

Product Number:	100464
Lot Number:	225612
Expiration:	28-Jan-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
Concentration	1.0 mg/mL

Review/Approval

Matt Smart

Name

QC

Title

M. RA 19 Sep 2022

Signature/Date