



## EXTENSION MEMO


Page 1 of 1

**Product Description(s):** SULFO-TAG™Anti-Human IgM Antibody

**Catalog Number(s) / Size(s):** D21ADD-3

**Lot Number:** D00V0011

Based on additional testing, the original date of expiration for this antibody has been extended by 12 months. The expiration date of D00V0011 has been extended to 31 AUG 2025.

Function	Name	Signature	Date
Quality	Karen Hamilla		25 JUN 2024

This memo supersedes all other statements regarding date of expiration for this product.



**CERTIFICATE OF CONFORMANCE**

Page 1 of 1

**Product Description:** SULFO-TAG™Anti-Human IgM Antibody

**Catalog Number:** D21ADD-3

**Lot Number:** D00V0011

**Expiration Date:** 31 AUG 2024

**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	<i>Karen Hamilla</i>	26 MAR 2021