

Product Number:	102648
Lot Number:	501627
Expiration:	19-Sep-2019
Platform(s):	HD-1

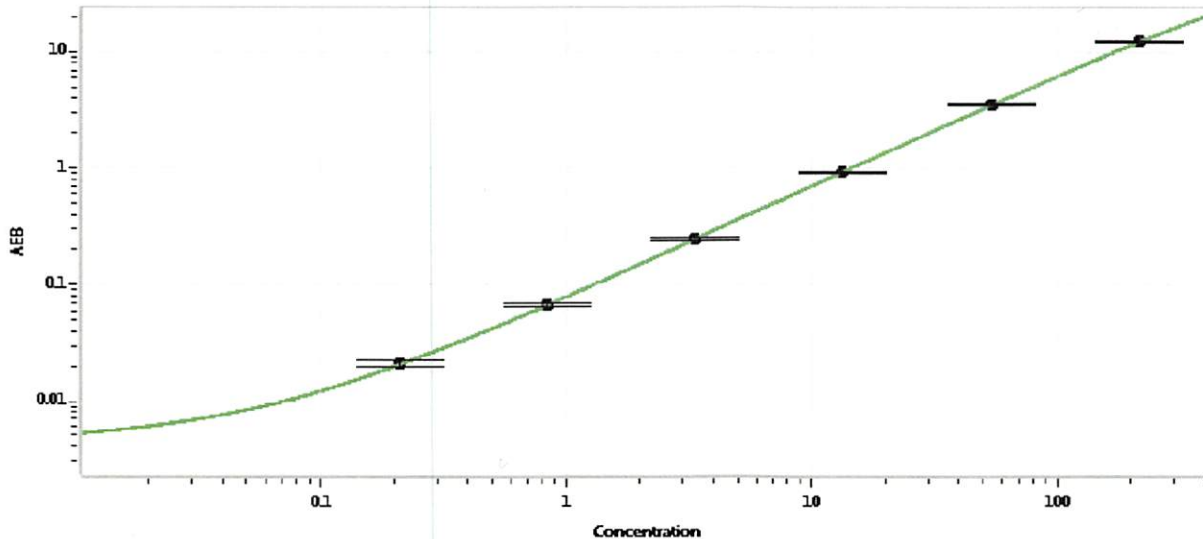


Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102641	901810
Detector Stock	102642	901908
Detector Diluent	102734	901907
Calibrator Concentrate	102645	902915
Calibrator Diluent	102643	901014
Sample Diluent	102644	826227

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 ng/ml
-------------------------------------	---------



Calibrator Levels (pg/mL)	
A	0.000
B	0.210
C	0.840
D	3.36
E	13.4
F	53.8
G	215

Review/Approval

Patrizia Stadler

QC Assoc. Director

*P. Stadler* 01 Feb 2019