

Product Number:	102233
Lot Number:	501390
Expiration:	21-Nov-18
Platform(s):	HD-1



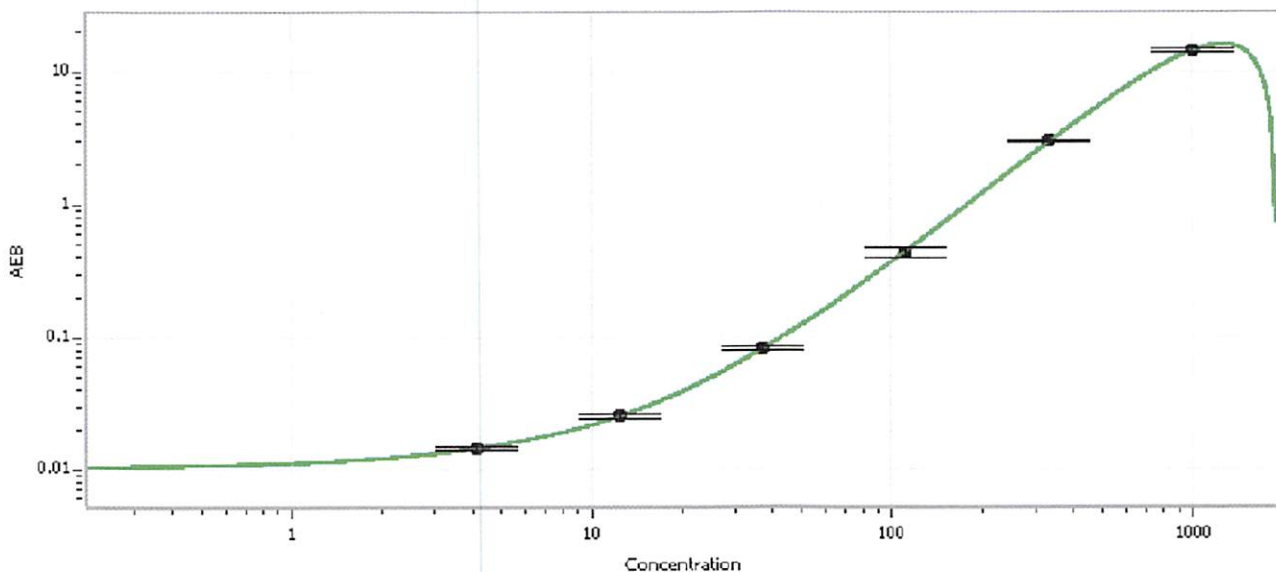
Data below represents results generated on the

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 Concentrate		101957
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

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date 31 AUG 2018