



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Diluent 39

Catalog Number: R5ABB-Series

Lot Number: M0390014

Expiration Date: 31 DEC 2022

Storage Temperature: 2-8°C

This product was manufactured according to MSD documents. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Xiomara Talavera	Quality	<i>Talavera</i>	27 APR 2021