

**Ac-DEVD-pNA \*CAS 189950-66-1\***

Catalog number: 13405

Unit size: 5 mg

Component	Storage	Amount
Ac-DEVD-pNA *CAS 189950-66-1*	Freeze (<-15 °C), Minimize light exposure	5 mg

**OVERVIEW**

Ac-DEVD-pNA is a colorimetric substrate for caspase 3, a protease that is rapidly activated when cells are exposed to apoptotic conditions and that cleaves poly(ADP-ribose) polymerase.

**AT A GLANCE**
**Important notes**

It is important to store at <-15 °C and should be stored in cool, dark place.

It can be used within 12 months from the date of receipt.

**PREPARATION OF STOCK SOLUTIONS**

Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20 °C after preparation. Avoid repeated freeze-thaw cycles.

*Ac-DEVD-pNA stock solution (100 mM):*

Add 80 µL of DMSO to 5 mg to make Ac-DEVD-pNA stock solution.

**Note** The stock solution should be used promptly; any unused solution should be aliquoted and refrozen at -20 °C. Avoid repeated freeze-thaw cycles and keep from light.

**PREPARATION OF WORKING SOLUTION**

*Ac-DEVD-pNA working solution (0.2 mM):*

Add assay buffer of your choice to Ac-DEVD-pNA stock solution to make working solution.

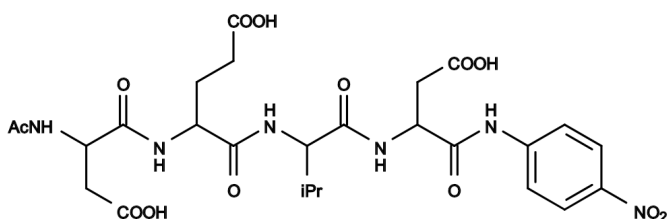
**Note** Buffer can be used such as 50 mM Tris-HCl, pH 7.5, 0.3% NP-40, 1.0 mM DTT)

**SAMPLE EXPERIMENTAL PROTOCOL**

1. Add 0.01 mL of enzyme solution (or cell lysate\*) into 0.09 mL assay solution (from Step 2) to make a final volume of 0.1 mL.

**Note** Lyse cells in 50 mM Tris-HCl, pH 7.5, 0.3% NP-40, 1.0 mM DTT, at a density of  $2 \times 10^6$ /mL).

2. Incubate the assay solution at 37 °C for 0-3 hr. Take periodic readings of absorbance at 405 nm.

**EXAMPLE DATA ANALYSIS AND FIGURES**


**Figure 1.** Chemical structure for Ac-DEVD-pNA \*CAS 189950-66-1\*

**DISCLAIMER**

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email [info@aatbio.com](mailto:info@aatbio.com) if you have any questions.