

Caspase-6 (Mch2) Assay Catalog #BML-SE170

Assay buffer

(50 mM HEPES, pH 7.4, 100 mM NaCl, 0.1% CHAPS, 1 mM EDTA, 10% glycerol, 10 mM DTT)

Caspase-6 (BML-SE170)

Dilute to 10 U/ μ l in assay buffer just before use.

Ac-VEID-pNA (BML-P437) colorimetric substrate

(400 μ M stock solution in Assay buffer; Store at -20°C . Warm to assay temperature before use.)

To 5 mg net peptide (MW=637.5) add 784 μ l Assay buffer (do not use DMSO), to prepare 10 mM stock. Dilute 10 mM stock to 400 μ M with Assay Buffer.

Reaction Conditions :

- 1) Add 45 μ l Assay buffer into 1/2 volume microtiter plate. Allow to equilibrate to assay temperature.
- 2) Add 5 μ l of caspase-6 (10 U/ μ l) to each appropriate well. Include 2 blanks (5 μ l assay buffer rather than caspase-6).
- 3) To start reaction, add 50 μ l Ac-VEID-pNA substrate (400 μ M in assay buffer) and mix well by pipetting up and down several times. Final substrate concentration=200 μ M.
- 4) Continuously monitor $A_{405\text{nm}}$.
- 5) Data analysis: Graph OD vs time and determine slope over the linear portion of the curve. Convert rates in OD/min to substrate/min using an extinction coefficient for p-nitroaniline of $10,500 \text{ M}^{-1} \text{ cm}^{-1}$, and adjusting for pathlength of sample (~ 0.5 cm for 100 μ l in well of a 1/2 volume, 96-well plate).

FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.