

Product Number: 103071
Lot Number: 502353
Expiration: 11-Dec-2020
Platform(s): HD-1/ HD-X & SR-X



Data below represents results generated on the

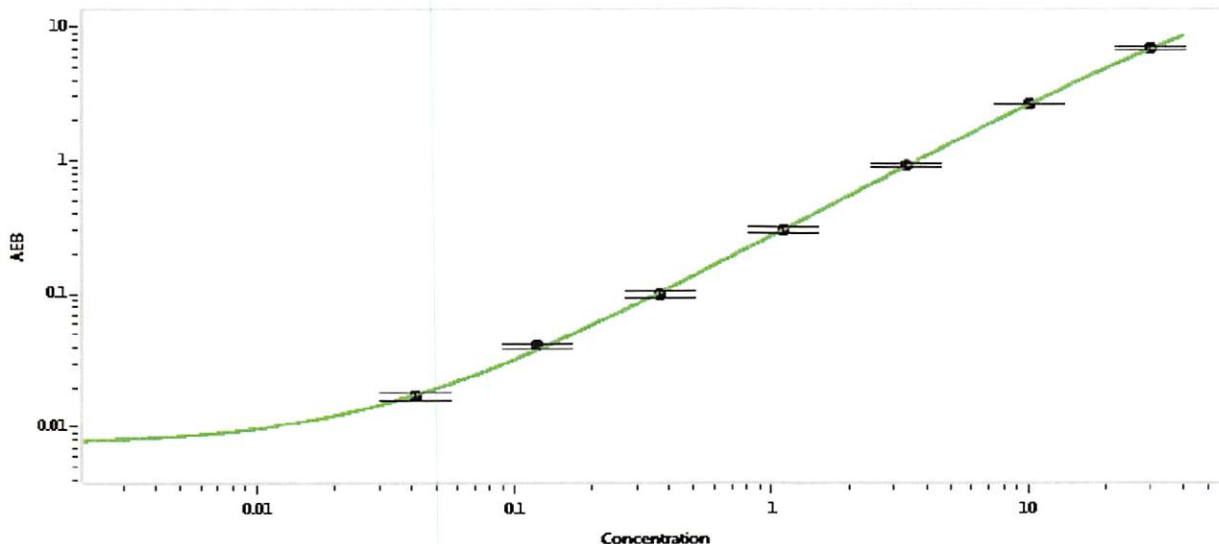
Component	Part Number	Lot Number
Bead Stock	103064	011704
Detector Stock	103065	011701
Calibrator Concentrate	103068	934309
Calibrator Diluent	103066	009210
Sample Diluent	103067	005227

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

600 pg/mL

Calibration Curve



Calibrator Levels
(pg/mL)

A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 04 May 2020
Signature/Date