

Product Information

Caspase 8 Assay Kit, Colorimetric

Product Code **CASP8C**

Storage Temperature $-20\text{ }^{\circ}\text{C}$

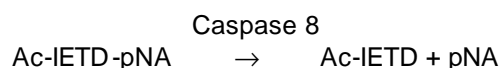
TECHNICAL BULLETIN

Product Description

Caspases (**C**ysteine-requiring **A**spartate proteases) belong to a highly conserved family of cysteine proteases with specificity for aspartic acid residues found in their substrates. Caspases play a central role in apoptosis.

Caspase 8, known also as Mch5,¹ MACH,² and FLICE,³ is localized at the top of the hierarchy of the caspase cascade and is a member of the “upstream” or initiator family of caspases. Caspase 8 exists in the cell as an inactive proenzyme of 55 kDa. It is converted to the active form, consisting of 18 and 12 kDa subunits, upon its recruitment to the cytoplasmic domain of activated death receptors such as Fas, via the adaptor protein FADD.⁴ The activation of the proenzyme is triggered by the protein’s aggregation, which leads to auto- or transprocessing. Caspase 8 activates downstream caspases (3, 6, and 7) that cleave key cellular substrates and lead to apoptotic death of the cells.^{5,6}

The assay is based on the hydrolysis of the peptide substrate Acetyl-Ile-Glu-Thr-Asp *p*-nitroaniline (Ac-IETD-pNA) by caspase 8 resulting in the release of a *p*-Nitroaniline (pNA) moiety.⁷ *p*-Nitroaniline has absorbance at 405 nm ($\epsilon^{\text{mM}} = 10.5$). The concentration of the pNA released from the substrate is calculated from the absorbance values at 405 nm or from a calibration curve performed with defined pNA solutions.



The Caspase 8 Colorimetric Assay Kit provides the reagents needed for a quick and efficient detection of caspase 8 activity in crude and purified preparations of caspase 8.

The Caspase 8 Colorimetric Assay Kit is designed for the fast and sensitive detection and measurement of caspase 8 activity in crude or purified preparations.

The assay can be performed in a 100 μl volume in a 96 well plate using a microplate fluorometer.

Reagents

The kit provides reagents sufficient for 100 tests of 100 μl in 96 well plates.

- 5' Lysis Buffer** 5 ml
 Product Code L2912
 250 mM HEPES, pH 7.4, 25 mM CHAPS,
 25 mM DTT
- 10' Assay Buffer** 5 ml
 Product Code A0344
 200 mM HEPES, pH 7.4, 1% CHAPS, 50 mM DTT,
 20 mM EDTA, 50% sucrose
- Caspase 8** 5 μg
 Product Code C6849
 Lyophilized powder. Reconstitution with 50 μl of
 17 M $\Omega\cdot\text{cm}$ water results in a solution of 100 $\mu\text{g}/\text{ml}$
 caspase 8 in 20 mM Tris-HCl, pH 8.0, 0.1%
 CHAPS, 20 mM 2-mercaptoethanol, 500 mM NaCl,
 2.5 mM EDTA, 150 mM imidazole and 10%
 sucrose.
 Specific Activity: >500 units per mg protein.
 Unit Definition: One unit is the amount of enzyme
 that will cleave 1.0 nmole of the substrate
 Ac-IETD-pNA per minute at pH 7.4 at 25 $^{\circ}\text{C}$.
- Acetyl-Ile-Glu-Thr-Asp-p-nitroaniline
Substrate** 0.1 ml
 Product Code A7332
 20 mM Ac-IETD-pNA in DMSO [MW 638.6]
- Acetyl-Ile-Glu-Thr-Asp-Aldehyde
Inhibitor** 0.1 ml
 Product Code A6464
 1 mM Ac-IETD-CHO in DMSO [MW 502.7]
- p*-Nitroaniline,** 1 ml
 Product Code N7773
 10 mM *p*-NA in DMSO [MW 138.1]
- Water (17 M $\Omega\cdot\text{cm}$)** 125 ml
 Product Code W3888

Reagents and Equipment Required but Not Provided

(Product Codes are given where appropriate)

- Apoptotic cells of interest
- BSA (optional) Product Code A8022
- ELISA reader, 96 well plate
- Polypropylene test tubes and microfuge tubes

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Preparation Instructions

Note: Use the provided 17 M Ω -cm water (Product Code W3888) in all the steps. Trace amounts of metal ions will inactivate caspase 8.

