

Ac-DEVD-CHO *CAS 169332-60-9*

Catalog number: 13403

Unit size: 1 mg

Component	Storage	Amount
Ac-DEVD-CHO *CAS 169332-60-9*	Freeze (<-15 °C), Minimize light exposure	1 mg

OVERVIEW

Ac-DEVD-CHO is a reversible inhibitor of IL-1 β converting enzyme (ICE); inhibits poly(ADP-ribose) polymerase cleavage by apopain (caspase 3).

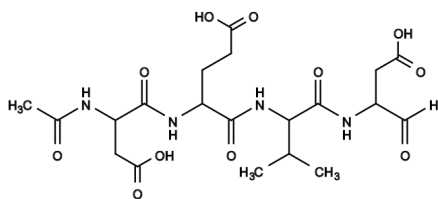
EXAMPLE DATA ANALYSIS AND FIGURES


Figure 1. Chemical structure for Ac-DEVD-CHO *CAS 169332-60-9*

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.