Glowing products for science

Ac-DEVD-CHO Caspase-3 Inhibitor

Ac-DEVD-CHO is a synthetic tetrapeptide competitive inhibitor for Caspase-3/7, and contains the amino acid sequence of the PARP cleavage site. Ac-DEVD-CHO can be used to inhibit Caspase-3/7 activity in apoptotic cells, and to study events downstream of Caspase-3/7 activation.



Product attributes

Call us: 800-304-5357

CAS number

184179-08-6

Email: btinfo@biotium.com

Product Description

Ac-DEVD-CHO is a synthetic tetrapeptide competitive inhibitor for Caspase-3/7, and contains the amino acid sequence of the PARP cleavage site. Ac-DEVD-CHO can be used to inhibit Caspase-3/7 activity in apoptotic cells, and to study events downstream of Caspase-3/7 activation.

- Soluble in DMSO
- Store at 4°C
- C₂₀H₃₀N₄O₁₁
- MW: 502.5

This datasheet was generated on November 7, 2025 at 04:44:29 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/ac-devd-cho-caspase-3-inhibitor/