



CERTIFICATE OF CONFORMANCE

Product Description: Diluent 100
Catalog Number: R50AA - Series
Lot Number: Y0270316
Expiration Date: 31 JAN 2025
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	Karen Hamilla	Quality	<i>Karen Hamilla</i>	07 SEP 2022