

Product Number:	100453
Lot Number:	328909
Expiration:	16-Oct-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	6.18	6.18 - 6.22	PASS

Review/Approval

Sonia Khillan
Name

Senior QC Analyst
Title

Sonia Khillan 13 NOV 2023
Signature/Date