

Product Number:	102209
Lot Number:	501659
Expiration:	11-Feb-2020
Platform(s):	HD-1 and SR-X



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

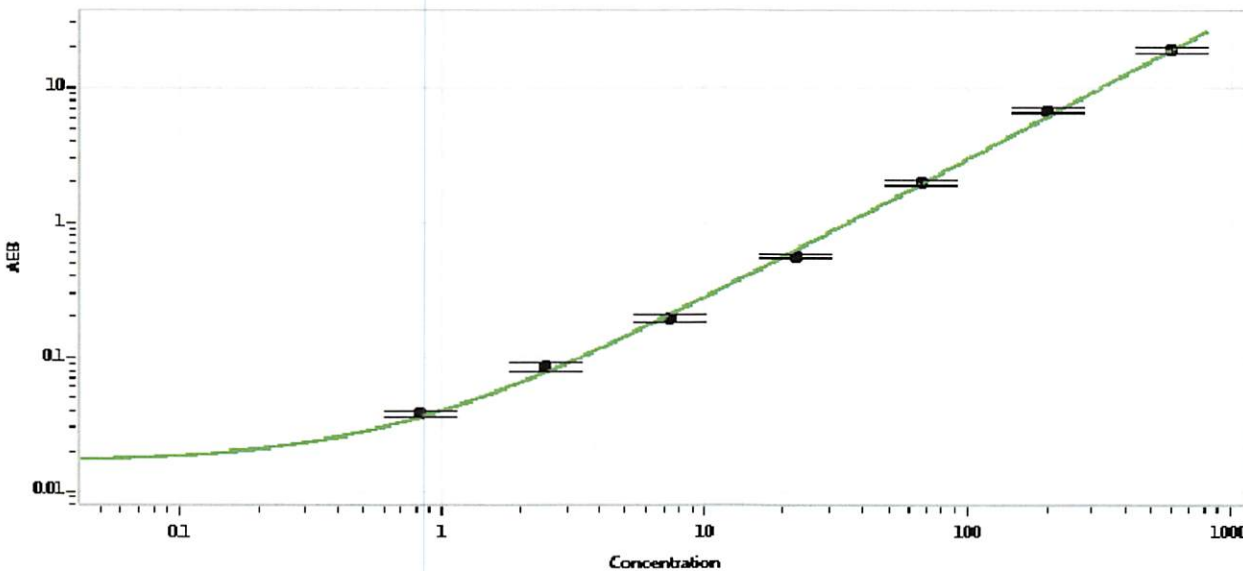
Component	Part Number	Lot Number
Bead Stock	102201	904307
Detector Stock	102202	904208
Calibrator Concentrate	102205	827108
Calibrator Diluent	102203	904210
Sample Diluent	102204	904211

NOTE: For use on the HD-1 platform - refer to the Package Insert (PI-0077).
For use on the SR-X platform - refer to the SR-X Kit Instructions (KI-0083).

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	12 ng/mL
--	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.823
C	2.47
D	7.41
E	22.2
F	66.7
G	200
H	600

Review/Approval

Patrizia stadler
Name

QC Assoc Director
Title

P. Stadler 28 Feb 2019
Signature/Date