

Product Number:	102233
Lot Number:	501759
Expiration:	1-Oct-19
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	102223	909317
Detector Stock	102224	909105
Calibrator Concentrate	102229	835303
Calibrator Diluent	102225	909104
Sample Diluent	102226	909103

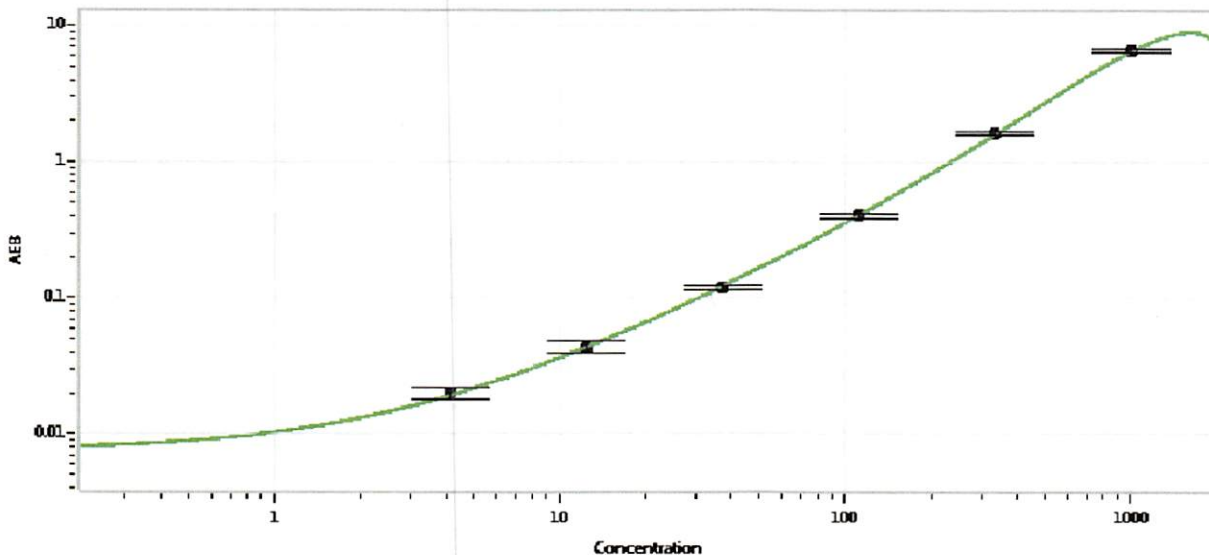


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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

Alpha-Synuclein Calibrator Concentrate Target Value	20 ng/mL
---	----------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	4.12
C	12.3
D	37.0
E	111
F	333
G	1000

Review/Approval

Patrizia Stadler Assoc. Director, QC

P. Stadler 10 Apr 2019

Name

Title

Signature/Date