

<b>Product Number:</b>	102907
<b>Lot Number:</b>	503688
<b>Expiration:</b>	17-Mar-2023
<b>Storage:</b>	2 - 8°C

Component	Part Number	Lot Number	Means of Analysis	Pass / Fail
Lysate Diluent A Reagent	102908	223408	pH meter	Pass
Lysate Diluent B Reagent	102910	210314		Pass
Lysate Diluent C Reagent	102912	207611		Pass
Lysate Diluent D Reagent	102914	215603		Pass

Review/Approval

*Matt Smart*

*QC*

*MAT 29 Aug 2022*

Name

Title

Signature/Date