

Product Data Sheet

Product Name: Caspase 8 Substrate 1, chromogenic

Catalog Number: AS-25258-5 (5 mg) Lot Number: See label on vial

Sequence: Ac-IETD-pNA (1-letter code)

Molecular Weight: 638.6

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Caspase 8 peptide is freely soluble in DMSO.

Storage: Caspase 8 is shipped at ambient temperature. Upon receipt, store lyophilized peptide at –20°C or lower. Reconstituted peptide can be aliquoted and stored at –20°C or lower.

Description: pNA (4-nitroaniline)-derived caspase substrates are widely used for the colorimetric detection of various caspase activities. Cleavage of pNA peptides by caspases generates pNA that is monitored colorimetrically at ~405 nm. pNA has maximum absorption around 408 nm. Caspase-8 substrate with Km = 66 uM; Substrate for granzyme B. Grutter MG *Curr Opin Struct Biol* 10, 649 (2000); Gastman BR *Head Neck* 23, 409 (2001); Grutter MG *Curr Opin Struct Biol* 10, 649 (2000); Stennicke HR and Salvesen GS *Cell Death Differ* 6, 1054 (1999); Stennicke HR and Salvesen GS *Biochim Biophys Acta* 1387, 17 (1998); Thornberry NA and Lazebnik *Science* 281, 1312 (1998); Talanian RV, et al. *J Biol Chem* 272, (1997); Fassy F, et al. *Eur J Biochem* 253, 76 (1998); Datta R, et al. *Blood* 88, 1936 (1996); Koeplinger KA, et al. *Protein Expr Purif* 18, 378 (2000).

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

Protein concentration was measured using the DC Protein Assay Kit (Bio-Rad, Hercules, CA) and equal amounts were incubated with 200 µmol/L caspase-8 substrate Ile-Glu-Thr-Asp-pNA (AnaSpec, San Jose, CA) for 4 hours at 37°C in water bath. Release of pNA was determined by reading the absorbance at 405 nm- Shao, R. et al. *Mol Cancer Res* 3, 219 (2005).

Published Citations:

Shao, R. et al. *Mol Cancer Res* **3**, 219 (2005). Tang, C-HA. et al. *J FEBS Letters* **579**, 265 (2005).

Related Products:

Name	Cat # S	ize
Caspase 8 Substrate 1m, fluorogenic	AS-25270-5	5 mg
Caspase 8 Substrate 1, chromogenic	AS-25258-5	5 mg

For Research Use Only