

Product Information

| | |
|---------------------------|---|
| Description | Boc-D-FMK is a Novel inhibitor of Caspase-3 |
| Target: | Apoptosis |
| Pathway: | Caspase |
| Chemical Name: | methyl 5-fluoro-3-[(2-methylpropan-2-yl)oxycarbonylamino]-4-oxopentanoate |
| Molecular Formula: | C ₁₁ H ₁₈ FNO ₅ |
| Supplied As: | Solid |
| Solubility: | insoluble in H ₂ O; ≥11.65 mg/mL in DMSO; ≥41.65 mg/mL in EtOH. For obtaining a higher solubility, please warm the tube at 37°C and shake it in the ultrasonic bath for a while. |
| Stability: | At least 6 months at -20°C. Avoid freeze/thaw cycles. |
| Storage: | -20°C |
| MW: | 263.26 Da |
| Cas No. | 187389-53-3, 634911-80-1 |
| Purity: | ≥98% by HPLC |

Chemical Structure

