

Product Number:	103896
Lot Number:	207207
Expiration:	13-Mar-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	pH Target	Result	Disposition
pH	7.35 - 7.45	7.42	Pass

Review/Approval

Matt Smart
Name

QC
Title

MSA 29542022
Signature/Date