



CERTIFICATE OF ANALYSIS

COA - 06464

Product Description: MSD Tris Lysis Buffer

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



Lot Number: Y0490304

Expiration Date: 30 NOV 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		08 FEB 2016
Review/Approval	Dawn Stollar	Quality Assurance		08 FEB 2016