



CERTIFICATE OF ANALYSIS

COA - 05353

Product Description: MSD Tris Wash Buffer (10x)

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



Lot Number: Y0500023

Expiration Date: 29 FEB 2016

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

pH: 7.5 ± 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		02 APR 2015
Review/Approval	Anastasia Belilovets	Quality Assurance		03 APR 2015