



## EXTENSION MEMO


Page 1 of 1

**Product Description(s):** SULFO-TAG™ Human ACE2 Protein

**Catalog Number:** D21ADG-3

**Lot Number:** D0081760

Based on additional testing, the original date of expiration for this antibody has been extended by 12 months. The expiration date of D0081760 has been extended to 31 JAN 2025.

Function	Name	Signature	Date
Quality	Karen Hamilla		05 OCT 2023

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



## CERTIFICATE OF CONFORMANCE

Page 1 of 1

**Product Description:** SULFO-TAG™ Human ACE2 Protein


**Catalog Number:** D21ADG-3

**Lot Number:** D0081760

**Expiration Date:** 31 JAN 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		17 FEB 2022