

Product Number:	103896
Lot Number:	202710
Expiration:	27-Jan-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	pH Target	Result	Disposition
pH	7.35-7.45	PASS	Pass

Review/Approval

Shivani Goel
Name

QC

Title

Shivani Goel 24 Feb 2022
Signature/Date