



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: MSD Phosphate Buffer (5X)


Catalog Number: R93SA - Series

Lot Number: Y0170119

Expiration Date: 31 AUG 2024

Storage Temperature: Room Temperature

This product was manufactured according to MSD documents. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|------------------------|---------------|-----------------|--|-------------|
| Review/Approval | Karen Hamilla | Quality |  | 20 MAY 2022 |