

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

Name

QC manager

Title

See 1261/c 26JUL2023

Signature/Date