
Product Manual

CytoSelect™ 48-Well Cell Adhesion Assay (Laminin-Coated, Fluorometric Format)

Catalog Number

CBA-057

48 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.
Creating Solutions for Life Science Research

Introduction

Cell adhesion is a complex process involved in embryogenesis, migration/invasion, tissue remodeling, and wound healing. To perform these processes, cells adhere to extracellular matrix components (via adhesion receptors), forming complexes with components of the cytoskeleton that ultimately affect cell motility, differentiation, proliferation, and survival. The Cell Biolabs CytoSelect™ Cell Adhesion Assay Kit provides a rapid, quantitative method for evaluating cell adhesion. The kit contains sufficient reagents for the evaluation of 48 samples (40 Mouse Laminin I-coated wells, 8 BSA-coated wells).

Assay Principle

The CytoSelect™ Cell Adhesion Assay Kit utilizes a Laminin-coated 48-well plate (see Adhesion Plate Layout). First, cells are seeded onto the coated substrate, where the adherent cells are captured. Next, unbound cells are removed with consecutive washes. Finally, the adherent cells are lysed and subsequently detected with CyQuant® GR Dye.

Related Products

1. CBA-050: CytoSelect™ 48-Well Cell Adhesion Assay, (Fibronectin-Coated, Colorimetric Format)
2. CBA-051: CytoSelect™ 48-Well Cell Adhesion Assay (Fibronectin-Coated, Fluorometric Format)
3. CBA-052: CytoSelect™ 48-Well Cell Adhesion Assay (Collagen I-Coated, Colorimetric Format)
4. CBA-053: CytoSelect™ 48-Well Cell Adhesion Assay (Collagen I-Coated, Fluorometric Format)
5. CBA-056: CytoSelect™ 48-Well Cell Adhesion Assay (Laminin-Coated, Colorimetric Format)
6. CBA-058: CytoSelect™ 48-Well Cell Adhesion Assay (Fibrinogen-Coated, Colorimetric Format)
7. CBA-059: CytoSelect™ 48-Well Cell Adhesion Assay (Fibrinogen-Coated, Fluorometric Format)
8. CBA-060: CytoSelect™ 48-Well Cell Adhesion Assay (Collagen IV-Coated, Colorimetric Format)
9. CBA-061: CytoSelect™ 48-Well Cell Adhesion Assay (Collagen IV-Coated, Fluorometric Format)
10. CBA-070: CytoSelect™ 48-Well Cell Adhesion Assay (ECM Array, Colorimetric Format)
11. CBA-071: CytoSelect™ 48-Well Cell Adhesion Assay (ECM Array, Fluorometric Format)

Kit Components

1. Laminin Adhesion Plate (Part No. 105601): One 48-well plate containing 40 Mouse Laminin I-coated wells and 8 BSA-coated wells (see layout below)
2. 4X Lysis Buffer (Part No. 10404): One Bottle – 10.0 mL
3. CyQuant® GR Dye (Part No. 105101): One tube – 50 µL

Adhesion Plate Layout

The following layout indicates the location of wells coated with Laminin and those coated with BSA.

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

Materials Not Supplied

1. Cell culture medium
2. Serum free medium, such as DMEM containing 0.5% BSA, 2 mM CaCl₂ and 2 mM MgCl₂
3. Cell culture incubator (37°C, 5% CO₂ atmosphere)
4. 1X PBS containing 2 mM CaCl₂ and 2 mM MgCl₂
5. Light microscope
6. 96-well plate suitable for a fluorescence plate reader
7. Fluorescence plate reader

Storage

Store all kit components at 4°C.

Assay Protocol

1. Under sterile conditions, allow the Laminin Adhesion Plate to warm up at room temperature for 10 minutes.
2. Prepare a cell suspension containing 0.1-1.0 x 10⁶ 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 μ L PBS.
6. Prepare sufficient 1X Lysis Buffer/CyQuant® GR dye solution for all samples by diluting the dye 1:300 in Lysis Buffer (for example, add 4 μ L dye to 300 μ L of 4X Lysis Buffer and 900 μ L of dH₂O).
7. Add 200 μ L of 1X Lysis Buffer/CyQuant® GR dye solution to each well containing cells. Incubate 20 minutes at room temperature with shaking.
8. Transfer 150 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.

Example of Results

The following figures demonstrate typical results with the CytoSelect™ 48-Well Cell Adhesion Assay Kit. Fluorescence measurement was performed on SpectraMax Gemini XS Fluorometer (Molecular Devices) with a 485/538 nm filter set and 530 nm cutoff. One should use the data below for reference only. This data should not be used to interpret actual results.

