



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog #/Size:** R92TD - Series

**Lot Number:** Y0060029

**Expiration Date:** 30 JUN 2021

Signal test, using controlled reagents and plates:

ECL, positive control (tolerance range): 72914 – 48609

Baseline ECL (tolerance range):  $\leq 158$

**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>	13 AUG 2018