

Product Number:	102233
Lot Number:	501469
Expiration:	21-Nov-18
Platform(s):	HD-1



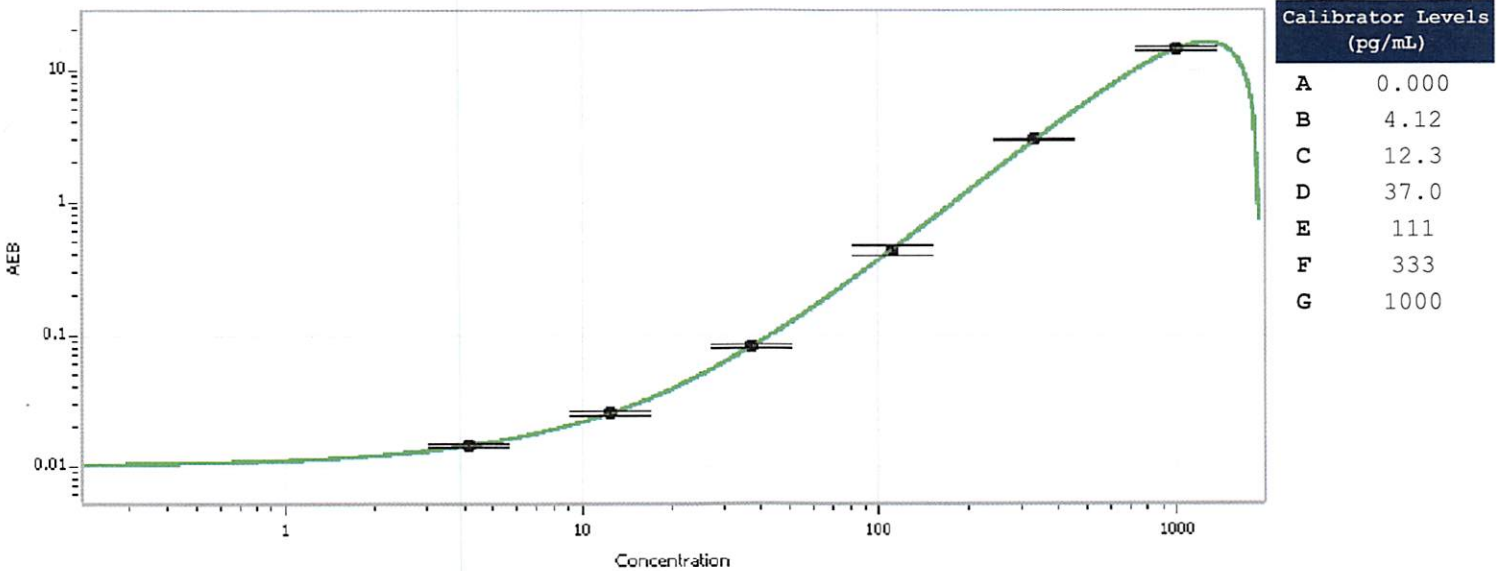
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102223	824003
Detector Stock	102224	822003
Calibrator Concentrate	102229	MF-112116-01
Calibrator Diluent	102225	822005
Sample Diluent	102226	822004

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] 24 Oct 2018
Signature/Date