



CERTIFICATE OF ANALYSIS

COA - 11211

Product Description: MSD Tris Wash Buffer (10X)

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



Lot Number: Y0500016

Expiration Date: 30 APR 2014

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		29 APR 2013
Review/Approval	Michael Eager	Quality Control		21 MAY 2013