



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: R92TC-2

Lot Number: Y0140403

Expiration Date: 30 APR 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		23 MAY 2022