



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Diluent 101

Catalog Number: R51AD-Series

Lot Number: M1010046

Expiration Date: 31 DEC 2024

Storage Temperature: $\leq -10^{\circ}\text{C}$

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Xiomara Talavera	Quality	<i>X Talavera</i>	26 JUL 2022