



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049-2
Lot Number: A0080381
Expiration Date: 30 SEP 2028
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1350	pg/mL
IL-1 β	649	pg/mL
IL-2	1560	pg/mL
IL-4	228	pg/mL
IL-6	738	pg/mL
IL-8	607	pg/mL
IL-10	389	pg/mL
IL-12p70	465	pg/mL
IL-13	556	pg/mL
TNF- α	378	pg/mL

Reconstitution Volume: 1000 μ L

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		09 NOV 2023