




### CERTIFICATE OF CONFORMANCE

**Product Description:** Diluent 43  
**Catalog Number:** R50AG-Series  
**Lot Number:** M0430052  
**Expiration Date** 31 DEC 2025  
**Storage Temperature:**  $\leq -10^{\circ}\text{C}$

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>	Xiomara Talavera	Quality		03 JAN 2024