



### CERTIFICATE OF CONFORMANCE

**Product Description:** SULFO-TAG™ Anti-Mouse IgM Antibody  
**Catalog Number:** D22AGR-Series  
**Lot Number:** D00V0031  
**Expiration Date:** 31 JAN 2027  
**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.

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Review/Approval</b>	Karen Hamilla	Quality		21 AUG 2023