

PRODUCT DATA SHEET



FAP (human, recombinant)
(EC 3.4.21.B28)

CATALOG NO.: SE-409

LOT NO.: temp

DESCRIPTION: MW=85.5 kDa by SDS-PAGE. Active recombinant human FAP (fibroblast activation protein α ; seprase, surface expressed protease; APCE, antiplasmin cleaving enzyme)¹ (GenBank Accession # NM_004460), residues Ser29-Asp760, representing a naturally-occurring cleaved (soluble) form, with a C-terminal purification tag. FAP is a homodimeric transmembrane serine prolyl peptidase that digests gelatin³, collagen, and α_2 -antiplasmin¹⁻⁴. Like DPPIV, it can be expressed as a soluble form³. Due to its implication in cancer and blood disorders, it is a target for drug discovery³⁻⁵. In addition, because this enzyme is related to DPPIV, it is useful for specificity screening of DPPIV inhibitors⁶.

PURITY: >90% by SDS-PAGE.

SPECIFIC ACTIVITY: 10.2 pmol/min/ μ g, using 5 μ M Ala-Pro-AMC in 50 mM Tris-HCl, pH 7.5, 10 mM MgCl₂, 1 mM MnCl₂, at room temperature.

USAGE: Study enzyme kinetics, cleave target substrates, and screen compounds.

SUPPLIED AS: 10 μ g at 0.6 μ g/ μ l in 25 mM Tris-Hcl pH 8.0, 100 mM NaCl, 0.05% Tween-20, 3 mM DTT, 50% 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. P. O'Brien and B.F. O'Connor *Biochim. Biophys. Acta* 2008 **1784** 1130
2. J.E. Park *et al. J. Biol. Chem.* 1999 **274** 36505
3. J.S. Rosenblum *et al. Curr. Opin. Chem. Biol.* 2003 **7** 496
4. K.N. Lee *et al. Blood.* 2006 **107** 1397
5. C.Y. Edosada *et al. J. Biol. Chem.* 2006 **281** 7437
6. P. van der Veken *et al. Curr. Top. Med. Chem.* 2007 **7** 621

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