



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: R92TD-Series

Lot Number: Y0060041

Expiration Date: 30 APR 2024

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 Select Product Description 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 # R0010698 and MSD GOLD™ 96-well High Bind Avidin SECTOR Plate Z0010674.

Concentration of BTI	Mean ECL Spec and Tolerance Range	Intraplate CV or SD spec	Pass/Fail
BTI with 4X	48609 - 72914	15%	Pass
Dil-100 with 4X	0 - 158	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		23 JUN 2021