



## Recombinant Human Active Caspase-6/Mch2

### PRODUCT ANALYSIS SHEET

<b>Catalog Number:</b>	PHZ0034
<b>Lot Number:</b>	L021104
<b>Quantity:</b>	25 units/vial x 4 vials.
<b>Formulation:</b>	Lyophilized, carrier free.
<b>Reconstitution:</b>	Reconstitute the contents of the vial in 25 µL PBS per vial.
<b>Purification:</b>	Recombinant human active caspase-6 is produced in <i>E. coli</i> and purified via sequential chromatography.
<b>Sterile:</b>	Filtered prior to lyophilization through a 0.22 micron sterile filter.
<b>Specificity:</b>	Caspase-6 (also known as Mch2) is a member of the Interleukin-1 $\beta$ Converting Enzyme (ICE) family of cysteine proteases. Pro-caspase-6 is a protein of approximately 34 kDa and has two alternatively spliced isoforms. The active caspase-6 is composed of two subunits and forms a signal amplification pathway with pro-caspase-3 during apoptosis. Caspase-6 activates pro-caspase-3; caspase-3 activates pro-caspase-6. Other downstream substrates include lamin A, PARP and nuclear mitotic apparatus protein (NuMA). Caspase-6, like caspase-3, is a major active caspase in apoptosis and is involved in nuclear apoptosis. The overexpression of the full-length cDNA of caspase-6, but not the shorter splice variant, results in apoptosis.
<b>Biological Activity:</b>	The enzyme activity of this product was analyzed using caspase-6 fluorometric (VEID-AFC) and colorimetric (VEID-pNA) substrates. A unit of the active recombinant caspase-6/Mch2 is defined as the enzyme activity that cleaves 1 nmole of the caspase substrate, VEID-pNA, per hour at 37°C at saturated substrate concentration.
<b>Applications:</b>	The active caspase-6 is suitable for study of caspase-6 inhibitors, in combination with caspase-6 enzyme activity assays. It can also be used as a positive control in caspase assays or in determining the specificity of substrates.
<b>Suggested Working Dilutions:</b>	For caspase-6 activity assay, use 0.5-1 units per test for fluorometric caspase assay and 2-5 units per test is recommended for colorimetric assay. The optimal concentration should be determined for each specific application.

**This product is for research use only. Not for use in diagnostic procedures.**

[www.invitrogen.com](http://www.invitrogen.com)

Manufactured under ISO 13485 Quality Standard

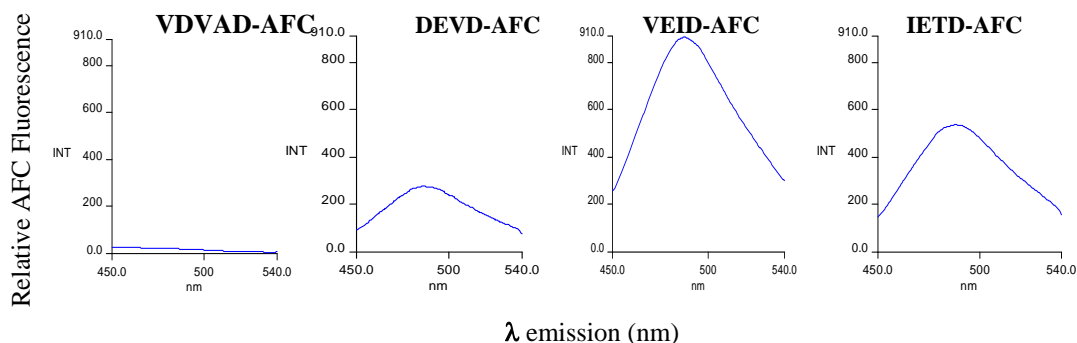
Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

PI PHZ0034

(Rev 1.0 ) (DCC-08-1232)

**Important Licensing Information** - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, [www.invitrogen.com](http://www.invitrogen.com)). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.





**Storage:** Store at  $-70^{\circ}\text{C}$ . Lyophilized powder: stable for 1 year at  $-70^{\circ}\text{C}$  and reconstituted solution: stable for 1 week at  $-70^{\circ}\text{C}$ . Keep freeze-thaw cycles to a minimum.

**References:**

Fernandes-Alnemri, T., et al. (1995) Mch2, a new member of the apoptotic Ced-3/Ice cysteine protease gene family. *Cancer Res.* 55(13):2737-42.

Alnemri, E. S., et al. (1996) Human ICE/CED-3 protease nomenclature. *Cell* 87(2):171.

Stennicke, H.R. and G.S. Salvesen, (1997) Biochemical characteristics of caspases-3, -6, -7, and -8. *J. Biol. Chem.* 272(41):25719-23.

Thornberry, N.A. and Y. Lazebnik, (1998) Caspases: enemies within. *Science* 281(5381):1312-6.

**This product is for research use only. Not for use in diagnostic procedures.**

[www.invitrogen.com](http://www.invitrogen.com)

Manufactured under ISO 13485 Quality Standard

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

PI PHZ0034

(Rev 1.0 ) (DCC-08-1232)

**Important Licensing Information** - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, [www.invitrogen.com](http://www.invitrogen.com)). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.

 invitrogen™