

Product Number:	100463
Lot Number:	303212
Expiration:	26-Oct-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
Concentration	5.1 mg/mL

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia Khillan 08FEB2023

Signature/Date