

**CERTIFICATE OF ANALYSIS**

Control Pack Description: Immune Response Panel 1 (human) Control Pack
Catalog Number: C4742-1
Lot Number: CP000175
Expiration Date: 31 AUG 2027


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Immune Response Panel 1 (human) Control 1	A00C0995	≤ -70°C	31 AUG 2027
Immune Response Panel 1 (human) Control 2	A00C0996	≤ -70°C	31 AUG 2027
Immune Response Panel 1 (human) Control 3	A00C0997	≤ -70°C	31 AUG 2027

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
GM-CSF	655	56.8	5.51	pg/mL
IFN- γ	664	53.2	4.62	pg/mL
IL-1 β	336	29.5	2.81	pg/mL
IL-2	1340	134	15.6	pg/mL
IL-6	459	33.4	2.42	pg/mL
IL-12p70	801	54.1	3.87	pg/mL
IL-10	485	37.8	2.70	pg/mL
IL-8	381	45.1	4.93	pg/mL
IL-17A	4920	378	30.0	pg/mL
TNF- α	1010	86.5	7.24	pg/mL

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		24 JUN 2025