



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: MSD® Blocker A
Catalog Number: R93BA - Series
Lot Number: Y0181068
Expiration Date: 31 JUL 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	Karen Hamilla	Quality	<i>Karen Hamilla</i>	14 JUL 2021