

# Caspase-14 (human), (recombinant)

Caspase-14 is considered a member of the caspase-1 subfamily and is the only subfamily member with a short (or absent) N-terminal prodomain. In mouse, caspase-14 is highly expressed in embryos, and in adult skin. Within the skin, caspase-14 is present only in the suprabasal layers of the epidermis. A deficiency in caspase-14 expression and processing is associated with psoriasis lesions, raising the possibility that low caspase-14 levels may be among the factors that cause the disease. Elevated caspase-14 expression has been reported in ductal carcinoma *in situ* and invasive breast cancers. Unlike other procaspases, which are processed by cleavage after aspartate residues, procaspase-14 processing occurs between Ile<sup>152</sup> and Lys<sup>153</sup>. In the presence of the kosmotropic salt, sodium citrate, both recombinant caspase-14, cut *in vitro* after Asp<sup>146</sup> by granzyme B9, and the endogenous form extracted from human epidermal stratum corneum, are activated to cleave tetrapeptide substrates incorporating the sequence WEHD.

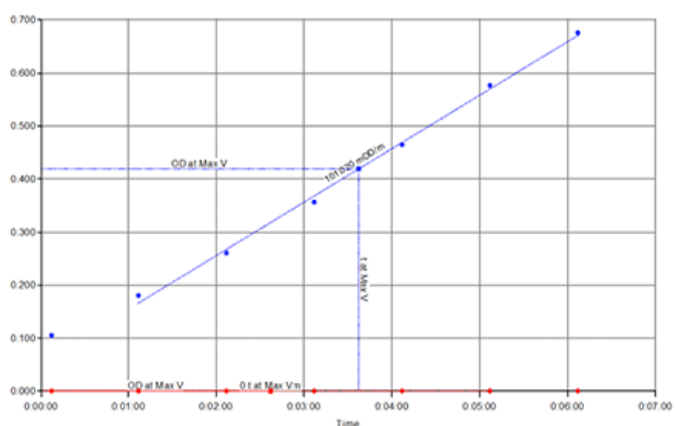
## Ordering Information

[Order Online »](#)

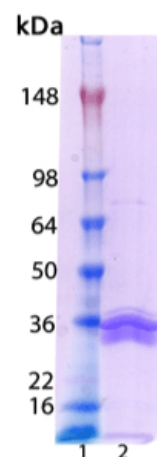
ENZ-PRT206-0020	20µg
-----------------	------

## Manuals, SDS & CofA

[View Online »](#)



Activity assay of Caspase-14 (Prod. No. ENZ-PRT206): Enzyme incubated with Ac-WEHD-pNA (Prod. No. ALX-260-082), 200  $\mu$ M, in Caspase-14 assay buffer (1.1 M sodium citrate, 100 mM HEPES/NaOH, pH 7.0, 60 mM NaCl, 0.01% CHAPS and 5 mM DTT) at 37C for 6 minutes, reading with 1 minute intervals. Blue: 2  $\mu$ L of enzyme; Red: Blank- no enzyme.



SDS-PAGE analysis: Lane 1: MW Marker, Lane 2: 2  $\mu$ g Caspase-14 (Prod. No. ENZ-PRT206).

## Handling & Storage

Handling	Centrifuge vial before opening. For prolonged storage, dilute to working aliquots and store at -80°C. Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application Notes	Useful tool to study the enzyme regulation and kinetics, cleave target substrates, screen for inhibitors.
Formulation	Lyophilized from sterile PBS, pH 7.4, 5% trehalose, 5% mannitol.
MW	~ 30kDa
Reconstitution	Reconstitute in sterile water.
Source	Produced in <i>E. coli</i> . Human caspase-14 (2-242 aa) fused to a polyhistidine tag at the N-terminus.
UniProt ID	P31944



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)