

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	MATRESIQUIMOD

To place an order or receive technical assistance

Order/Customer Service:
[SigmaAldrich.com/order](https://www.sigmaaldrich.com/order)

Technical Service:
[SigmaAldrich.com/techservice](https://www.sigmaaldrich.com/techservice)

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

MilliporeSigma
400 Summit Drive
Burlington, MA 01803

The life science business of Merck KGaA, Darmstadt, Germany operates as MilliporeSigma in the U.S. and Canada.

© 2022 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved. MilliporeSigma, the vibrant M, Millipore and PyroMAT are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources. Lit. No. MS_PF2111EN Ver. 3.0 04/2022