



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD Read Buffer T (4x), Surfactant Free  
**Catalog Number:** R92TD-1-Series  
**Lot Number:** Y0060047  
**Expiration Date:** 31 JUL 2026

**Statement:** MSD Read Buffer T (4x), Surfactant Free is verified to ensure that consistent signals are obtained across lots. Details on the quality control testing of MSD Read Buffer T (4x), Surfactant Free are provided in the product insert available at [www.mesoscale.com](http://www.mesoscale.com).

MSD Read Buffer T (4x), Surfactant Free was tested using SULFO-TAG™ Conjugated Biotinylated IgG (BTI) lot # R0010830 and MSD GOLD™ 96-well Streptavidin SECTOR Plate lot # Z0021899

Concentration of BTI	Mean ECL Spec and Tolerance Range	Intraplate CV or SD spec	Pass/Fail
BTI with 4X	27429 - 41131 counts	15 %	PASS
Dil-100 with 4X	0 - 240 counts	15 counts	PASS

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality	<i>Karen Hamilla</i>	13 SEP 2023