

**Z-DEVD-ProRed™ 620**

Catalog number: 13433

Unit size: 1 mg

Component	Storage	Amount
Z-DEVD-ProRed™ 620	Freeze (<-15 °C), Minimize light exposure	1 mg

**OVERVIEW**

ProRed™-derived protease substrates are colorless and non-fluorescent. Cleavage of blocking protease-cleavable peptide residue by caspases generates the strongly red fluorescent ProRed™ that can be monitored fluorimetrically at ~620 nm with excitation of ~530 nm. ProRed™-derived caspase substrates are the most sensitive red indicators for the fluorimetric detection of various caspase activities. This DEVD-ProRed™ substrate is specific for detecting caspases 3 and 7.

**AT A GLANCE**

**Important** 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.*

**1. Z-DEVD-ProRed™ 620 stock solution (10 mM):**

Add 65 µL of DMSO into the vial of 1 mg Z-DEVD-ProRed™ 620 to make 10 mM stock solution.

**PREPARATION OF WORKING SOLUTION****Caspase 3/7 assay solution (2X):**

Mix 50 µL Z-DEVD-ProRed™ 620 stock solution (10 mM), 100 µL DTT (1M), 400 µL EDTA (100 mM) and 10 mL Tris Buffer (20 mM), pH =7.4.

**SAMPLE EXPERIMENTAL PROTOCOL**

1. Mix equal volume of the caspase 3/7 standards or samples with 2X caspase 3/7 assay reaction solution and incubate at room temperature for at least 1 hour.
2. Monitor the fluorescence increase at Ex/Em = 535/620 nm.

**EXAMPLE DATA ANALYSIS AND FIGURES**

Example data analysis and images of this product can be found on the web at:

<https://www.aatbio.com/products/z-devd-prored-620>

**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.