



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Influenza H5 Bridging Serology Assay Kit


Catalog Number: K150AXJU-Series

Lot Number: K0082716

Expiration Date: 31 OCT 2027

Component Description	Lot Number	Storage Temperature	Expiration Date
BIOTIN-H5 Conjugated Protein	538534	$\leq -70^{\circ}\text{C}$	31 OCT 2027
SULFO-TAG H5 Protein	538533	$\leq -70^{\circ}\text{C}$	31 OCT 2027
H5 Control 1.1	A00C1010	$\leq -70^{\circ}\text{C}$	30 APR 2035
H5 Control 1.2	A00C1011	$\leq -70^{\circ}\text{C}$	30 APR 2035
H5 Control 1.3	A00C1012	$\leq -70^{\circ}\text{C}$	30 APR 2035
MSD GOLD 96-Well Small Spot Streptavidin plate	Not kit lot specific	2-8°C	See lot specific certificate
Diluent 100	Not kit lot specific	2-8°C	See lot specific certificate
Blocker A	Not kit lot specific	Room Temperature	See lot specific certificate
MSD GOLD Read Buffer B	Not kit lot specific	Room Temperature	See lot specific certificate

This product was manufactured according to approved MSD documents. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		23 MAY 2025