

Product Number:	100463
Lot Number:	235308
Expiration:	26-Oct-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
Concentration	5.1 mg/mL

Review/Approval

Sania Khillan QC Manager Sania Khillan / 19 JAN 2023

Name

Title

Signature/Date