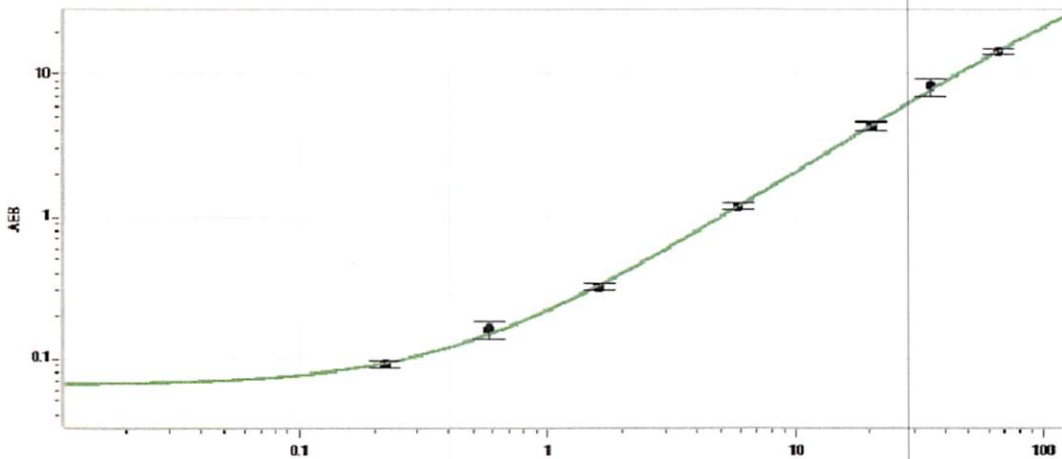


Product Number:	103644
Lot Number:	026813
Expiration:	24-Dec-2020
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Bead Concentration	Multisizer	1.35×10^9 Beads/mL
Functional	HD-X Analyzer	PASS



Review/Approval

Patrizia Stadler Assoc. Director, QC
Name

P. Stadler 01 OCT 2020
Signature/Date