



## CERTIFICATE OF ANALYSIS

COA - 04511

**Product Description:** MSD Tris Lysis Buffer

**Catalog #/Size:** R60TX-3 / 50 mL; R60TX-2 / 200 mL



**Lot Number:** Y0490300

**Expiration Date:** 31 MAR 2015

**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
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		12 JUN 2014
<b>Review/Approval</b>	Michael Eager	Quality Control		29Jun2014