

Product Number:	103710
Lot Number:	212909
Expiration:	9-May-2024
Storage:	-80 °C

Laboratory Analysis

Means of Analysis	Characteristic	(ug/mL)
HD-X Analyzer	Concentration	1.0

Review/Approval

Sonia Khillan QC Manager  18 APR 2023

Name	Title	Signature/Date
Sonia Khillan	QC Manager	 18 APR 2023