

Product Number:	100464
Lot Number:	235301
Expiration:	28-Jan-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
Concentration	1.0 mg/mL

Review/Approval

Sonia Khillan QC Manager *Sonia Khillan* 1/03/JAN/2023

Name Title Signature/Date