

PRODUCT DATA SHEET



DPPII/DPP7 (human, recombinant)
(EC 3.4.14.2)

CATALOG NO.: SE-564

LOT NO.: temp

DESCRIPTION: MW=52.5 kDa by SDS-PAGE. Active recombinant human DPPII/DPP7/QPP (dipeptidyl peptidase II, quiescent cell proline dipeptidase) (GenBank Accession # NM_013379, amino acids 26-492) with a C-terminal purification tag. DPPII is a circulating homodimeric serine protease that cleaves short bioactive peptides such as substance P, glucagon, and bradykinin¹⁻³. Its physiological function is unknown, but it is implicated in cell differentiation and the degradation of short fragments of collagen. This enzyme is related to DPPIV and is useful for specificity screening of DPPIV inhibitors⁴.

PURITY: >80% by SDS-PAGE.

SPECIFIC ACTIVITY: 1882 pmol/min/μg using 20 μM AP-AMC in 10 mM Tris-HCl, pH 7.4, 10 mM MgCl₂, 1 mM MnCl₂ at room temperature.

USAGE: Study enzyme kinetics, cleave target substrates, and screen compounds^{1,5}.

SUPPLIED AS: 10 μg at 0.6μg/μl in 25 mM Tris-Hcl pH 8.0, 130 mM NaCl, 0.05% Tween-20, 10% glycerol.

STORAGE: -70°C. The enzyme is stable on ice for at least several hours. However, it is recommended that thawing and dilution of the enzyme be done within as short a time as possible before start of the assay. After initial defrost, aliquot product into individual tubes and refreeze at -70°C. Avoid repeated freeze/defrost cycles.

NOTE: When stored under the above conditions, this enzyme is stable at the concentration supplied, in its current storage buffer. Procedures such as dilution of the enzyme followed by refreezing could lead to loss of activity.

REFERENCES:

1. M.-B. Maes *et al.* *Biochem J.* 2005 **386** 315
2. M.-B. Maes *et al.* *Clin. Chem. Acta* 2007 **380** 31
3. J.S. Rosenblum and J.W. Kozarich *Curr. Opin. Chem. Biol.* 2003 **7** 496
4. B. Lee *et al.* *Eur. J. Pharmacol.* 2008 **589** 306
5. P. Van der Veken *et al.* *J. Med. Chem.* 2007 **50** 5568

FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.