



CERTIFICATE OF ANALYSIS

Product Description: MSD Read Buffer T (4x) Surfactant Free
CatalogNumber/Size: R92TD-Series
Lot Number: Y0060037
Expiration Date: 31 MAY 2023
Storage Temperature: Room Temperature

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 # R0010605 and MSD GOLD™ 96-well High Bind Avidin SECTOR Plate lot # Z0010663

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

Statement: This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		15 MAY 2020