

**CERTIFICATE OF ANALYSIS**

Control Pack Description: Cytolytic Panel 1 (human) Control Pack
Catalog Number: C4741-1
Lot Number: CP000170
Expiration Date: 30 JUN 2027

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Cytolytic Panel 1 (human) Control 1	A00C0986	≤ -70°C	30 JUN 2027
Cytolytic Panel 1 (human) Control 2	A00C0987	≤ -70°C	30 JUN 2027
Cytolytic Panel 1 (human) Control 3	A00C0988	≤ -70°C	30 JUN 2027

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
GM-CSF	608	50.1	4.71	pg/mL
Granzyme B	803	54.1	3.88	pg/mL
IFN- γ	671	52.2	4.51	pg/mL
IL-2	1320	132	14.4	pg/mL
IL-6	542	37.9	3.19	pg/mL
Perforin	64900	4640	376	pg/mL
TNF- α	904	78.1	6.74	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		01 AUG 2025