

Product Number:	102907
Lot Number:	503689
Expiration:	19-Apr-2023
Storage:	2 - 8°C

Component	Part Number	Lot Number	Means of Analysis	Pass / Fail
Lysate Diluent A Reagent	102908	223608	pH meter	Pass
Lysate Diluent B Reagent	102910	223612		Pass
Lysate Diluent C Reagent	102912	210924		Pass
Lysate Diluent D Reagent	102914	223615		Pass

Review/Approval

Matt Smart

QC

MAA 31 Aug 2022

Name

Title

Signature/Date