


**CERTIFICATE OF CONFORMANCE**

**Product Description:** MSD Read Buffer T (4X) With Surfactant  
**Catalog Number:** R92TC-Series  
**Lot Number:** Y0140426  
**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.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		04 JUN 2024