



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (rat) Control Pack

Catalog Number: C4044-1

Lot Number: CP000056

Expiration Date: 28 FEB 2022

Concentration of Materials Tested:

Units = pg/mL	Assay								
	Rat IFN- γ	Rat IL-1 β	Rat IL-4	Rat IL-5	Rat IL-6	Rat KC/GRO	Rat IL-10	Rat IL-13	Rat TNF- α
Control 1 (A00C0473)	5011	13 088	995	3974	8425	1033	30 052	1502	1200
Control 2 (A00C0474)	1543	3516	345	1834	2819	550	8605	485	357
Control 3 (A00C0475)	477	872	94.2	935	835	284	2946	170	102

Reconstitution Volume: 250 μ L

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (rat) Control 1	A00C0473	2-8°C	28 FEB 2022
Proinflammatory Panel 1 (rat) Control 2	A00C0474	2-8°C	28 FEB 2022
Proinflammatory Panel 1 (rat) Control 3	A00C0475	2-8°C	28 FEB 2022

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	14 JUL 2017