



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD® Read Buffer T(1X) with Surfactant  
**Catalog Number:** R92TA-Series  
**Lot Number:** Y0110096  
**Expiration Date:** 31 DEC 2025

**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 #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 1X	25565 – 47479 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
<b>Review/Approval</b>	Xiomara Talavera	Quality		08 FEB 2023