



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Read Buffer T (1x) with Surfactant

Catalog Number: R92TA-Series

Lot Number: Y0110090

Expiration Date: 30 JUN 2024

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.

MSD Read Buffer T (1x) with Surfactant was tested using SULFO-TAG™ Conjugated Biotinylated IgG (BTI) lot # R0010698 and MSD GOLD™ 96-well Streptavidin SECTOR Plate 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

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		28 JUL 2021