



**CERTIFICATE OF CONFORMANCE**

**Product Description:** Diluent 49  
**Catalog Number:** R50AM-Series  
**Lot Number:** M0490017  
**Expiration Date:** 28 FEB 2026  
**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

		<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Review/Approval</b>	Karen Hamilla	Quality	<i>Karen Hamilla</i>	03 OCT 2023