

eBioscience™ Annexin V Apoptosis Detection Kit APC

Catalog Number: 88-8007

For Research Use Only. Not for use in diagnostic procedures.

Product Information

REF

Contents: eBioscience™ Annexin V
Apoptosis Detection Kit APC
Catalog Number: 88-8007
Concentration: 5 µL/test



LOT



Formulation: aqueous buffer, 0.09% sodium azide, may contain carrier protein/stabilizer
Temperature Limitation: Store at 2-8°C. Do not freeze. Light sensitive material.
Batch Code: Refer to vial
Use By: Refer to vial
Contains sodium azide

Description

Annexins are a family of calcium-dependent phospholipid-binding proteins that preferentially bind phosphatidylserine (PS). Under normal physiologic conditions, PS is predominantly located in the inner leaflet of the plasma membrane. Upon initiation of apoptosis, PS loses its asymmetric distribution across the phospholipid bilayer and is translocated to the extracellular membrane leaflet marking cells as targets of phagocytosis. Once on the outer surface of the membrane, PS can be detected by fluorescently labeled Annexin V in a calcium-dependent manner.

In early-stage apoptosis, the plasma membrane excludes viability dyes such as propidium iodide (PI), 7-AAD, or Fixable Viability Dyes such as eFluor® 660 or eFluor® 780. These cells will stain with Annexin V but not a viability dye, thus distinguishing cells in early apoptosis. However, in late stage apoptosis, the cell membrane loses integrity thereby allowing Annexin V to also access PS in the interior of the cell. A viability dye can be used to resolve these late-stage apoptotic and necrotic cells (Annexin V, viability dye-positive) from the early-stage apoptotic cells (Annexin V positive, viability dye-negative).

Note: Fixable Viability Dye eFluor® 450 is not recommended for use with Annexin V Apoptosis Detection Kits.

Components

Cat. No. 88-8007-72:

10X Binding Buffer (cat. 00-0055): 30 mL, store at 2-8°C.

Annexin V APC (cat. 17-8007): 5 µL/test, 50 tests, store at 2-8°C. Protect from light.

Propidium Iodide Staining Solution (cat. 00-6990): 5 µL/test, 100 tests, store at 2-8°C. Protect from light.

Cat. No. 88-8007-74:

10X Binding Buffer (cat. 00-0055): 100 mL, store at 2-8°C.

Annexin V APC (cat. 17-8007): 5 µL/test, 200 tests, store at 2-8°C. Protect from light.

Propidium Iodide Staining Solution (cat. 00-6990): 5 µL/test, 2x100 tests, store at 2-8°C. Protect from light.

Applications Reported

The Annexin V Apoptosis Detection Kit APC has been reported for use in flow cytometric analysis.

Applications Tested

The Annexin V Apoptosis Detection Kit APC has been pre-titrated and tested on mouse thymocytes cultured overnight in medium (to induce apoptosis) or on Jurkat cells treated with 10 µM Camptothecin for 4 hours. Both the Annexin V APC and the Propidium iodide components can be used at 5 µL per test. A test is defined as the amount (µg) of antibody that will stain a cell sample in a final volume of 100 µL. Cell number should be determined empirically but can range from 10⁵ to 10⁸ cells/test.

Special Notes

The Annexin V Apoptosis Detection Kits are compatible with intracellular staining. Please refer to the Best Protocols: Annexin V Staining Protocol, Protocol C for details.

Not for further distribution without written consent.

Copyright © 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Tel: 888.999.1371 or 858.642.2058 • Fax: 858.642.2046 • thermofisher.com/ebioscience •

info@ebioscience.com

eBioscience™ Annexin V Apoptosis Detection Kit APC

Catalog Number: 88-8007

For Research Use Only. Not for use in diagnostic procedures.

References

Andree HA, Reutelingsperger CP, Hauptmann R, Hemker HC, Hermens WT, Willems GM. Binding of vascular anticoagulant alpha (VAC alpha) to planar phospholipid bilayers. *J Biol Chem.* 1990; 265(9):4923-4928

Koopman G, Reutelingsperger CP, Kuijten GA, Keehnen RM, Pals ST, van Oers MH. Annexin V for flow cytometric detection of phosphatidylserine expression on B cells undergoing apoptosis. *Blood.* 1994; 84(5):1415-1420

Vermes I, Haanen C, Steffens-Nakken H, Reutelingsperger C. A novel assay for apoptosis. Flow cytometric detection of phosphatidylserine expression on early apoptotic cells using fluorescein labelled Annexin V. *J Immunol Methods.* 1995; 184(1):39-51

Related Products

00-6993 eBioscience™ 7-AAD Viability Staining Solution

65-0864 eBioscience™ Fixable Viability Dye eFluor™ 660

65-0865 eBioscience™ Fixable Viability Dye eFluor™ 780

88-8005 eBioscience™ Annexin V Apoptosis Detection Kit FITC

88-8006 eBioscience™ Annexin V Apoptosis Detection Kit eFluor™ 450

88-8008 eBioscience™ Annexin V Apoptosis Detection Kit PerCP-eFluor™ 710

88-8102 eBioscience™ Annexin V Apoptosis Detection Kit PE

Legal

Pat. No. EP 181 465 B2, EP 0509 026, USP 5,066,787

Not for further distribution without written consent.

