



## CERTIFICATE OF ANALYSIS

COA - 07898

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

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

**Lot Number:** Y0140341

**Expiration Date:** 31 JAN 2020

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

pH:  $7.8 \pm 0.1$

Signal test, using controlled reagents and plates:

ECL, positive control (mean  $\pm$  tolerance): 1X:  $36\,522 \pm 30\%$

2X:  $50\,830 \pm 15\%$

Baseline ECL: 1X:  $31.5 \pm 40\%$

2X:  $63 \pm 40\%$

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

	Name	Function	Signature	Date
<b>Preparer</b>	Xiomara Talavera	Quality Control	<i>X Talavera</i>	18 FEB 2017
<b>Review/Approval</b>	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	22 FEB 2017