



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog #/Size: R92TD - Series

Lot Number: Y0060028

Expiration Date: 31 OCT 2020

Signal test, using controlled reagents and plates:

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

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	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	19 DEC 2017