

Product Number:	101736
Lot Number:	502453
Expiration:	13-May-2021
Storage:	2-8°C

Component	Part Number	Lot Number
RGP Reagent	103159	01340

Means of Analysis	PASS/FAIL
HD-1/HD-X	Pass

Review/Approval

Patrizia Stadler Assoc. Director, QC
Name Title

P. Stadler 30 Jun 2020
Signature/Date