



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD Read Buffer T (4X) With Surfactant

**Catalog Number:** R92TC-Series

**Lot Number:** Y0140366

**Expiration Date:** 31 MAR 2022

Signal test, using controlled reagents and plates:


ECL, positive control (tolerance range): 1X: 25 565 – 47 479

2X: 43 205 – 58 454

Baseline ECL (tolerance range): 1X: 19 – 44

2X: 38 – 88

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Review/Approval</b>	Dawn Stollar	Quality		22 APR 2019