



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD[®] Read Buffer T(1X) with Surfactant
Catalog Number: R92TA-Series
Lot Number: Y0110097
Expiration Date: 31 JAN 2026

Statement: MSD[®] Read Buffer T(1X) with Surfactant is verified to ensure that consistent signals are obtained across lots. Details on the quality control testing of MSD[®] Read Buffer T(1X) with Surfactant are provided in the product insert available at www.mesoscale.com.

MSD[®] Read Buffer T(1X) with Surfactant was tested using SULFO-TAG[™] Conjugated Biotinylated IgG (BTI)lot #R0010769 and MSD GOLD[™] 96-well Streptavidin 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 1X	20733 – 38503 counts	15%	PASS
Dil-100 with 1X	19 – 44 counts	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		09 MAR 2023