



**CERTIFICATE OF CONFORMANCE**

**Product Description:** Diluent 100  
**Catalog Number:** R50AA-Series  
**Lot Number:** Y0270332  
**Expiration Date:** 31 MAR 2026  
**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.

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Review/Approval</b>	Karen Hamilla	Quality	<i>Karen Hamilla</i>	25 OCT 2023