

Product Number:	102648
Lot Number:	502373
Expiration:	21-Feb-2021
Platform(s):	HD-1/HD-X & SR-X



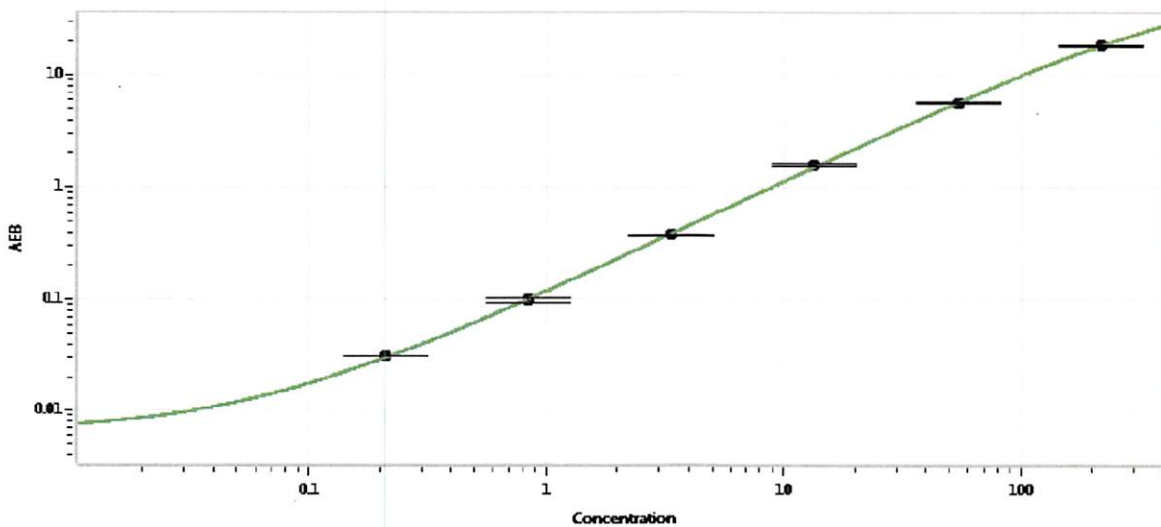
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102641	013102
Detector Stock	102642	013105
Calibrator Concentrate	102645	924602
Calibrator Diluent	102643	010813
Sample Diluent	102644	005231
Detector Diluent	102734	013103

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	3,000 pg/mL
-------------------------------------	-------------

Calibration Curve



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

Shivani Goel
Name

QC Manager
Title

Shivani Goel 15 May 2020
Signature/Date