



### CERTIFICATE OF CONFORMANCE

**Product Description:** Diluent 57  
**Catalog Number:** R50BZ -Series  
**Lot Number:** M0570011  
**Expiration Date:** 31 JUL 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>	Karen Hamilla	Quality	<i>Karen Hamilla</i>	14 MAR 2023