



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** R92TC-Series

**Lot Number:** Y0140412

**Expiration Date:** 31 JAN 2026

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

MSD Read Buffer T (4x) with Surfactant was tested using SULFO-TAG™ Conjugated Biotinylated IgG (BTI) lot # R0010769 and MSD GOLD™ 96-well Streptavidin SECTOR Plate Z0021899. Specifications are BTI lot specific.

Concentration of BTI	Mean ECL Spec and Tolerance Range	Intraplate CV or SD spec	Pass/Fail
BTI with 4X	31137 – 46705 counts	10 %	Pass
Dil-100 with 4X	66 – 156 counts	15 counts	Pass
BTI with 2X	34665 – 51998 counts	10%	PASS
Dil-100 with 2X	38 – 88 counts	15 counts	PASS
BTI with 1X	19533 – 36278 counts	10%	PASS
Dil-100 with 1X	19 – 44 counts	15 counts	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 MAR 2023