

## Caspase-3/7 Inhibitor Ac-DEVD-CHO

Table 1. Contents and Storage Information

Catalog No.	Material	Amount	Storage	Stability
A043A	Ac-DEVD-CHO	1 mg	-20 °C	The product is stable for 1 year when stored as directed.
A043B	Ac-DEVD-CHO	5 mg		

### Product Description

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

### Specifications:

- Substrate name: Ac-DEVD-CHO
- Molecular weight: 502.5
- Peptide sequence: DEVD
- Soluble in DMSO

### Application

Enzyme assay.