



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (mouse) Control Pack

Catalog Number: C4048-1

Lot Number: CP000072

Expiration Date: 30 APR 2023

Storage Temperature: 2-8°C

Concentration of Materials Tested:

Units = pg/mL	Assay									
	ms IFN- γ	ms IL-1 β	ms IL-2	ms IL-4	ms IL-5	ms IL-6	ms KC/GRO	ms IL-10	ms IL-12p70	ms TNF- α
Reconstituted Control 1 (A00C0579)	467	741	919	480	368	1050	695	1160	13 500	166
Reconstituted Control 2 (A00C0580)	32.2	53.8	54.8	62.0	55.5	167	116	345	3380	29.9
Reconstituted Control 3 (A00C0581)	2.15	4.48	5.84	15.9	8.58	31.5	17.7	136	806	4.93

Reconstitution Volume: 250 μ L

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (mouse) Control 1	A00C0579	2-8°C	30 APR 2023
Proinflammatory Panel 1 (mouse) Control 2	A00C0580	2-8°C	30 APR 2023
Proinflammatory Panel 1 (mouse) Control 3	A00C0581	2-8°C	30 APR 2023

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

	Name	Function	Signature	Date
Review/Approval	Jessica Drolet	Quality Assurance		17 May 2018