

Product Number:	101956
Lot Number:	205008
Expiration:	19-Feb-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval	Shivani Goel	QC Manager	<i>Shivani Goel</i> 06 Apr 2022
	Name	Title	Signature/Date