

Product Number:	100464
Lot Number:	303213
Expiration:	22-Aug-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
Concentration	1.1 mg/mL

Review/Approval

Sonia Khillan

QC Manager

Sonia Khillan 10 FEB 2023

Name

Title

Signature/Date