



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: SULFO-TAG™ Anti-Human IgA Antibody

Catalog Number: D21ADE-Series

Lot Number: D00V0020

Expiration Date: 30 JUN 2025

Storage Temperature: 2-8°C

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. 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>	22 APR 2022