



CERTIFICATE OF CONFORMANCE

Product Description: SULFO-TAG™ Anti-Rat IgG Antibody
Lot Number: D00V0028
Expiration Date: 31 Dec 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	Xiomara Talavera-Herrera	Quality	<i>X Talavera</i>	08 JUL 2022