



### CERTIFICATE OF ANALYSIS

**Product Description:** Immune Response Panel 1 (human) Calibrator Blend  
**Catalog Number:** C0742-2  
**Lot Number:** A0080421  
**Expiration Date:** 31 AUG 2027  
**Storage Temperature:** ≤ -70°C

**Concentration Assignment:**

Analyte	Stock Concentration	Unit of Measure
GM-CSF	6850	pg/mL
IFN-γ	7450	pg/mL
IL-1β	3645	pg/mL
IL-2	10800	pg/mL
IL-6	5150	pg/mL
IL-12p70	8300	pg/mL
IL-10	5200	pg/mL
IL-8	4200	pg/mL
IL-17A	53500	pg/mL
TNF-α	10250	pg/mL

**Dilution to Top of Curve** 5 X

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>	Michelle Likiam	Quality		20 JUN 2025