



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD Read Buffer T(4X) with Surfactant
CatalogNumber/Size: R92TC-Series
Lot Number: Y0140382
Expiration Date: 30 SEP 2023
Storage Temperature: Room Temperature

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 MSD Read Buffer T(4X) with Surfactant are provided in the product insert available at www.mesoscale.com.

T(4X) with Surfactant was tested using SULFO-TAG™ Conjugated Biotinylated IgG (BTI) lot # R0010605 and MSD GOLD™ 96-well Streptavidin SECTOR Plate lot # Z0021486.

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

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

	Name	Function	Signature	Date
Review/Approval	Carissa Hubbard	Quality		16 SEP 2020