

ab325586 – Leukocyte-endothelium Adhesion Assay

A robust system for the quantitative determination of leukocyte-endothelium interactions.
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit: www.abcam.com/ab325586

Storage and Stability: 500X Tracker Solution and TNF-alpha should be removed from the kit and stored at -20°C immediately. Store all other components at 4°C. Refer to list of materials supplied for storage conditions of individual components.

Materials Supplied

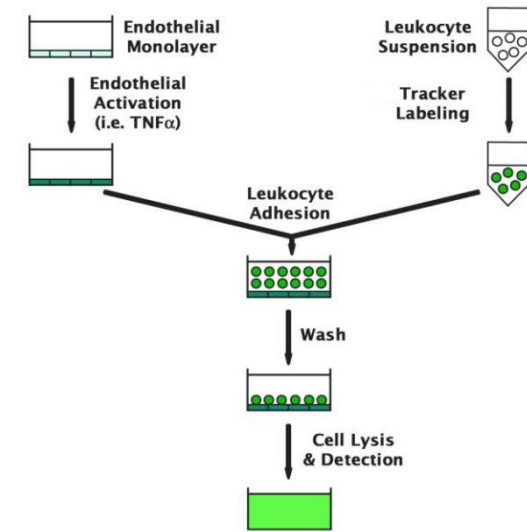
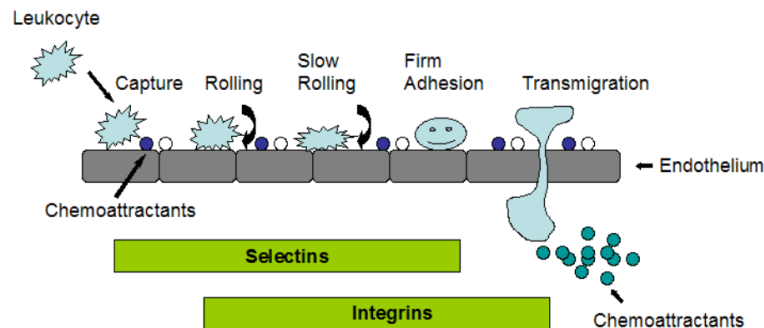
Item	Quantity 100 Tests	Storage Condition
4X Lysis Solution	10 mL	+4°C
500X Tracker Solution	100 µL	-20°C
Gelatin Solution	12 mL	+4°C
10X Wash Buffer	20 mL	+4°C
TNF-alpha	100 µL	-20°C

Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

- Endothelial cells and cell culture medium
- Leukocytes
- 96-well or 48-well tissue culture plate
- Serum free medium, such as DMEM containing 0.5% BSA, 2 mM CaCl₂ and 2 mM MgCl₂
- Sterile 1X PBS
- Cell culture incubator (37°C, 5% CO₂ atmosphere)
- Light microscope
- 96-well plate suitable for a fluorescence plate reader
- Fluorescence plate reader

The Leukocyte-Endothelium Adhesion Cascade



Preparation of Reagents

- **1X Wash Buffer:** Prepare a 1X Wash Buffer by diluting the provided 10X stock 1:10 in deionized water. Store the diluted solution at room temperature.
- **1X Lysis Buffer:** Prepare a 1X Lysis Buffer by diluting the provided 4X stock 1:4 in deionized water. Store the diluted solution at room temperature.

Gelatin Coating

1. Under sterile conditions, add 200 µL of the Gelatin Solution to each well of a 48-well tissue culture treated plate, or 100 µL of the Gelatin Solution to each well of a 96-well tissue culture treated plate.
2. Incubate for 60 min at 37°C in a cell culture incubator.
3. Wash twice with sterile 1X PBS. Aspirate the final wash before use.

Assay Protocol

1. Add 50,000-100,000 endothelial cells/well to the Gelatin-coated 48-well or 96-well plate.
2. Culture cells for 48-72 hrs until the endothelial cells form a monolayer.
3. Treat endothelial cell monolayer or leukocyte with desired activator or inhibitor for 6-12 hrs.
4. Harvest leukocytes and prepare a cell suspension at 1.0×10^6 cells/ml in serum free media. Add Tracker to a final concentration of 1X (for example, add 2 µL of 500X Tracker solution to 1.0 mL of leukocyte cell suspension).
5. Incubate for 60 min at 37°C in a cell culture incubator. Spin down cells at 1000 rpm for 2 minutes, aspirate the medium and wash cell pellet with serum free media. Repeat the wash twice. Resuspend the cell pellet at $0.25 - 1.0 \times 10^6$ cells/ml in serum free media.

6. Aspirate endothelial culture media and wash once with serum free media. Add 200 μ L of the cell suspension to each well already containing the endothelial monolayer.
7. Incubate for 30-90 min in a cell culture incubator.
8. Carefully discard or aspirate the media from each well (**Δ Note:** Do not allow wells to dry). Gently wash each well 3 times with 250 μ L 1X Wash Buffer.
9. (Optional) Count the adherent leukocytes under an inverted fluorescence microscope; average at least three separate fields per well.
10. Aspirate the final wash and add 150 μ L of 1X Lysis Buffer to each well containing cells. Incubate 5 minutes at room temperature with shaking.
11. Transfer 100 μ L of the mixture to a 96-well plate suitable for fluorescence measurement. Read fluorescence with a fluorescence plate reader at 480 nm/520 nm.

Technical Support

Copyright © 2026 Abcam, All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

Version 1 | 2026-03-11

For all technical or commercial enquiries please go to:

<https://www.abcam.com/en-us/contact-us>

<https://www.abcam.cn/contact-us> (China)

<https://www.abcam.co.jp/contact-us> (Japan)