# Z-DEVD-ProRed<sup>™</sup> 620

Ordering Information

Product Number: 13433 (1 mg)

#### **General Properties**

Molecular Weight: 1565.50 Maximum excitation: 534 nm Maximum Emission: 619 nm Solvents: DMSO

## **Biological Applications**

Caspases play important roles in apoptosis and cell signaling. The activation of Caspase 3/7 (CPP32/apopain) is important for the initiation of apoptosis. Caspase 3/7 is also identified as a drug-screening target. Caspase inhibitors have anti-cancer and other pharmalogical potentials. It has been proven that Caspase 3/7 has substrate selectivity for the peptide sequence Asp-Glu-Val-Asp (DEVD). ProRed<sup>TM</sup>-derived protease substrates are colorless and non-fluorescent. Cleavage of blocking protease-cleavable peptide residue by caspases generates the strongly red fluorescent ProRed<sup>TM</sup> that can be monitored fluorimetrically at ~620 nm with excitation of ~535 nm. ProRed<sup>TM</sup>-derived caspase substrates are the most sensitive red indicators for the fluorimetric detection of various caspase activities. This DEVD-ProRed<sup>TM</sup> substrate is specific for detecting caspases 3 and 7.

## **Sample Protocol**

- 1. Make a 10 mM stock solution by adding 65 µL of DMSO into the vial of 1 mg Z-DEVD-ProRed<sup>™</sup> 620
- Prepare 2X Caspase 3/7 assay solution as the following: 50 μL (Z-DEVD)<sub>2</sub>-R110 stock solution (10 mM) 100 μL DTT (1M) 400 μL EDTA (100 mM) 10 mL Tris Buffer (20 mM), pH =7.4
- **3.** Mix equal volume of the caspase 3/7 standards or samples with 2X caspase 3/7 assay solution, and incubate at room temperature for at least 1 hour.
- 4. Monitor the fluorescence increase at Ex/Em = 535/620 nm.

## **References**

- 1. N. A. Thornberry and Y. Lazebnik, *Science* 281, 1312-1316 (1998).
- 2. J. C. Reed, J.Clin.Oncol. 17, 2941-2953 (1999).
- 3. Y. A. Lazebnik, S. H. Kaufmann, S. Desnoyers, G. G. Poirier, W. C. Earnshaw, Nature 371, 346-347 (1994).
- 4. P. Villa, S. H. Kaufmann, W. C. Earnshaw, Trends Biochem. Sci. 22, 388-393 (1997).
- 5. Y. Liu et al., Anal.Biochem. 267, 331-335 (1999).
- 6. M. Fennell, H. Chan, A Wood. J Biomol Screen, 11, 296 (2006)
- 7. X. Wu, J. Simone, D. Hewgill, R. Siegel, PE. Lipsky, L. He. Cytometry A, 69, 477 (2006)

**Disclaimer:** This product is for research use only and is not intended for therapeutic or diagnostic applications. Please contact our technical service representative for more information.

Storage Conditions

Store at -20 °C