



CERTIFICATE OF CONFORMANCE

Page 1 of 1

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


Catalog Number: D21ADE-3

Lot Number: D00V0012

Expiration Date: 28 FEB 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		22 SEP 2021