

Product Number:	100457
Lot Number:	220606
Expiration:	25-Jul-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Matt Smart
Name

QC
Title

MSA 15 Aug 2022
Signature/Date