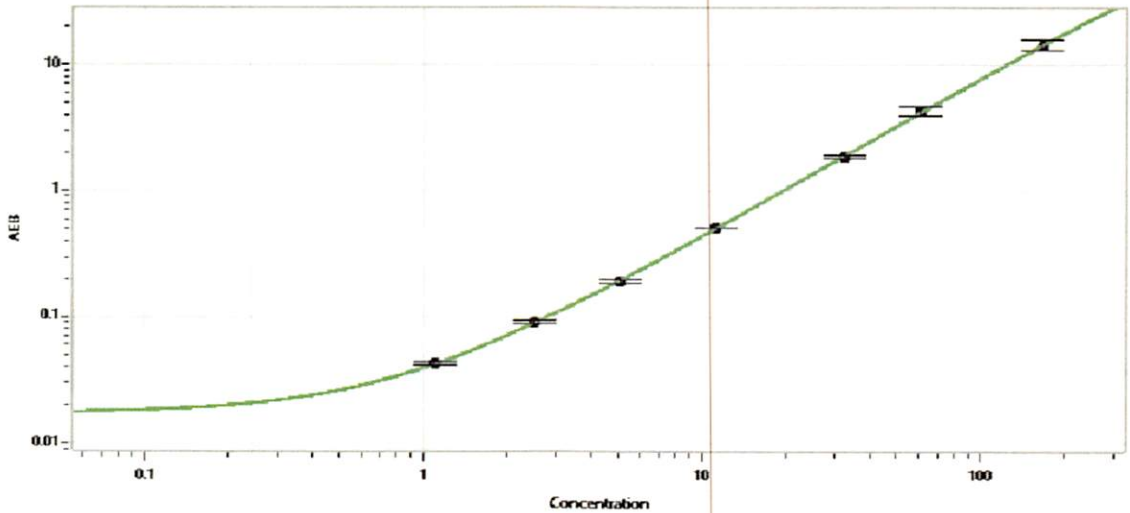


Product Number:	103643
Lot Number:	028717
Expiration:	13-Jan-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.11 x 10 ⁹ 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	20.7	10.4-24.3	
Control 2	013902	280	161-376	
Panel 1	003802	72.9	44.5-103	
Panel 2	003802	35.4	25.1-51.1	
Panel 3	003802	88.3	46.9-127	

Review/Approve

Shivani Goel

QC Manager

Shivani Goel 27 Oct 2020

Name

Title

Signature/Date