

Product Number:	100463
Lot Number:	122803
Expiration:	26-Oct-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
Concentration	5.1 mg/mL

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 18 Aug 2021
Signature/Date