



CERTIFICATE OF CONFORMANCE

Product Description: SULFO-TAG™ Anti-Human IgG Antibody
Catalog Number: D21ADF-Series
Lot Number: D00V0027
Expiration Date: 30 NOV 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>	30 JUN 2022