

PyroMAT® System Monocyte Activation Test (MAT)

PyroMAT® kit – Ref. PYROMATKIT

Product description

The kit contains reagents for the incubation with the cells and ELISA reagents and plate for Interleukin-6 detection.

Note: the PyroMAT® cells are not included and should be ordered separately, as they have different shipment and storage conditions.

Reagents for Step 1: "PREPARATION & INCUBATION WITH PyroMAT® CELLS"

- Pyrogen-free water
- 96-wells plate
- RPMI

Reagents for Step 2: "DETECTION OF IL-6 WITH ELISA"

- IL-6 microplate
- Assay diluent
- IL- 6 conjugate
- Wash buffer
- Substrate Solution
- Stop solution
- Plate sealers
- IL-6 control

The positive controls (Endotoxin, and Non-Endotoxin Pyrogen positive controls) are to be ordered separately from the kit.

Storage condition

Store unopened kit at 2-8 °C

Quick Procedure

For use of the PyroMAT® kit reagents, refer to the user guide available on the website.

PyroMAT® system products

Product	Catalog number
PyroMAT® cells	PYROMATCELLS
PyroMAT® kit	PYROMATKIT
Reference Standard Endotoxin	1.44161.0001
NEP Control Flagellin	MATFLAGELLIN
NEP Control HKSA	MATHKSA
NEP Control FSL-1	MATFSL1
NEP Control PAM3CSK4	MATPAM3CSK4
NEP Control Resiquimod	MATRESIQUIMOD

To place an order or receive technical assistance

Order/Customer Service:
SigmaAldrich.com/order

Technical Service:
SigmaAldrich.com/techservice

[SigmaAldrich.com/PyroMAT](https://sigmaaldrich.com/PyroMAT)

Merck KGaA
Frankfurter Strasse 250
64293 Darmstadt, Germany