



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD Read Buffer T (4x), Surfactant Free

**Catalog Number:** R92TD-Series

**Lot Number:** Y0060045

**Expiration Date:** 31 JUL 2025

**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 # R0010769 and MSD GOLD™ 96-well High Bind Avidin SECTOR Plate Z0010674. Specifications are BTI lot specific.

Concentration of BTI	Mean ECL Spec and Tolerance Range	Intraplate CV or SD spec	Pass/Fail
BTI with 4X	48609 - 72914 counts	15 %	PASS
Dil-100 with 4X	0 - 158 counts	50 %	PASS

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		26 AUG 2022