



**CERTIFICATE OF ANALYSIS**

Page 1 of 1

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

**Catalog #/Size:** R92TC-Series

**Lot Number:** Y0140364

**Expiration Date:** 31 DEC 2021

**Statement:** This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

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 Assurance | <i>Dawn Stollar</i> | 05 FEB 2019 |