

Product Number:	101736
Lot Number:	504273
Expiration Date:	17-Oct-2025



Component	Part Number	Lot Number
RGP Reagent	103159	41080

Means of Analysis	PASS / FAIL
HD-X	PASS

Review/Approval

Robert Somogyi	QC Manager		24 MAY 2024
Name	Title	Signature/Date	