

Product Number:	101736
Lot Number:	504650
Expiration Date:	15-Nov-2026

Component	Part Number	Lot Number
RGP Reagent	103159	51350

Means of Analysis	PASS / FAIL
HD-X	PASS



Review/Approval

Phurmi Patel

Name

DOC Control II, QA

Title

SDP / *03 Jul 2025*

Signature/Date