



**CERTIFICATE OF ANALYSIS**

**Product Description:** MSD® Read Buffer T(1X) with Surfactant  
**CatalogNumber/Size:** R92TA-Series  
**Lot Number:** Y0110076  
**Expiration Date:** 31 Jul 2022  
**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™ Conjugated Biotinylated IgG (BTI) lot # R0010605 and MSD GOLD™ 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		15 AUG 2019