



**CERTIFICATE OF ANALYSIS**

**Product Description:** MSD Read Buffer T (4x) Surfactant Free  
**CatalogNumber/Size:** R92TD-Series  
**Lot Number:** Y0060033  
**Expiration Date:** 31 DEC 2022  
**Storage Temperature:** Room Temperature

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

MSD Read Buffer T (4x) Surfactant Free was tested using SULFO-TAG™ Conjugated Biotinylated IgG (BTI) lot # R0010605 and MSD GOLD™ 96-well High Bind Avidin SECTOR Plate lot # Z0010663

Concentration of BTI	Mean ECL spec and tolerance range	Intraplate CV or SD spec	Pass/Fail?
BTI with 4X	48609 - 72914 counts	15 %	PASS
Dil-100 with 4X	0 - 158 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		11 DEC 2019