

Product Number:	101955
Lot Number:	331416
Expiration:	10-Nov-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.39	7.30 - 7.50	PASS

Review/Approval

Sonia Khilkin

Name

Senior QC Analyst

Title

Sonia Khilkin 30 NOV 2023

Signature/Date