Copyright © 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Tel: 888.999.1371 or 858.642.2058 • Fax: 858.642.2046 • thermofisher.com/ebioscience •

info@ebioscience.com

Annexin V Staining Protocols

Protocol: Annexin V Staining

Note: Due to the calcium dependence of the Annexin V:PS interaction, it is critical to avoid buffers containing EDTA or other calcium chelators during Annexin V experiments. Annexin V can only be used as a marker of apoptosis in cells where the plasma membrane is intact because destroying the integrity of the plasma membrane will allow non-specific binding of Annexin V to PS inside the cell.

Experimental Procedure

1. Dilute 10X Binding Buffer to 1X using distilled water (1 mL 10X Binding Buffer + 9 mL dH₂O).
2. Wash cells once in PBS, then once in 1X Binding Buffer.
3. Resuspend cells in 1X Binding Buffer at 1-5x10⁶/mL.
4. Add 5 µL of fluorochrome-conjugated Annexin V to 100 µL of the cell suspension.
5. Incubate 10-15 minutes at room temperature.
6. Wash cells in 1X Binding Buffer and resuspend in 200 µL of 1X Binding Buffer.
7. Add 5 µL of Propidium Iodide Staining Solution (cat. 00-6990) or 7-AAD Viability Staining Solution (cat. 00-6993).
8. Analyze by flow cytometry within 4 hours, storing at 2-8°C in the dark

Protocol: Annexin V Staining with Fixable Viability Dyes

Note: Due to the calcium dependence of the Annexin V:PS interaction, it is critical to avoid buffers containing EDTA or other calcium chelators during Annexin V experiments. Annexin V can only be used as a marker of apoptosis in cells where the plasma membrane is intact because destroying the integrity of the plasma membrane will allow non-specific binding of Annexin V to PS inside the cell.

Materials Needed

- PBS without sodium azide
- Fixable Viability Dyes
 - Fixable Viability Dye eFluor™ 455UV (Cat. No. 65-0868)
 - Fixable Viability Dye eFluor™ 450 (Cat. No. 65-0863)
 - Fixable Viability Dye eFluor™ 520 (Cat. No. 65-0867)
 - Fixable Viability Dye eFluor™ 660 (Cat. No. 65-0864)
 - Fixable Viability Dye eFluor™ 780 (Cat. No. 65-0865)
- Distilled water
- Flow Cytometry Staining Buffer (Cat. No. 00-4222)

Experimental Procedure

1. Choose an appropriate viability stain that has an emission profile compatible with the Annexin V- conjugate to be used.
Note: Fixable Viability Dye eFluor™ 450 is not recommended for use with the Annexin V Apoptosis Detection Kits.
2. Follow the staining protocol for the chosen Fixable Viability Dye to stain late-apoptotic/dead cells.
Refer to the Best Protocols webpage (Viability Staining Protocol C in the Resources tab of the home page).
3. After staining with Fixable Viability Dye, be sure to wash cells twice with a protein-containing buffer such as Flow Cytometry Staining Buffer (cat. 00-4222).
4. Dilute 10X Binding Buffer to 1X using distilled water (1 mL 10X Binding Buffer + 9 mL dH₂O).
5. Wash cells once with the 1X Binding Buffer.
6. Resuspend cells in 1X Binding Buffer at 1-5x10⁶/mL.
7. Add 5 µL of fluorochrome-conjugated Annexin V to 100 µL of the cell suspension.
8. Incubate 10-15 minutes at room temperature, protected from light.
9. Wash cells in 1X Binding Buffer and resuspend in 200 µL of 1X Binding Buffer.
10. Analyze by flow cytometry within 4 hours, storing at 2-8°C in the dark.

Documentation and support

Customer and technical support

Visit thermofisher.com/support for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
 - Product FAQs
 - Software, patches, and updates
- Order and web support
- Product documentation, including:
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at thermofisher.com/support.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Corporate entity: Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. All other trademarks are properties of their respective owners.

For support visit thermofisher.com/support or email techsupport@lifetech.com

thermofisher.com

23 January 2017

ThermoFisher
SCIENTIFIC