



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (rat) Calibrator Blend
Catalog Number: C0044-2
Lot Number: A0080405
Expiration Date: 31 AUG 2029
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	5190	pg/mL
IL-1 β	12400	pg/mL
IL-4	1090	pg/mL
IL-5	9510	pg/mL
IL-6	14300	pg/mL
KC/GRO	2720	pg/mL
IL-10	21100	pg/mL
IL-13	1470	pg/mL
TNF- α	1160	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		10 OCT 2024