



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (mouse) Calibrator Blend
Catalog Number: C0048-2
Lot Number: A0080373
Expiration Date: 31 AUG 2028
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	930	pg/mL
IL-1 β	1700	pg/mL
IL-2	2580	pg/mL
IL-4	1870	pg/mL
IL-5	923	pg/mL
IL-6	5160	pg/mL
KC/GRO	2150	pg/mL
IL-10	3770	pg/mL
IL-12p70	27000	pg/mL
TNF- α	610	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		29 SEP 2023