

Product Number:	101353
Lot Number:	322620
Expiration:	14-Aug-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.40	7.35 - 7.45	PASS

Review/Approval

Sonia Khilkin

Name

QC Manager

Title

Sonia 12/11/2023

Signature/Date