



CERTIFICATE OF ANALYSIS

COA - 07242

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

Catalog #/Size: R92TD-3 / 50 mL; R92TD-2 / 200 mL; R92TD-1 / 1000 mL

Lot Number: Y0060024

Expiration Date: 31 MAY 2019

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

pH: 7.8 ± 0.1

Signal test, using controlled reagents and plates:

ECL, positive control (mean \pm tolerance): $44\ 332 \pm 20\%$

Baseline ECL: ≤ 158

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

	Name	Function	Signature	Date
Preparer	Xiomara Talavera	Quality Control	<i>Xiomara Talavera</i>	26 AUG 2016
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	29 AUG 2016