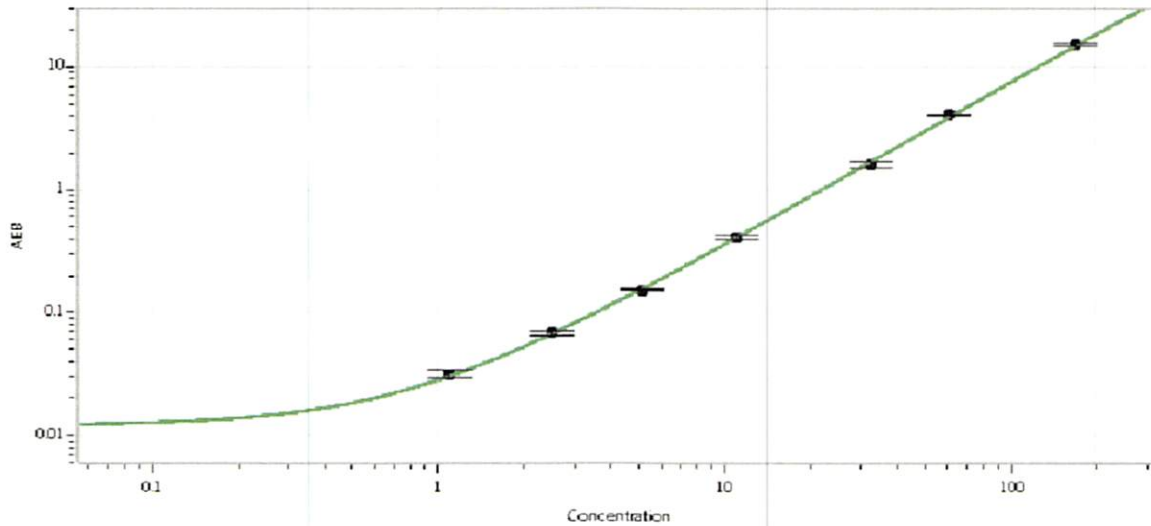


<b>Product Number:</b>	103643
<b>Lot Number:</b>	020308
<b>Expiration:</b>	21-Oct-2020
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.27 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	013902	17.0	10.4-24.3	
Control 2	013902	273	161-376	
Panel 1	003802	52.8	44.3-107	
Panel 2	003802	27.7	25.5-52.5	
Panel 3	003802	65.9	45.2-132	

Review/Approva	Shivani Goel	QC Manager	Shivani Goel 12 Aug 2020
Name	Title	Signature/Date	