

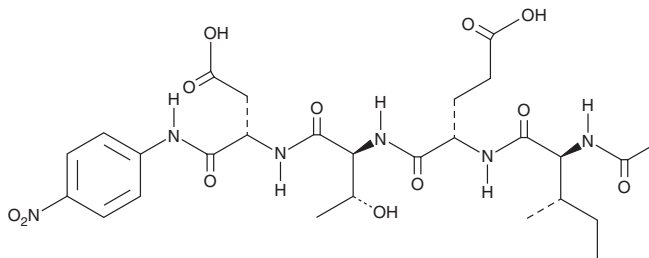
# PRODUCT INFORMATION



## Ac-IETD-pNA

Item No. 24568

**CAS Registry No.:** 219138-21-3  
**Formal Name:** N-acetyl-L-isoleucyl-L- $\alpha$ -glutamyl-L-threonyl-N-(4-nitrophenyl)-L- $\alpha$ -asparagine  
**Synonyms:** Ac-Ile-Glu-Thr-Asp-pNA, Caspase-8 Chromogenic Substrate I  
**MF:** C<sub>27</sub>H<sub>38</sub>N<sub>6</sub>O<sub>12</sub>  
**FW:** 638.6  
**Purity:**  $\geq$ 95%  
**Supplied as:** A lyophilized powder  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Ac-IETD-pNA is supplied as a lyophilized powder. A stock solution may be made by dissolving the Ac-IETD-pNA in the solvent of choice. Ac-IETD-pNA is soluble in the organic solvent formic acid, which should be purged with an inert gas, at a concentration of approximately 1 mg/ml.

### Description

Ac-IETD-pNA is a substrate for caspase-8.<sup>1</sup> Caspase-8 preferentially binds to and cleaves the Ile-Glu-Thr-Asp (IETD) peptide sequence to release *p*-nitroaldehyde, which can be quantified by colorimetric detection at 405 nm as a measure of enzyme activity.

### Reference

1. Smith, G.K., Barrett, D.G., Blackburn, K.J., *et al.* Expression, preparation, and high-throughput screening of caspase-8: Discovery of redox-based and steroid diacid inhibition. *Arch. Biochem. Biophys.* **399**(2), 195-205 (2002).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/01/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM