



## CERTIFICATE OF CONFORMANCE

Page 1 of 1

**Product Description:** MSD Phosphate Buffer (5X)


**Catalog Number:** R93SA-Series

**Lot Number:** Y0170117

**Expiration Date:** 31 DEC 2023

**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.

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Review/Approval</b>	Karen Hamilla	Quality		27 JUL 2021