

Product Number: 102462
Lot Number: 501904
Expiration: 30-Mar-2020
Platform(s): HD-1/ HD-X

Component	Part Number	Lot Number
Bead Stock	102455	911208
Detector Stock	102456	910904
Calibrator Concentrate	102459	911304
Calibrator Diluent	102457	910903
Sample Diluent	102458	910902

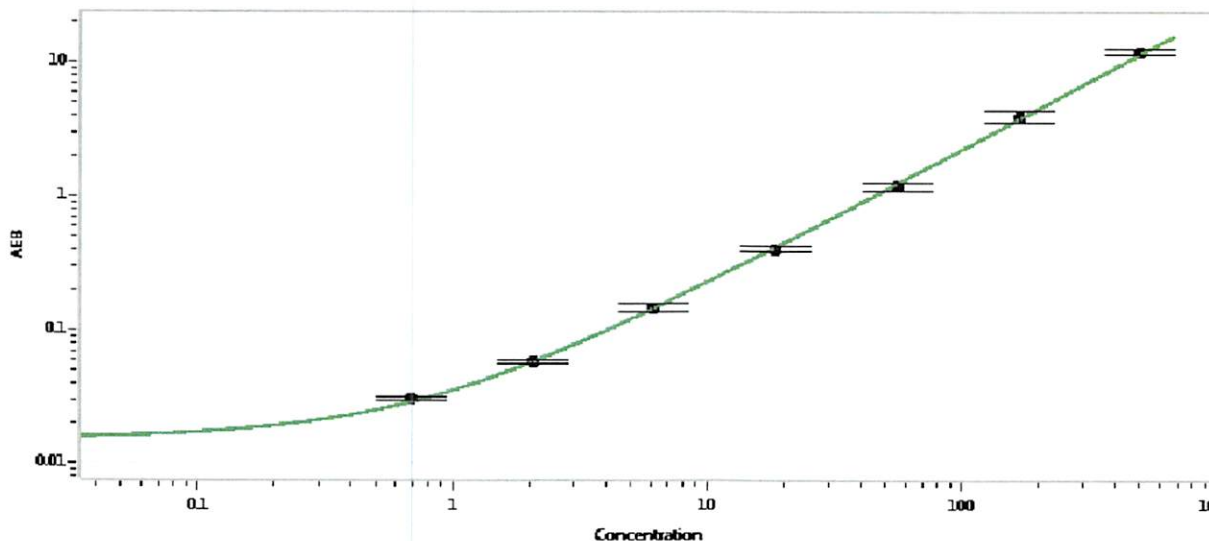


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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 10 ng/mL

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.686
C	2.06
D	6.17
E	18.5
F	55.6
G	167
H	500

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 09 Jul 2019
Signature/Date