

# PRODUCT INFORMATION



## Suc-YVAD-pNA

Item No. 28135

CAS Registry No.: 208264-84-0

Formal Name: N-(3-carboxy-1-oxopropyl)-L-tyrosyl-L-valyl-L-alanyl-N-(4-nitrophenyl)-L- $\alpha$ -asparagine

Synonyms: Caspase-1 and Caspase-4 Chromogenic Substrate, Suc-Tyr-Val-Ala-Asp-p-nitroanilide

MF:  $C_{31}H_{38}N_6O_{12}$

FW: 686.7

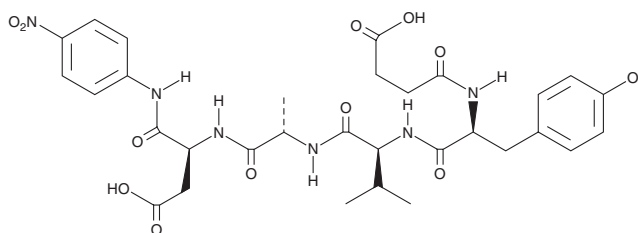
Purity:  $\geq 95\%$

UV/Vis.:  $\lambda_{max}$ : 223, 317 nm

Supplied as: A solid

Storage:  $-20^{\circ}C$

Stability:  $\geq 4$  years



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

### Laboratory Procedures

Suc-YVAD-pNA is supplied as a solid. A stock solution may be made by dissolving the Suc-YVAD-pNA in the solvent of choice, which should be purged with an inert gas. Suc-YVAD-pNA is soluble in the organic solvent acetic acid (50%).

### Description

Suc-YVAD-pNA is a colorimetric substrate for caspase-1/interleukin-1 $\beta$ -converting enzyme (ICE) and caspase-4.<sup>1</sup> Caspase-1 and caspase-4 preferentially bind to and cleave the Tyr-Val-Ala-Asp (YVAD) peptide sequence to release p-nitroanilide (pNA), which can be quantified by colorimetric detection at 405 nm as a measure of caspase-1 and caspase-4 activity.

### Reference

1. Margolin, N., Raybuck, S.A., Wilson, K.P., et al. Substrate and inhibitor specificity of interleukin-1 $\beta$ -converting enzyme and related caspases. *J. Biol. Chem.* **272**(11), 7223-7228 (1997).

#### 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, 10/04/2022

#### 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