




### CERTIFICATE OF CONFORMANCE

**Product Description:** MSD Read Buffer T (4X) Surfactant Free  
**Catalog Number:** R92TD-Series  
**Lot Number:** Y0060048  
**Expiration Date:** 30 APR 2027  
**Storage Temperature:** Room Temperature

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		16 MAY 2024