

Product Number:	100464
Lot Number:	204107
Expiration:	28-Jan-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
Concentration	1.0 mg/mL

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 15 Feb 2022

Signature/Date