



## CERTIFICATE OF CONFORMANCE

Page 1 of 1

**Product Description:** MSD® Human Myeloperoxidase (MPO) Kit

**Catalog Number:** K151EEC- Series

**Lot Number:** K0050855

**Expiration Date:** N/A (See expiration date on specific components below)

### Plate Specifications:

| Coating Uniformity             | Specifications |
|--------------------------------|----------------|
| Mean Intraplate CV             | ≤ 15%          |
| Mean Interplate CV             | ≤ 18%          |
| No more than one plate with CV | > 18%          |
| No plate with CV               | > 20%          |

### Components:

| Description                       | Lot Number | Storage Temperature | Expiration Date |
|-----------------------------------|------------|---------------------|-----------------|
| Human Myeloperoxidase (MPO) Plate | Z0056539   | 2-8°C               | 30 NOV 2021     |
| Human MPO Calibrator              | A0020251   | ≤ -70°C             | 30 APR 2021     |
| Anti-hMPO Antibody                | D0050797   | 2-8°C               | 30 NOV 2022     |

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        |
|------------------------|------------------|----------|-----------|-------------|
| <b>Review/Approval</b> | Xiomara Talavera | Quality  |           | 07 APR 2020 |