

Product Number:	100463
Lot Number:	103516
Expiration:	18-Jul-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
Concentration	5.0 mg/mL

Review/Approval

Junpil Park

Supervisor, QC

 05 May 2021

Name

Title

Signature/Date