



CERTIFICATE OF ANALYSIS

COA - 06006

Product Description: MSD Tris Lysis Buffer

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



Lot Number: Y0490303

Expiration Date: 31 MAY 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		25 SEP 2015
Review/Approval	Monika Jain	Quality Assurance		25 SEP 2015