

Product Number:	100463
Lot Number:	103512
Expiration:	28-Jan-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
Concentration	5.1 mg/mL

Review/Approval

Junpil Park

Supervisor, QC

 18 Feb 2024

Name

Title

Signature/Date