

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"

- Endotoxin-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*
- Color reagent A
- Color reagent B
- Stop solution
- Plate sealers
- *Wash buffer may turn yellow over time. This has no impact on the test.

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	Pyr0MATCELLS
PyroMAT® kit	Pyr0MATKIT
Reference Standard Endotoxin	1.44161.0001
IL-6 Control	Pyr0MATIL6
NEP Control Flagellin	MATFLAGELLIN
NEP Control HKSA	MATHKSA
NEP Control FSL-1	MATFSL1
NEP Control PAM3CSK4	MATPAM3CSK4
NEP Control Resiquimod	MATRESIQUIM0D

To place an order or receive technical assistance

Order/Customer Service: SigmaAldrich.com/order

Technical Service:

SigmaAldrich.com/techservice

SigmaAldrich.com/PyroMAT

Merck KGaA Frankfurter Strasse 250 64293 Darmstadt, Germany