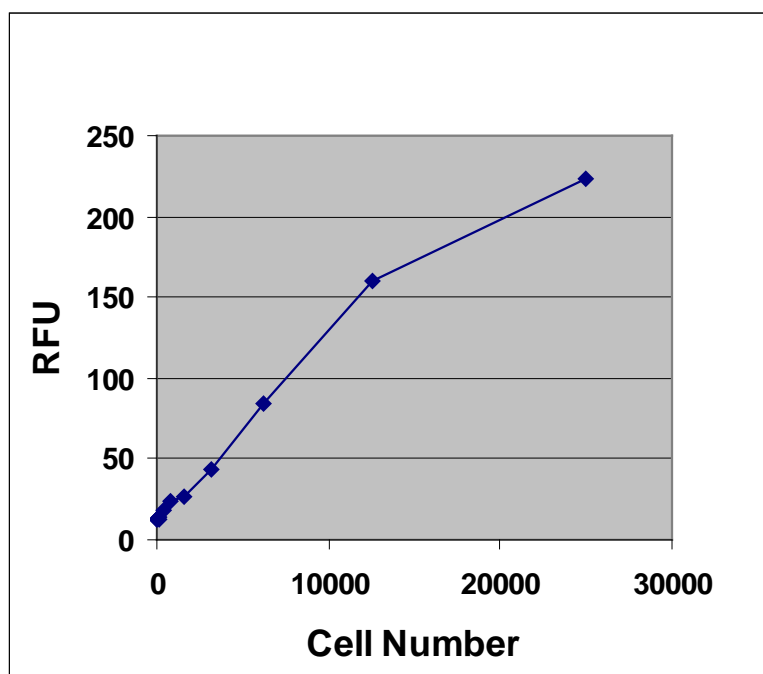


Figure 1: Quantitation of Human HT-1080. HT-1080 cells were titrated in 1X PBS, then subsequently lysed and detected with 1X Lysis Buffer/CyQuant® GR Dye (225 μ L of cell suspension was mixed with 75 μ L of 4X Lysis Buffer and 1 μ L of CyQuant® GR dye).

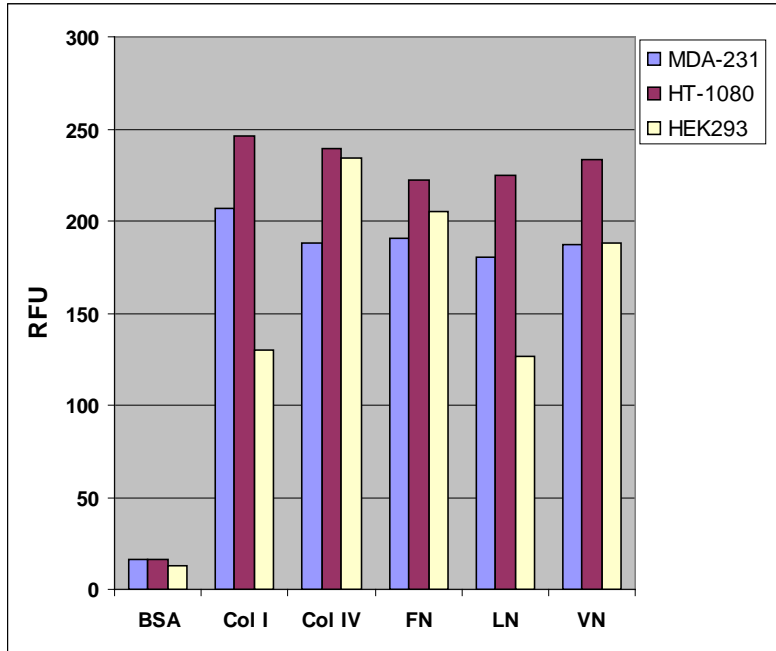


Figure 2. ECM-mediated Cell Adhesion. Serum starved cells were allowed to attach to ECM-coated 48-well plate for 1 hr at 100,000 cells/well. Adherent cells were lysed and quantified by CyQuant® GR Dye as described in the Assay Protocol.

References

1. Hynes, R. O. (1992) *Cell* **69**, 11-25.
2. Schwartz, M. A., Schaller, M. D. and Ginsberg, M. H. (1995) *Annu. Rev. Cell Dev. Biol.* **11**, 549-599.

Recent Product Citation

Lee, J. et al. (2013). Selective inhibition of prostaglandin E2 receptors EP2 and EP4 inhibits adhesion of human endometriotic epithelial and stromal cells through suppression of integrin-mediated mechanisms. *Biol Reprod.* **88**:77.

License Information

This product is provided under an intellectual property license from Life Technologies Corporation. The purchase of this product conveys to the buyer the non-transferable right to use the purchased product and components of the product only in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The sale of this product is expressly conditioned on the buyer not using the product or its components, or any materials made using the product or its components, in any activity to generate revenue, which may include, but is not limited to use of the product or its components: (i) in manufacturing; (ii) to provide a service, information, or data in return for payment; (iii) for therapeutic, diagnostic or prophylactic purposes; or (iv) for resale, regardless of whether they are resold for use in research. For information on purchasing a license to this product for purposes

other than as described above, contact Life Technologies Corporation, 5791 Van Allen Way, Carlsbad CA 92008 USA or outlicensing@lifetech.com.

Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS' sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products

Contact Information

Cell Biolabs, Inc.
7758 Arjons Drive
San Diego, CA 92126
Worldwide: +1 858-271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2004-2016: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.