



CERTIFICATE OF CONFORMANCE

Product Description: Diluent 41
Catalog Number: R50AH-Series
Lot Number: M0410045
Expiration Date: 31 JUL 2024
Storage Temperature: ≤ -10°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>Talavera</i>	23 SEP 2022