



## Product Data Sheet

---

**Product Name:** Caspase 3 (Apopain) Substrate 1r-z, fluorogenic  
**Catalog Number:** AS-60304-5 (5 mg)      **Lot Number:** See label on vial  
**Sequence:** (Z-Asp-Glu-Val-Asp)2Rh110 (3-letter code)  
(Z-DEVD)2Rh110 (1-letter code)  
**Molecular Weight:** 1515.5  
**% Peak Area by HPLC:** ≥95  
**Appearance:** Lyophilized yellow powder

**Peptide Reconstitution:** Use fresh concentrated Anhydrous DMSO as the solvent. Do not use DMSO diluted with aqueous solvents, as this is not an effective solvent. Do not add aqueous solvent to the peptide prior to DMSO addition as this may prevent proper solubilization. Add concentrated DMSO directly to the lyophilized peptide powder to obtain a final concentration of approximately 0.5mg/mL to 1mg/mL. Gently vortex to mix.

**Storage:** Peptide is shipped at ambient temperature. Upon receipt, store lyophilized powder at –20°C or lower. Reconstituted peptide should be aliquoted into several freezer vials and stored at –20°C or lower. Do not freeze thaw.

**Description:** Rh110 (Rhodamine 110)-derived caspase substrates are probably the most sensitive indicators widely used for the fluorimetric detection of various caspase activities. Cleavage of Rh110 peptides by caspases generates strongly fluorescent Rh110 that is monitored fluorimetrically at 510-530 nm with excitation of 488 nm, the most common excitation light source used in fluorescence instruments. Ref: Gurtu V, et al. *Anal Biochem* **251**, 98 (1997); Martins LM, et al. *J Biol Chem* **272**, 7421 (1997).

**Published Citations:**

[Sanchez-Gonzalez, PD. et al. \*Toxicol Lett\* \*\*203\*\*, 154 \(2011\).](#)

***For Research Use Only***