



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD Read Buffer T (4x) with Surfactant

Catalog Number: R92TC-Series

Lot Number: Y0140414

Expiration Date: 31 MAR 2026

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

MSD Read Buffer T (4x) 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 4X	31137 – 46705 counts	10 %	Pass
Dil-100 with 4X	66 – 156 counts	15 counts	Pass
BTI with 2X	34665 – 51998 counts	10%	Pass
Dil-100 with 2X	38 – 88 counts	15 counts	Pass
BTI with 1X	19533 – 36278 counts	10%	Pass
Dil-100 with 1X	19 – 44 counts	15 counts	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		25 APR 2023