

Product Number:	102648
Lot Number:	502014
Expiration:	30-Apr-2020
Platform(s):	HD-1/HD-X & SR-X



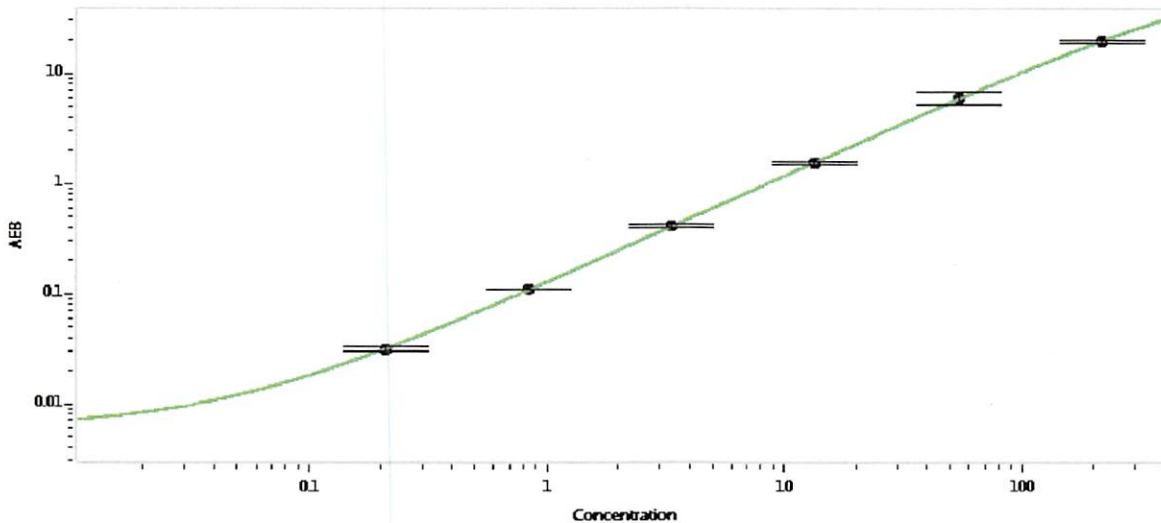
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102641	924609
Detector Stock	102642	924903
Calibrator Concentrate	102645	924602
Calibrator Diluent	102643	924007
Sample Diluent	102644	912024

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	3,000 pg/mL
-------------------------------------	-------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.21
C	0.84
D	3.36
E	13.4
F	53.8
G	215

Review/Approval

Patrizia Stadler  
Name

Assoc. Director, QC  
Title

*P. Stadler 13 Sep 2019*  
Signature/Date