



CERTIFICATE OF CONFORMANCE

Product Description: Diluent 100
Catalog Number: R50AA-Series
Lot Number: Y0270327
Expiration Date: 30 SEP 2025
Storage Temperature: 2-8°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	Karen Hamilla	Quality	<i>Karen Hamilla</i>	18 APR 2023