



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: MSD[®] Human ProInflammatory 9 Plex Tissue Culture Kit

Catalog Number: K15007B-Series

Lot Number: K0035623

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 ProInflammatory 9-Plex Plate	Z0046681	2-8°C	30 JUN 2020
Anti-hProInflammatory 9-Plex Detection Antibody Blend	D0033821	2-8°C	30 NOV 2019
Human Tissue Culture Calibrator 2	A0032316	≤ -70°C	30 NOV 2019
Human IFN-γ Calibrator	A0032308	≤ -70°C	30 SEP 2019

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		28 MAY 2020