



## CERTIFICATE OF ANALYSIS

COA - 07794

**Product Description:** MSD Phosphate Buffer (5x)

**Catalog #/Size:** R93SA - Series

**Lot Number:** Y0170102

**Expiration Date:** 30 APR 2019

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

pH: 7.0 ± 0.1

**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>Preparer</b>	Xiomara Talavera	Quality Control	<i>X Talavera</i>	19 JAN 2017
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	23 JAN 2017