

Product Number:	103896
Lot Number:	102528
Expiration:	25-Jan-2022
Storage:	2 - 8°C


Laboratory Analysis

Characteristic	pH Target	Result	Disposition
pH	7.35-7.45	7.44	Pass

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 01Mar2021
Signature/Date