

## ab325574 – Fibronectin Cell Adhesion Assay (Fluorometric)

A rapid, quantitative method for evaluating cell adhesion.  
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit: [www.abcam.com/ab325574](http://www.abcam.com/ab325574)

**Storage and Stability:** Store kit at 2-8°C immediately upon receipt. Refer to list of materials supplied for storage conditions of individual components.

### Materials Supplied

Item	Quantity 48 Tests	Storage Condition
Fibronectin Adhesion Plate	1 unit	+4°C
4X Lysis Solution	10 mL	+4°C
GR Dye	50 µL	+4°C

### Adhesion Plate Layout

The layout below indicates the location of wells coated with Fibronectin and those coated with BSA.

	1	2	3	4	5	6	7	8
A	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin
B	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin
C	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin
D	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin
E	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin	Fibronectin
F	BSA	BSA	BSA	BSA	BSA	BSA	BSA	BSA

### Materials Required, Not Supplied

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

- Cell culture medium
- Serum free medium, such as DMEM containing 0.5% BSA, 2 mM CaCl<sub>2</sub> and 2 mM MgCl<sub>2</sub>
- Cell culture incubator (37°C, 5% CO<sub>2</sub> atmosphere)
- 1X PBS containing 2 mM CaCl<sub>2</sub> and 2 mM MgCl<sub>2</sub>
- Light microscope
- 96-well plate suitable for a fluorescence plate reader
- Fluorescence plate reader

### Assay Protocol

1. Under sterile conditions, allow the Fibronectin Adhesion Plate to warm up at room temperature for 10 minutes.
2. Prepare a cell suspension containing 0.1-1.0 x 10<sup>6</sup> cells/mL in serum free media. Agents that inhibit or stimulate cell adhesion can be added directly to the cell suspension.
3. Add 150 µL of the cell suspension to the inside of each well (BSA-coated wells are provided as a negative control).
4. Incubate for 30-90 min in a cell culture incubator.
5. Carefully discard or aspirate the media from each well (Note: Do not allow wells to dry). Gently wash each well 4-5 times with 250 µL PBS.
6. Prepare sufficient 1X Lysis Buffer/GR dye solution for all samples by diluting the dye 1:300 in Lysis Buffer (for example, add 4 µL dye to 300 µL of 4X Lysis Buffer and 900 µL of dH<sub>2</sub>O).
7. Add 200 µL of 1X Lysis Buffer/GR dye solution to each well containing cells. Incubate 20 minutes at room temperature with shaking.
8. Transfer 150 µL 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 | 2026-03-10

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)