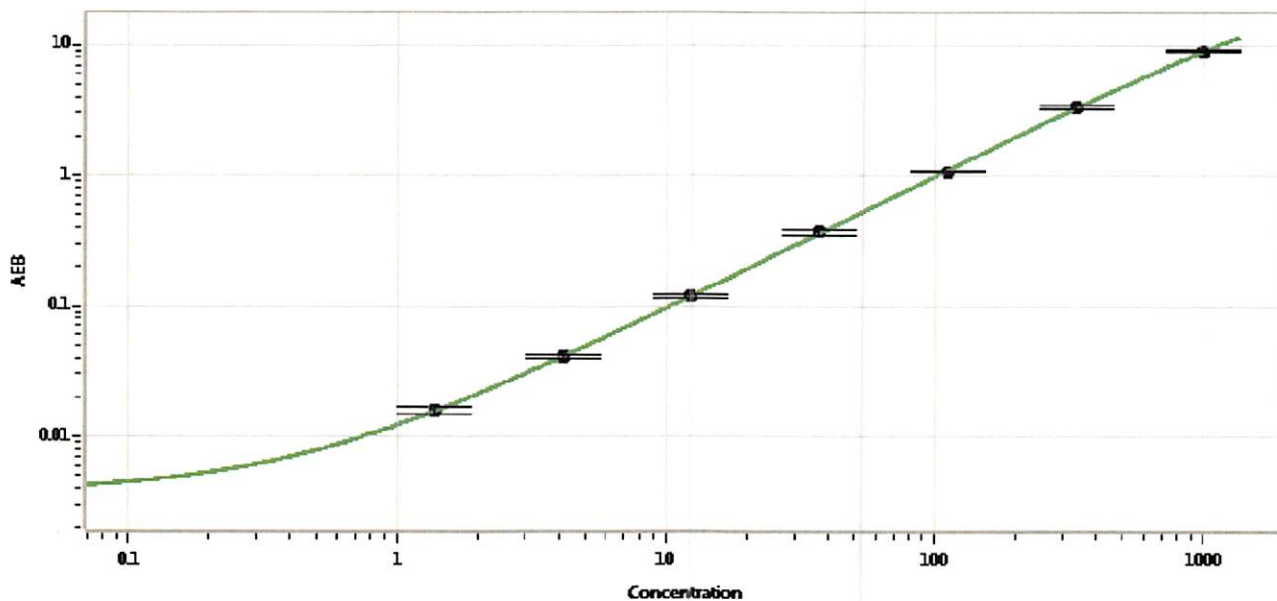


Product Number:	102331
Lot Number:	029306
Expiration:	19-Oct-2021
Storage:	2-8°C

Laboratory Analysis

Means of Analysis	Characteristic	Result
HD-1 Analyzer	Functional Test	PASS

Calibration Curve



Review/Approval

Shivani Goel	QC Manager	<i>Shivani Goel</i> 08 Feb 2021
Name	Title	Signature/Date