



## CERTIFICATE OF ANALYSIS

COA - 05907

**Product Description:** MSD Tris Wash Buffer (10x)

**Catalog #/Size:** R61TX-1 / 1000 mL; R61TX-2 / 200 mL



**Lot Number:** Y0500024

**Expiration Date:** 30 JUN 2016

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

pH:  $7.5 \pm 0.1$

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

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		25 SEP 2015
Review/Approval	Monika Jain	Quality Assurance		25 SEP 2015