

Product Number:	101473
Lot Number:	319914
Control Number:	319914A
Expiration:	18-Jul-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.39	7.35 - 7.45	PASS

Review/Approval

ROBERT SOMOGYI

QC, MANAGER



25AUG2023

Name

Title

Signature/Date