



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Blocker S1
Catalog Number: R93AG-1
Lot Number: R0010757
Expiration Date: 31 JAN 2024
Storage Temperature: ≤ -10°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	Xiomara Talavera	Quality	<i>Talavera</i>	15 NOV 2021