

Data Sheet

FAP, His-Tag (Mouse)

Mouse, Recombinant, C-terminal His-tag

Catalog #: 100226

Lot #: 181023

Conc.: 1.25 mg/ml

Description: Mouse FAP, GenBank Accession No. NM_007986, amino acids 26-762 (end), with C-terminal His-tag, MW=88 kDa, expressed in HEK293 expression system.

Formulated in: 8mM phosphate, 110mM NaCl, 2.2mM KCl, pH7.4, and 20% glycerol

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

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).

Quality Assurance

4-20% SDS-PAGE

Coomassie Staining

Lane 1:

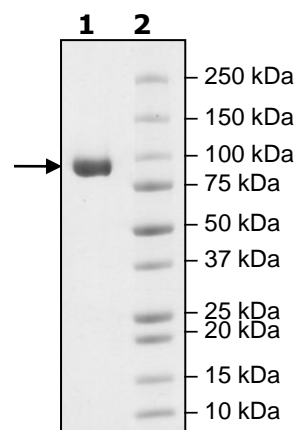
4 μg FAP

Lane 2:

Protein Marker

MW: 88 kDa

Purity: 90%



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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com