

Product Number:	103896
Lot Number:	114419
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 Supervisor
Title

Shivani Goel 12 Jul 2021
Signature/Date