## DPPIII (human), (recombinant)

DPPIII is a metalloproteinase that cleaves biologically active peptides such as enkephalins, angiotensins, and endomorphins, and has been implicated in cancer, pain, and inflammation.

## **Ordering Information**

Order Online »

BML-SE529-0010

10µg

Manuals, SDS & CofA

**View Online** »

## **Handling & Storage**

**Long Term Storage** -80°C

**Shipping** Dry Ice

Regulatory Status RUO - Research Use Only

**Product Details** 

**Alternative Name** Dipeptidyl Peptidase III, DPP3

**Application Notes** Useful tool to study enzyme kinetics, cleave target

substrates, and screen for inhibitors.

MW 109 kDa

Source Produced in insect cells. Active recombinant human DPPIII

with an N-terminal purification tag. Produced in a

baculovirus expression system.

**UniProt ID Q9NY33** 

Last modified: May 29, 2024



info-

eu@enzolifesciences.com