

Product Number:	103400
Lot Number:	502059
Expiration:	15-Apr-2020
Platform(s):	HD-1/HD-X and SR-X

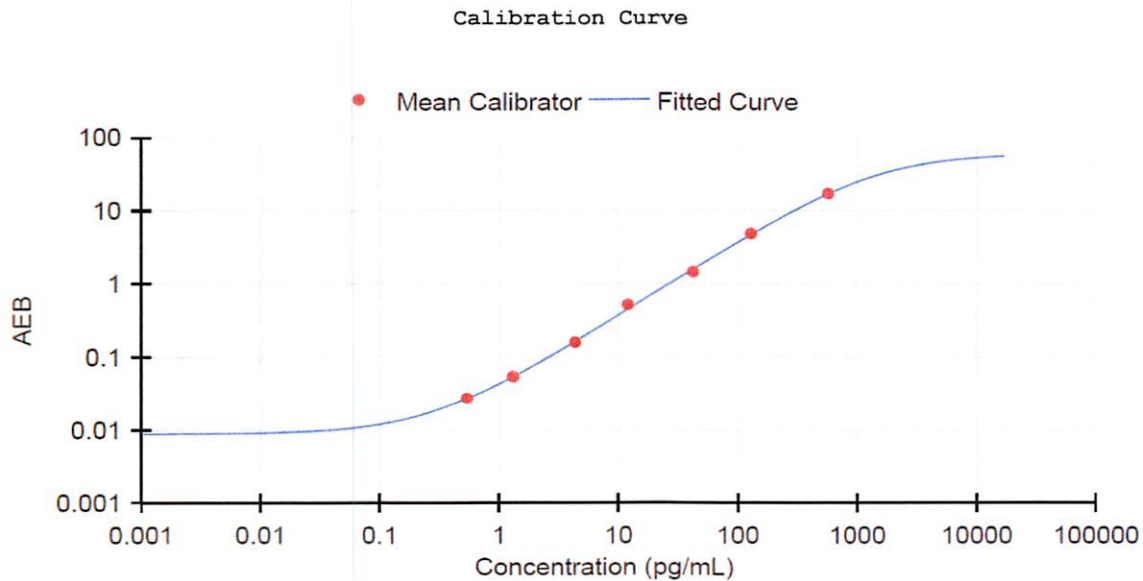


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

Component	Part Number	Lot Number
Bead Reagent	102246	927001
Detector Reagent	102248	926801
SBG Reagent	102250	926608
Sample Diluent	103399	910505
Calibrators	102255	925601
Controls	102256/102257	925901
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific.

Release Materials				Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	925901	2.95	2.58-4.81	
Control 2	925901	114	99-173	
Panel 1	922601	18.6	14.1-21.1	
Panel 2	922602	6.12	4.94-7.41	
Panel 3	922603	39.0	28.4-42.6	
Panel 4	922604	155	129-204	



Calibrator Levels (pg/mL)
A 0.000
B 0.542
C 1.31
D 4.34
E 12.0
F 41.8
G 129
H 567

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

p. Stadler 11 Oct 2019
Signature/Date