

Annexin V-FITC Apoptosis detection Kit

PRODUCT INFORMATION & MANUAL

Catalog Number  *BMS500FI/20CE, BMS500FI/100CE, BMS500FI/300CE*



For in-vitro diagnostic use. Not for therapeutic procedures.

Information in this document is subject to change without notice.

DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC 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

Trademarks

All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. All other trademarks are the property of their respective owners.

© 2017 Thermo Fisher Scientific Inc. All rights reserved.



Bender MedSystems GmbH Campus Vienna Biocenter 2 1030 Vienna, Austria

TABLE OF CONTENTS

1 Intended Use	3
2 Summary	3
3 Principles of the Test	3
4 Reagents Provided	3
5 Storage and Stability	3
6 Materials Required But Not Provided	4
7 Precautions for Use	4
8 Preparation of Reagents	5
9 Test Protocol	5
10 Limitations	5
Documentation and support.....	6

1 Intended Use

The Annexin V-FITC Kit can be used to detect phosphatidylserine on the outer leaflet of the cell membrane using flow cytometry.

The Annexin V-FITC Kit is for in vitro diagnostic use. Not for use in therapeutic procedures.

2 Summary

Annexins are a family of calcium-dependent phospholipid-binding proteins. They are abundant in eukaryotic organisms belonging to a family of ubiquitous cytoplasmic proteins involved in signal transduction. All annexins have been shown to have a putative binding site for protein kinase C (PKC) but only annexin V would possess a potential pseudo- substrate site. Thus annexin V seems to modulate the activity of some PKCs on their substrates.

Annexin V was found to play a major role in matrix vesicle-initiated cartilage calcification as a collage-regulated calcium channel. Annexin V binds to procoagulant phospholipids (Vascular anticoagulant alpha) with high affinity.

Annexin V's preferential binding partner is phosphatidylserine (PS). PS is predominantly located in membrane leaflets, which face the cytosol. However, recent findings show that each cell type has the molecular machinery to expose PS at its cell surface. This machinery is activated during the execution of apoptosis. Once PS is exposed at the cell surface it exhibits procoagulant and proinflammatory activities. Annexin V will bind to the PS-exposing apoptotic cell and can inhibit the procoagulant and proinflammatory activities of the dying cell.

For literature update refer to www.thermofisher.com.

3 Principles of the Test

Annexin V exhibits anti-phospholipase activity and binds to phosphatidylserine. FITC labelling allows simple direct detection by FACS analysis. Counterstaining by propidium iodide allows the discrimination of apoptotic cells.

4 Reagents Provided

- **rh Annexin V-FITC**
- **Binding Buffer (4x)**
- **Propidium Iodide (20 µg/ml)**

	Catalog Number		
	BMS500FI/20CE for 20 tests	BMS500FI/100CE for 100 tests	BMS500FI/300CE for 300 tests
rh Annexin V-FITC	1 vial (100µl)	1 vial (500µl)	1 vial (1.5 ml)
Binding Buffer (4x)	1 bottle (50 ml)	1 bottle (50 ml)	1 bottle (50 ml)
Propidium Iodide	1 vial (1.6 ml)	1 vial (1.6 ml)	2 vials (1.6 ml)

5 Storage and Stability

Store kit reagents between 2° and 8°. Immediately after use remaining reagents should be returned to cold storage (2° to 8°C). Expiry of the kit and reagents is stated on labels.

Expiry of the kit components can only be guaranteed if the components are stored properly, and if, in case of repeated use of one component, this reagent is not contaminated by the first handling.

6 Materials Required But Not Provided

- 5 ml and 10 ml graduated pipettes
- 5 µl to 1000 µl adjustable single channel micropipettes with disposable tips
- Beakers, flasks, cylinders necessary for preparation of reagents
- Glass-distilled or deionized water
- Bench top centrifuge
- Flow Cytometer
- PBS (for 1 liter: 8.00 g NaCl, 0.20 g KCl, 1.78 g Na₂HPO₄ x 2 H₂O, 0.27 g KH₂PO₄)

7 Precautions for Use

- All reagents should be considered as potentially hazardous. We therefore recommend that this product is handled only by those persons who have been trained in laboratory techniques and that it is used in accordance with the principles of good laboratory practice. Wear suitable protective clothing such as laboratory overalls, safety glasses and gloves. Care should be taken to avoid contact with skin or eyes. In the case of contact with skin or eyes wash immediately with water. See material safety data sheet(s) and/or safety statement(s) for specific advice.
- Reagents are intended for in vitro diagnostic use only and are not for use in therapeutic procedures.
- Do not mix or substitute reagents with those from other lots or other sources.
- Do not use kit reagents beyond expiration date on label.
- Do not expose kit reagents to strong light during storage or incubation.
- Do not pipette by mouth.
- Do not eat or smoke in areas where kit reagents or samples are handled.
- Avoid contact of skin or mucous membranes with kit reagents or specimens.
- Rubber or disposable latex gloves should be worn while handling kit reagents or specimens.
- Avoid splashing or generation of aerosols.
- In order to avoid microbial contamination or cross-contamination of reagents which may invalidate the test use disposable pipette tips and/or pipettes.
- Glass-distilled water or deionized water must be used for reagent preparation.
- Decontaminate and dispose specimens and all potentially contaminated materials as they could contain infectious agents. The preferred method of decontamination is autoclaving for a minimum of 1 hour at 121.5°C.
- Liquid wastes not containing acid and neutralized waste may be mixed with sodium hypochlorite in volumes such that the final mixture contains 1.0% sodium hypochlorite. Allow 30 minutes for effective decontamination. Liquid waste containing acid must be neutralized prior to the addition of sodium hypochlorite.

8 Preparation of Reagents

Dilute **Binding Buffer** (4x) 1:4 in distilled water (50 ml binding buffer and 150 ml distilled water).

9 Test Protocol

1. Wash cells in **PBS** by gentle shaking or pipetting up and down.
2. Resuspend cells in 200 μ l **Binding Buffer** (1x); cell density should be 2-5x10⁵/ml.
3. Add 5 μ l **Annexin V-FITC** to 195 μ l cell suspension.
4. Mix and incubate for 10 min at room temperature.
5. Wash cells with 200 μ l Binding Buffer (1x) and resuspend in 190 μ l Binding buffer (1x).
6. Add 10 μ l **Propidium Iodide** (20 μ g/ml)
7. Perform FACS analysis.

10 Limitations

- Bacterial or fungal contamination of either screen samples or reagents or cross-contamination between reagents may cause erroneous results.
- Disposable pipette tips, flasks or glassware are preferred, reusable glassware must be washed and thoroughly rinsed of all detergent before use.

Documentation and support

Obtaining support

Technical support	<p>For the latest services and support information for all locations, visit www.thermofisher.com.</p> <p>At the website, you can:</p> <ul style="list-style-type: none">• Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities• Search through frequently asked questions (FAQs)• Submit a question directly to Technical Support (thermofisher.com/support)• Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents• Obtain information about customer training• Download software updates and patches
Safety Data Sheets (SDS)	<p>Safety Data Sheets (SDSs) are available at thermofisher.com/support.</p>
Limited product warranty	<p>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 www.thermofisher.com/support.</p>

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

thermofisher.com

23 January 2017

ThermoFisher
SCIENTIFIC