

Product Number:	103896
Lot Number:	125609
Expiration:	13-Mar-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	pH Target	Result	Disposition
pH	7.35-7.45	7.44	Pass

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 06 Oct 2021

Signature/Date