

FAP (human), (recombinant) (His- tag)

FAP is a homodimeric serine prolyl peptidase that digests gelatin, collagen, and α 2-antiplasmin. Due to its implication in cancer and blood disorders, it is a target for drug discovery. In addition, because this enzyme is related to DPPIV, it is useful for specificity screening of DPPIV inhibitors.

Ordering Information

[Order Online »](#)

BML-SE409-0010	10 μ g
----------------	------------

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	Fibroblast activation protein α , Seprase, Surface expressed protease, APCE, Antiplasmin cleaving enzyme
Application Notes	Useful tool to study enzyme kinetics, cleave target substrates, and screen for inhibitors.
Formulation	Liquid. In 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 80 mM Imidazole, 3 mM DTT, and 20% glycerol.
MW	86 kDa
Source	Produced in a baculovirus expression system. Active recombinant human FAP (aa 29-760), representing a naturally-occurring cleaved (soluble) form fused at the C-terminus to a His-tag.
UniProt ID	Q12884

Last modified: July 3, 2024



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)