



CERTIFICATE OF ANALYSIS

COA - 06703

Product Description: MSD Tris Wash Buffer (10x)

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



Lot Number: Y0500025

Expiration Date: 30 SEP 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 |  | 29 MAR 2016 |
| Review/Approval | Dawn Stollar | Quality Assurance |  | 29 MAR 2016 |