



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: MSD Blocker D-R


Catalog Number: R93BR - Series

Lot Number: R0010641

Expiration Date: 28 FEB 2021

Storage Temperature: $\leq -10^{\circ}\text{C}$

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		09 JUL 2019