



### CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** R92TC-Series

**Lot Number:** Y0140390

**Expiration Date:** 30 APR 2024

**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 Select Product Description are provided in the product insert available at [www.mesoscale.com](http://www.mesoscale.com).

MSD Read Buffer T (4x) with Surfactant was tested using SULFO-TAG™ Conjugated Biotinylated IgG (BTI) lot # R0010698 and MSD GOLD™ 96-well Streptavidin SECTOR Plate Z0021486.

Concentration of BTI	Mean ECL Spec and Tolerance Range	Intraplate CV or SD spec	Pass/Fail
BTI with 4X	36622 – 54933	15%	Pass
Dil-100 with 4X	66 - 156	50%	Pass
BTI with 2X	43205 - 58454	15%	Pass
Dil-100 with 2X	38 - 88	50%	Pass
BTI with 1X	25565 - 47479	15%	Pass
Dil-100 with 1X	19 - 44	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		28 MAY 2021