



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD<sup>®</sup> Read Buffer T(1X) with Surfactant  
**CatalogNumber/Size:** R92TA-Series  
**Lot Number:** Y0110084  
**Expiration Date:** 31 JAN 2024  
**Storage Temperature:** Room Temperature

**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<sup>™</sup> Conjugated Biotinylated IgG (BTI) lot # R0010698 and MSD GOLD<sup>™</sup> 96-well Streptavidin SECTOR Plate lot # Z0021486.

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

**Statement:** This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		24 FEB 2021