



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

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

CAS number	184179-08-6
------------	-------------

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
- $C_{20}H_{30}N_4O_{11}$
- MW: 502.5

This datasheet was generated on June 7, 2026 at 03:26:51 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/ac-devd-cho-caspase-3-inhibitor/>