

Product Number:	103896
Lot Number:	220914
Expiration:	28-Jul-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	pH Target	Result	Disposition
pH	7.35 - 7.45	7.41	Pass

Review/Approval

Sonia Khillar

Name

QC

Title

Sonia Khillar / 24 AUG 2022

Signature/Date