

<b>Product Number:</b>	102382
<b>Lot Number:</b>	501635
<b>Expiration:</b>	10-Dec-2019
<b>Platform(s):</b>	HD-1



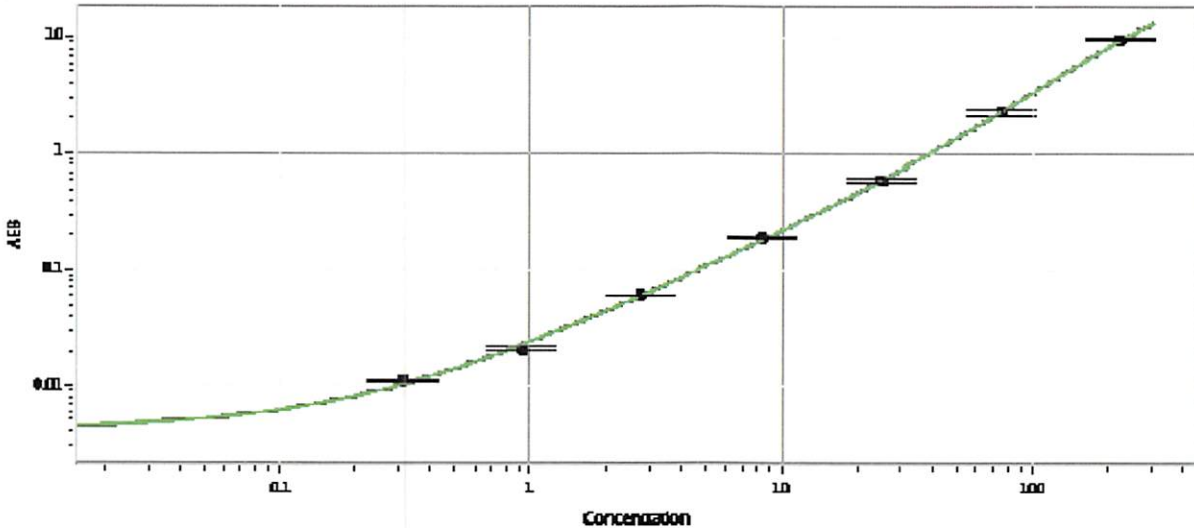
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102374	903301
Detector Stock	102375	834404
Calibrator Concentrate	102379	828408
Calibrator Diluent	102376	834408
Sample Diluent	102378	834407

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

<b>Calibrator Concentrate Target Value</b>	4500 pg/mL
--	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.309
C	0.926
D	2.78
E	8.33
F	25.0
G	75.0
H	225

Review/Approval

Patrizia Stadler  
Name

Assoc. Director, QC  
Title

*P. Stadler 14 Feb 2019*  
Signature/Date