



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog #/Size: R92TD - Series

Lot Number: Y0060025

Expiration Date: 30 NOV 2019

Signal test, using controlled reagents and plates:

ECL, positive control (tolerance range): 53198 – 35466

Baseline ECL (tolerance range): ≤ 158

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

	Name	Function	Signature	Date
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	18MAY2017