



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: MSD® STAT Panel Whole Cell Lysate Set


Catalog Number: C1202-1

Lot Number: CL0010111 / CL0010112

Expiration Date: 31 MAR 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	Karen Hamilla	Quality		10 MAY 2021