

Product Number:	102233
Lot Number:	501558
Expiration:	10-Jun-19
Platform(s):	HD-1



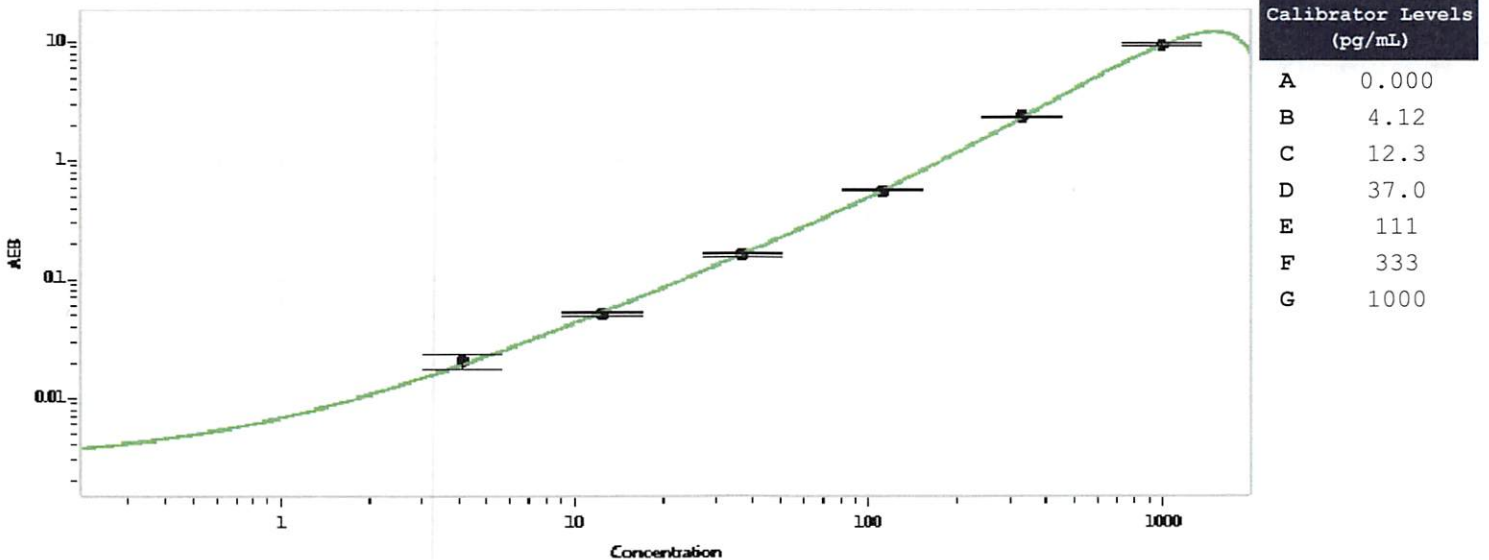
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102223	834804
Detector Stock	102224	834411
Calibrator Concentrate	102229	835002
Calibrator Diluent	102225	834410
Sample Diluent	102226	834409

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



Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date
20 DEC 2018