

Product Number:	100877
Lot Number:	219213
Expiration:	11-Jul-2024
Storage:	-80°C

Laboratory Analysis

Means of Analysis	Characteristic	Result (ug/mL)
HD-X Analyzer	Concentration	43

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 20 JUN 2023
Signature/Date