

Product Number:	100463
Lot Number:	105804
Expiration:	26-Oct-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
Concentration	5.1 mg/mL

Review/Approval

Junpil Park

Name

QC Supervisor

Title

 04Mar2021

Signature/Date