

Product Number:	101736 (103159)
Lot Number:	502300 (00570)
Expiration:	26-Feb-2021
Storage:	2-8°C

	PASS/FAIL
Release Criteria	Pass

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P-Stadler

25 MAR 2020

Name

Title

Signature/Date