



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: MSD® Phosphate Buffer (5X)


Catalog Number: R93SA-Series

Lot Number: Y0170114

Expiration Date: 30 JUN 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.

	Name	Function	Signature	Date
Review/Approval	Carissa B. Hubbard	Quality		06 APR 2021