



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: R92TC-Series

Lot Number: Y0140400

Expiration Date: 28 FEB 2025

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 Z0021725. 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	22325 - 33488 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		28 MAR 2022