



CERTIFICATE OF CONFORMANCE

Product Description: Diluent 100
Catalog Number: R50AA-Series
Lot Number: Y0270322
Expiration Date: 30 JUN 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>	01 FEB 2023