



## CERTIFICATE OF CONFORMANCE

Page 1 of 1

**Product Description:** MSD® JAK-STAT Whole Cell Lysate Set


**Catalog Number:** C10KQ-1

**Lot Number:** CL0010123/CL0010124

**Expiration Date:** 30 APR 2026

**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

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	Quality		19 AUG 2021