



CERTIFICATE OF CONFORMANCE

Product Description: MSD Read Buffer T (4X) With Surfactant
Catalog Number: R92TC-Series
Lot Number: Y0140421
Expiration Date 31 JAN 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	<i>Karen Hamilla</i>	14 Feb 2024