

Product Number: 103071
Lot Number: 501853
Expiration: 1-Apr-2020
Platform(s): HD-1 & SR-X



Data below represents results generated on the

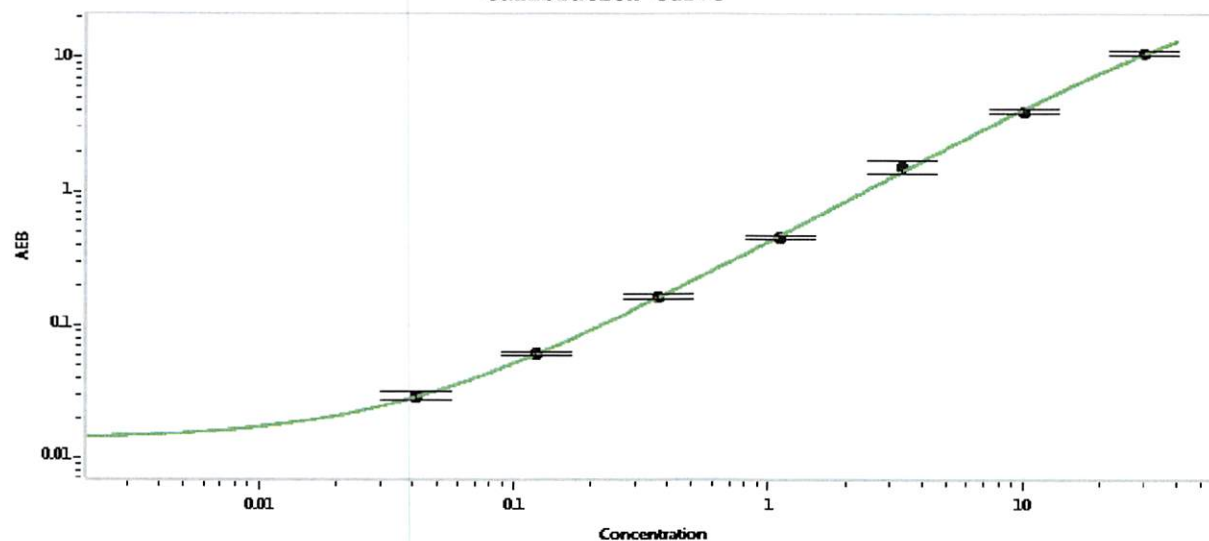
Component	Part Number	Lot Number
Bead Stock	103064	913502
Detector Stock	103065	913309
Calibrator Concentrate	103068	915704
Calibrator Diluent	103066	912017
Sample Diluent	103067	908911

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

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 21 JUN 2019
Signature/Date