



**CERTIFICATE OF ANALYSIS**

Page 1 of 1

**Product Description:** MSD Read Buffer T (4X) With Surfactant  
**Catalog #/Size:** R92TC-Series  
**Lot Number:** Y0140345  
**Expiration Date:** 31 MAY 2020

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>	08 JUN 2017