

Product Number:	101362
Lot Number:	132509
Expiration:	21-Nov-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 29 Nov 2021
Signature/Date