



## CERTIFICATE OF CONFORMANCE

Page 1 of 1

**Product Description:** MSD® Blocker A  
**Catalog Number:** R93BA - Series  
**Lot Number:** Y0181054  
**Expiration Date:** 31 AUG 2022  
**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	<i>Karen Hamilla</i>	14 OCT 2020