

Product Number:	101472
Lot Number:	230509
Expiration:	1-Nov-2023
Storage:	2-8 °C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia Khillan / 07NOV2022

Signature/Date