



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07855

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049 -2
Lot Number: A0080109
Expiration Date: 31 JAN 2022
Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay									
	hu IFN- γ	hu IL-1 β	hu IL-2	hu IL-4	hu IL-6	hu IL-8	hu IL-10	hu IL-12p70	hu IL-13	hu TNF- α
Reconstituted Calibrator (pg/mL)	1550	599	1660	263	779	657	380	488	506	371

Reconstitution Volume: 1.0 mL

Statement: The above product was manufactured and tested using MSD procedures and it conforms to MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	03 FEB 2017