

Product Number:	100463
Lot Number:	119303
Expiration:	26-Oct-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
Concentration	5.1 mg/mL

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 15 Jul 2021

Signature/Date