

Product Number:	103359
Lot Number:	329403
Expiration:	21-Oct-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	8.01	7.95 - 8.05	PASS

Review/Approval

Sonia Khillan
Name

Senior QC Analyst
Title

Sonia Khillan 25 Oct 2023
Signature/Date