- **1 \times Assay Buffer** - 20 mM HEPES, pH 7.4, 2 mM EDTA, 0.1% CHAPS, 5 mM DTT, 5% sucrose. Dilute the 10 \times Assay Buffer (Product Code A0344) 10-fold with the supplied 17 M Ω -cm water.
- **Caspase 8 Colorimetric Substrate** (Ac-IETD-pNA) - Dilute the caspase substrate solution (Product Code A7332) 10-fold with 1 \times Assay Buffer for the microwell assay to yield 2 mM solution.
- **Caspase 8 Inhibitor** (Ac-IETD-CHO) - Dilute an aliquot of the caspase inhibitor solution (Product Code A6464) 40-fold with 1 \times Assay Buffer to yield 25 μ M.
- **Caspase 8 Positive Control.** Reconstitute the vial (5 μ g) with 50 μ l of the supplied 17 M Ω -cm water (~100 μ g/ml). Store in aliquots at -70 $^{\circ}$ C. Just before use, dilute 10-fold with 1 \times Assay Buffer to 10 μ g/ml in a polypropylene tube and use a sample of 10 μ l in the assay.
Note: Caspase 8 may adsorb to glass surfaces, therefore all dilutions should be prepared in polypropylene tubes.
- **p-Nitroaniline Standard Solution** - For the 96 well plate assay, dilute the p-Nitroaniline solution (Product Code N7773) 50-fold with 1 \times Assay Buffer to 200 μ M and build a calibration curve as shown in Table 2.

- **1 \times Lysis Buffer** - Dilute the 5 \times Lysis Buffer (Product Code L2912) 5-fold with the supplied 17 M Ω -cm water.

Note: In order to protect the cell lysate caspases from non-specific proteolysis, protease inhibitors that do not include cysteine protease inhibitors (E-64 and leupeptin) may be added.

Storage/Stability

Store the kit at -20 $^{\circ}$ C.

Procedure

When determination of caspase 8 activity in cell lysates is desired, lyse the cells with 1 \times Lysis Buffer with appropriate protease inhibitors.

The reaction schemes for assays of unknown samples and controls are shown in Table 1. Three controls are recommended for each caspase 8 colorimetric assay:

- Inhibitor-treated caspase 8 (for measuring the nonspecific hydrolysis of the substrate)
- Caspase 8 positive control.
- Reagent blank (negative control)

Equipment Required

- Flat bottomed 96 well plate for colorimetric assay
 - ELISA plate reader
1. Place 10 μ l of the diluted Caspase 8 Positive Control or x μ l of the unknown sample in a well.
 2. Add the appropriate amount of 1 \times Assay Buffer (see Table 1).
 3. Add the Caspase 8 Inhibitor to the appropriate wells. Ensure that all the cells are mixed gently to avoid bubble formation and let sit for 5 minutes at room temperature.
 4. Start the reaction with the addition of 10 μ l of the prepared Caspase 8 Colorimetric Substrate using a multichannel pipette.
 5. Place in the ELISA reader and measure the initial absorbance at 405 nm (t = 0) and then read at 5 minute intervals for t minutes (where t can be from 20-60 minutes or even longer for very dilute samples).

Table 1.Reaction Scheme for 96 Well Plate Microassay Method (Total volume = 100 μ l)

	1 st Assay Buffer	Caspase 8 10 μ g/ml	Unknown Sample	Caspase 8 Inhibitor Ac-IETD-CHO 25 μ M	Caspase 8 Substrate Ac-IETD-pNA 2 mM
Reagent blank	90 μ l	----	----	----	10 μ l
Caspase 8 Positive Control	80 μ l	10 μ l	----	----	10 μ l
Caspase 8 Positive Control + Inhibitor	78 μ l	10 μ l	----	2 μ l	10 μ l
Unknown Sample	90 -x μ l	----	x μ l	----	10 μ l
Unknown Sample + Inhibitor	88 -x μ l	----	x μ l	2 μ l	10 μ l

Table 2.*p*-Nitroaniline Calibration Curve

nmole pNA per well	pNA standard solution μ l per well	1x Assay buffer μ l per well
0	0	100
1	5	95
2	10	90
5	25	75
10	50	50
15	75	25
20	100	0

References

1. Srinivasula, S.M., *et al.*, Proc. Natl. Acad. Sci. USA, **93**, 14486-14491 (1996).
2. Boldin, M.P., *et al.*, Cell, **85**, 803-815 (1996).
3. Madema, J.P., EMBO J., **16**, 2794-2804 (1997).
4. Stennicke, H.R., *et al.*, J. Biol. Chem., **273**, 27084-27090 (1998).
5. Cohen, G.M., Biochem. J., **326**, 1-16 (1997).
6. Green, D.R., Cell, **94**, 695-698 (1998)
7. Villa P., *et al.*, Trends Biol. Sci., **22**, 388-392 (1997).

LPG,AC,MAM 10/05-1

Calculation

Calculate the caspase 8 activity as nmoles of *p*-nitroaniline released per minute per ml for the unknown sample or per mg protein for the positive control.

$$\text{Activity (nmole/min/ml)} = \frac{(A_t - A_0) \times d}{(A_{1 \text{ nmole}}) \times t \times v}$$

v = volume of sample in ml

d = dilution factor

t = reaction time in minutes

 $A_{1 \text{ nmole}}$ = absorbance of 1 nmole of pNA in the well A_t = absorbance at time t minutes A_0 = absorbance at zero time

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.