

Product Number:	100457
Lot Number:	210318
Expiration:	13-Apr-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 / 07 JUL 2022
Signature/Date