

Product Number:	100464
Lot Number:	320612
Expiration:	22-Aug-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
Concentration	1.1 mg/mL


Review/Approval

ROBERT SOMOSYI

Name

QC, MAJAGR

Title

 08AUG2023

Signature/Date