

Product Contents

CaspACE™ FITC-VAD-FMK In Situ Marker:

Part No.	Size
G746A	50µl
G746B	125µl

Description: CaspACE™ FITC-VAD-FMK In Situ Marker is a fluoroisothiocyanate (FITC) conjugate of the cell permeable caspase inhibitor VAD-FMK. This structure allows delivery into the cell where it binds to activated caspase, serving as an in situ marker for apoptosis. The bound marker is localized by fluorescence detection.

Formulation: CaspACE™ FITC-VAD-FMK In Situ Marker is supplied as a 5mM solution in DMSO.

Storage Conditions: Store aliquots at -20°C protected from light and moisture, preferably in a desiccator. Avoid multiple freeze-thaw cycles and exposure to frequent temperature changes.

Usage Notes: Concentration gradients may form in frozen products and should be dispersed upon thawing. Mix well prior to use.

Usage Information

Protocol for use of CaspACE™ FITC-VAD-FMK In Situ Marker

1. Seed Jurkat cells at 1×10^5 cells/ml and grow in RPMI-1640 + 10% FBS in a 37°C, 5% CO₂ incubator for 2–3 days, until they reach a density of 5×10^5 cells/ml.
2. Prepare poly-L-lysine-coated slides. A 0.01% poly-L-lysine solution is used to coat each chamber on multi-chamber slides. When partially dry, the slides are rinsed in NANOpure® water and then allowed to air-dry. Poly-L-lysine-coated slides can be prepared in advance and stored at 4°C for up to 7 days prior to use.
3. Add anti-Fas mAb (Clone CH-11, PanVera Cat. #SY-100) to a final concentration of 0.1µg/ml. Do not add to controls. Incubate for 3–4 hours at 37°C.
4. Add CaspACE™ FITC-VAD-FMK In Situ Marker to the Jurkat cells at a final concentration of 10µM. Protect cells from light and incubate for 20 minutes in the incubator. Keep cells protected from light for the remaining steps.
5. Centrifuge at 300 ×g for 5 minutes.
6. Wash cells with PBS then centrifuge at 300 ×g for 5 minutes.
7. Suspend cells in PBS to 1.5×10^6 cells/ml.
8. Add cells to poly-L-lysine-coated slides and incubate at room temperature for 5 minutes to allow the cells to adhere to the poly-L-lysine.
9. Fix in 10% buffered formalin for 30 minutes at room temperature (protected from light).
10. Rinse 3 × 5 minutes in PBS.
11. Add mounting medium and cover slips to the slides. Analyze under a fluorescence microscope.

Part# 9PIG746

Revised 10/03



Promega

Promega Corporation

2800 Woods Hollow Road	
Madison, WI 53711-5399	USA
Telephone	608-274-4330
Toll Free	800-356-9526
Fax	608-277-2516
Internet	www.promega.com

PRODUCT USE LIMITATIONS, WARRANTY, DISCLAIMER

Promega manufactures products for a number of intended uses. Please refer to the product label for the intended use statements for specific products. Promega products contain chemicals which may be harmful if misused. Due care should be exercised with all Promega products to prevent direct human contact.

Each Promega product is shipped with documentation stating specifications and other technical information. Promega products are warranted to meet or exceed the stated specifications. Promega's sole obligation and the customer's sole remedy is limited to replacement of products free of charge in the event products fail to perform as warranted. Promega makes no other warranty of any kind whatsoever, and SPECIFICALLY DISCLAIMS AND EXCLUDES ALL OTHER WARRANTIES OF ANY KIND OR NATURE WHATSOEVER, DIRECTLY OR INDIRECTLY, EXPRESS OR IMPLIED, INCLUDING, WITHOUT LIMITATION, AS TO THE SUITABILITY, PRODUCTIVITY, DURABILITY, FITNESS FOR A PARTICULAR PURPOSE OR USE, MERCHANTABILITY, CONDITION, OR ANY OTHER MATTER WITH RESPECT TO PROMEGA PRODUCTS. In no event shall Promega be liable for claims for any other damages, whether direct, incidental, foreseeable, consequential, or special (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including negligence) or strict liability arising in connection with the sale or the failure of Promega products to perform in accordance with the stated specifications.

CaspACE is a trademark of Promega Corporation.

© 1999–2003 Promega Corporation. All Rights Reserved.

All specifications are subject to change without prior notice.

NANOpure is a registered trademark of Barnstead/ThermoLyne Corporation

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.

Part# 9PIG746
Revised 10/03.