

Product Number:	103896
Lot Number:	114420
Expiration:	24-Nov-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	pH Target	Result	Disposition
pH	7.35-7.45	7.45	Pass

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 14 Jul 2021

Signature/Date