POP/PREP (human), (recombinant)

POP/PREP is a cytoplasmic, membrane-associated, or serum serine protease that cleaves peptide hormones and neuropeptides such as such as substance P, oxytocin, neurotensin, angiotensins, and bradykinin. This enzyme is related to DPPIV and is useful for specificity screening of DPPIV inhibitors.

Ordering Information

Order Online »

BML-SE545-0010

10µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Prolyl oligopeptidase, Prolyl endopeptidase, Post-proline

cleaving enzyme

Application NotesUseful tool to study enzyme kinetics, cleave target

substrates, and screen for inhibitors.

MW 81.6 kDa

Source Produced in E. coli.

UniProt ID P48147

Last modified: May 29, 2024



eu@enzolifesciences.com