



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: R92TD-Series

Lot Number: Y0060046

Expiration Date: 31 DEC 2025

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.

MSD® Read Buffer T (4x), Surfactant Free was tested using SULFO-TAG™ Conjugated Biotinylated IgG (BTI) lot # R0010769 and MSD GOLD™ 96-well High Bind Avidin 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	48609 - 72914 counts	15 %	PASS
Dil-100 with 4X	0 - 158 counts	50 %	PASS

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

	Name	Function	Signature	Date
Review/Approval	Nithya Santhanam	Quality		13 Feb 2023