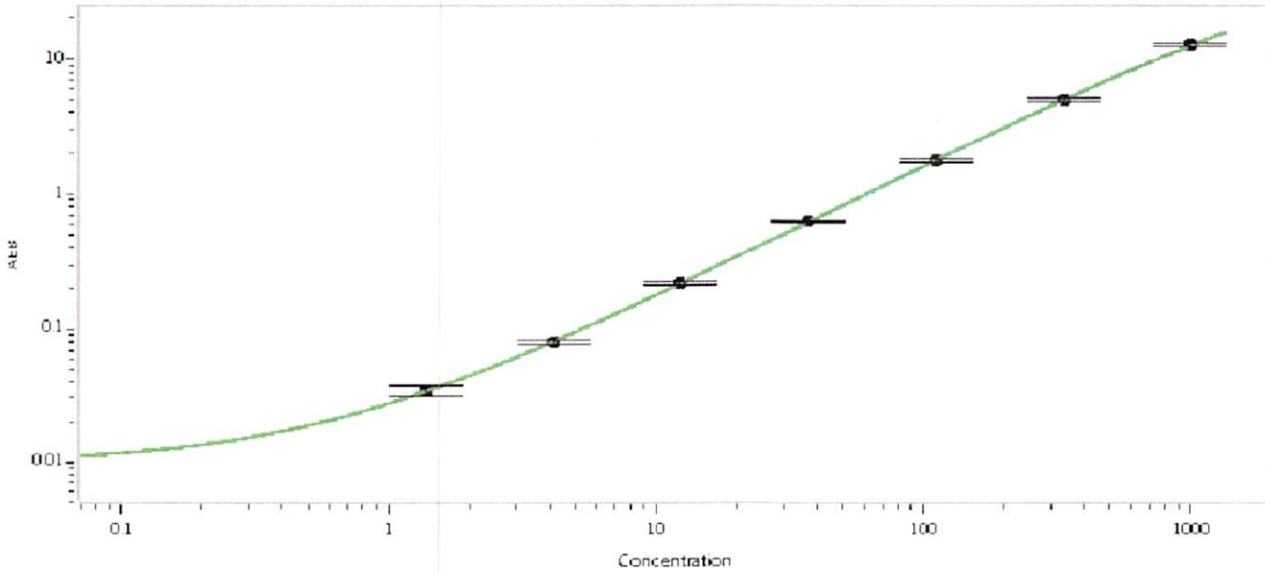
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102176	005109
Detector Stock	102177	005104
Calibrator Concentrate	102180	005605
Calibrator Diluent	102178	003514
Sample Diluent	102179	933820

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	1.37
C	4.12
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

p. Stadler 28 Feb 2020
Signature/Date