

Data Sheet

FAP, His-Tag

Human, Recombinant, C-terminal His-tag

Catalog #: 80100

Lot #: 190607

Conc.: 0.63 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200mM Imidazole, 3 mM DTT and 20% glycerol.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

References:

1. Edosada, C.Y. *et al.*, *J. Biol. Chem.* **281** (11), 7437-7444 (2006).
2. Edosada, C.Y. *et al.*, *FEBS Lett.* **580** (6), 1581-1586 (2006).

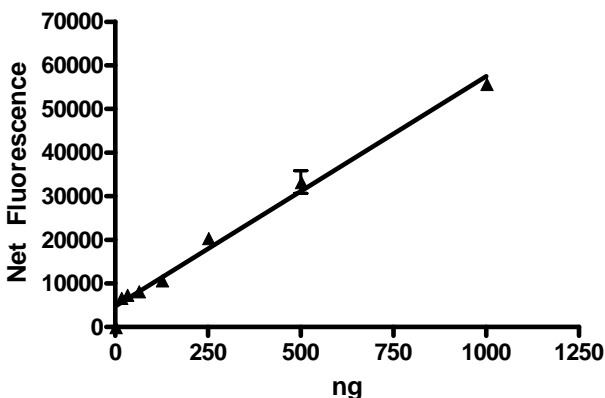
Description: Human FAP, also known as Fibroblast Activation Protein, Integral Membrane Serine Protease, and Surface-Expressed Protease, GenBank Accession No. NM_004460, a.a. 29-760 (end) with C-terminal His-tag, expressed in a Baculovirus infected Sf9 cell expression system. MW = 86 kDa.

Specific Activity: ≥ 83 pmol/min/ μg
Assay Conditions: Assay was done according to BPS Fluorogenic FAP Assay Kit #80210 with FAP enzyme titrated from 1000 ng/rxn – 0 ng/rxn.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

Specific Activity

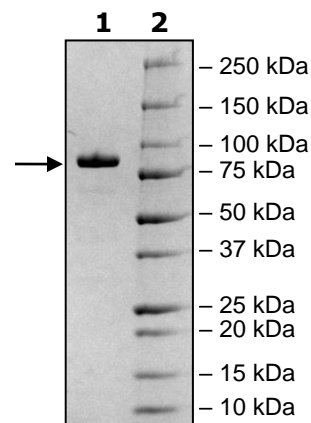


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2 μg FAP
Lane 2:
Protein Marker

MW: 86 kDa

Purity: $\geq 90\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com