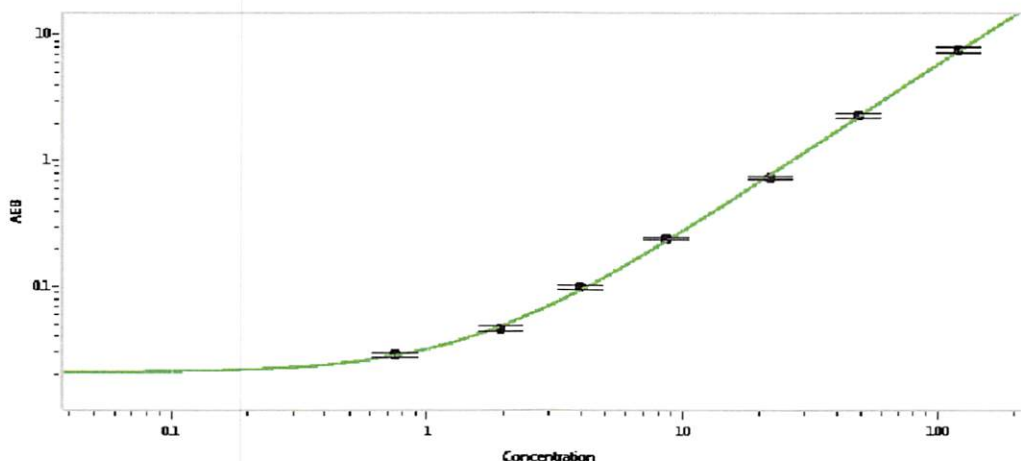


<b>Product Number:</b>	103643
<b>Lot Number:</b>	005008
<b>Expiration:</b>	19-Aug-2020
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.47 x 10 <sup>9</sup> Beads/mL

**Functional Testing**



	Lot	Result (pg/mL)	Mean Range (pg/mL)	* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
Control 1	923503	10.7	6.72-14.2	
Control 2	923503	171	119-217	
Panel 1	003802	61.6	44.3-107	
Panel 2	003802	31.8	25.5-52.5	
Panel 3	003802	71.4	45.2-132	

Review/Approve: Patrizia Stadler Assoc. Director, QC *P. Stadler 30X082020*  
Name Title Signature/Date