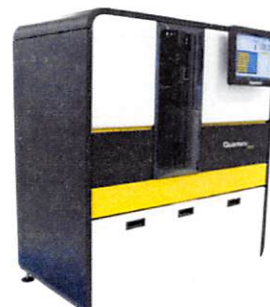


Product Number:	102475
Lot Number:	502451
Expiration:	22-Jan-2021
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	102468	002210
Detector Stock	102469	016908
Calibrator Concentrate	102472	004511
Calibrator Diluent	102470	015115
Sample Diluent	102471	015418

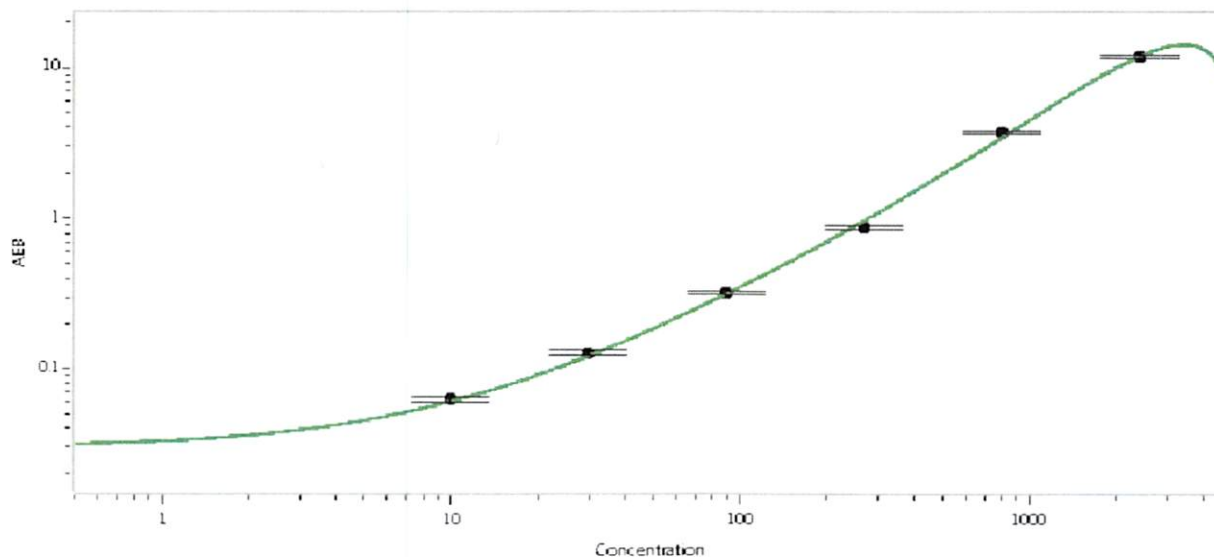
Generic Components - Included With Kit But Not Lot Number Specific		
Discovery Bead Diluent		101954
Discovery Detector Diluent		101955
Discovery SBG Diluent		101956
Discovery SBG Stock		103297
RGP Reagent		103159



Data below represents results generated on the Simoa™ HD-X Analyzer.

Calibrator Concentrate Target Value	60 ng/mL
--	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	9.88
C	29.6
D	88.9
E	267
F	800
G	2400

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler 30 Jun 2020

Name

Title

Signature/Date