

Product Number:	100463
Lot Number:	220604
Expiration:	26-Oct-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
Concentration	5.1 mg/mL

Review/Approval

Matt Smart

QC

MAT 15 Aug 2022

Name

Title

Signature/Date