

Product Number:	103400
Lot Number:	502045
Expiration:	25-Feb-2020
Platform(s):	HD-1/HD-X and SR-X



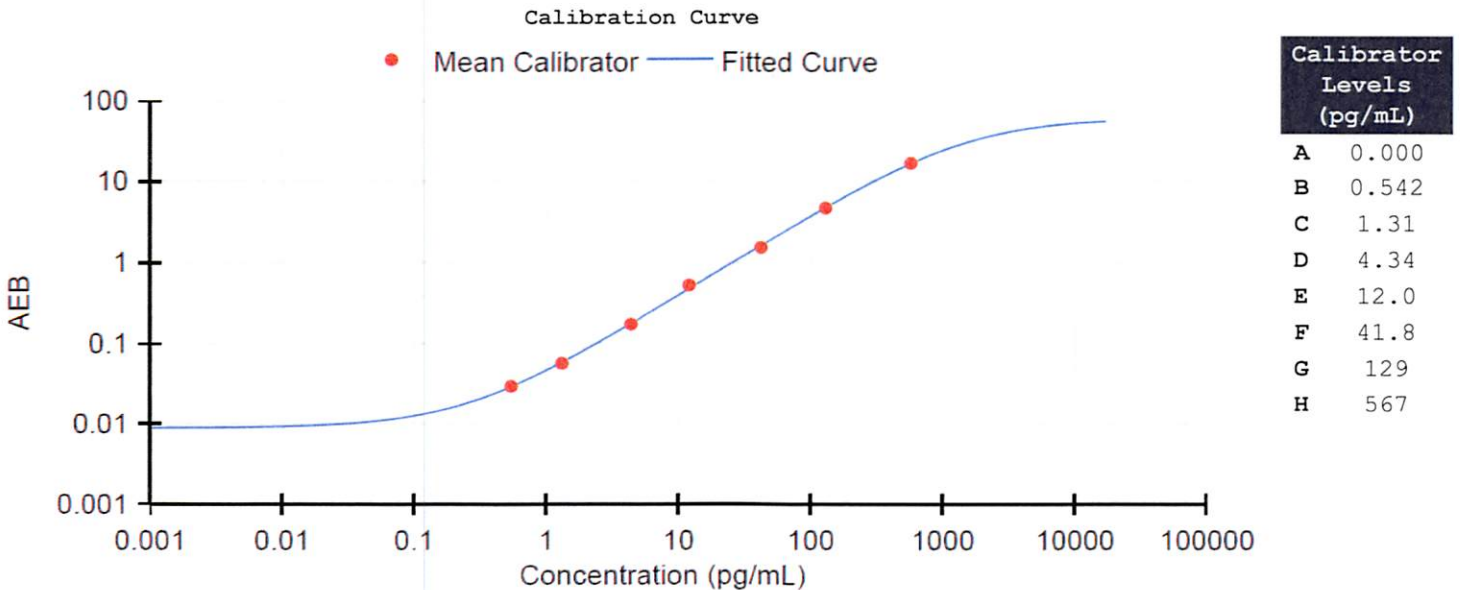
Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

Component	Part Number	Lot Number
Bead Reagent	102246	905611
Detector Reagent	102248	913005
SBG Reagent	102250	921313
Sample Diluent	103399	905601
Calibrators	102255	925601
Controls	102256/102257	925901
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific.

Release Materials			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	925901	4.13	3.07-4.61
Control 2	925901	167	111-170
Panel 1	922601	18.2	14.0-21.0
Panel 2	922602	6.23	4.95-7.42
Panel 3	922603	37.0	28.1-42.1
Panel 4	922604	151	131-205

Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.



Review/Approval	Patrizia Stadler Name	Assoc. Director, QC Title	<i>p. Stadler</i> Signature	27 Sep 2019 Date
-----------------	--------------------------	------------------------------	--------------------------------	---------------------