

<b>Product Number:</b>	101736
<b>Lot Number:</b>	504296
<b>Expiration Date:</b>	10-Nov-2025



Component	Part Number	Lot Number
RGP Reagent	103159	41310

Means of Analysis	PASS / FAIL
HD-X	PASS

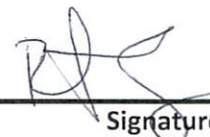
Review/Approval

Robert Somogyi

QC Manager

Name

Title

 24 JUN 2024  
Signature/Date