



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD Read Buffer T (1x) with Surfactant

**Catalog Number:** R92TA-Series

**Lot Number:** Y0110101

**Expiration Date:** 31 DEC 2026

**Statement:** MSD Read Buffer T (1x) with Surfactant is verified to ensure that consistent signals are obtained across lots. Details on the quality control testing of MSD Read Buffer T (1x) with Surfactant are provided in the product insert available at [www.mesoscale.com](http://www.mesoscale.com).

MSD Read Buffer T (1x) with Surfactant was tested using SULFO-TAG™ Conjugated Biotinylated IgG (BTI) lot # R0010830 and MSD GOLD™ 96-well Streptavidin SECTOR Plate Z0022134. Specifications are BTI lot specific.

Concentration of BTI	Mean ECL Spec and Tolerance Range	Intraplate CV or SD spec	Pass/Fail
BTI with 1X	20733 – 38503 counts	15 %	Pass
Dil-100 with 1X	19– 44 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		09 Feb 2024