

## Product Information

|                           |  |
|---------------------------|--|
| <b>Description</b>        | Z-VEID-FMK is the specific recognition sequence for caspase-6/Mch2.  |
| <b>Target:</b>            | Apoptosis  |
| <b>Pathway:</b>           | Caspase  |
| <b>Chemical Name:</b>     | methyl (4S)-5-[[[(2S)-1-[[[(3S)-5-fluoro-1-methoxy-1,4-dioxopentan-3-yl]amino]-3-methyl-1-oxopentan-2-yl]amino]-4-[[[(2S)-3-methyl-2-(phenylmethoxycarbonylamino)butanoyl]amino]-5-oxopentanoate                             |
| <b>Molecular Formula:</b> | C <sub>31</sub> H <sub>45</sub> FN <sub>4</sub> O <sub>10</sub>  |
| <b>Supplied As:</b>       | Solid  |
| <b>Solubility:</b>        | Insoluble in H <sub>2</sub> O; ≥113.4 mg/mL in DMSO; ≥3.01 mg/mL in EtOH with gentle warming and ultrasonic For obtaining a higher solubility, please warm the tube at 37°C and shake it in the ultrasonic bath for a while. |
| <b>Stability:</b>         | At least 6 months at -20°C. Avoid freeze/thaw cycles.  |
| <b>Storage:</b>           | -20°C  |
| <b>MW:</b>                | 652.71 Da  |
| <b>Purity:</b>            | ≥98%   |

## Chemical Structure

