

Data Sheet

PrCP

Human, recombinant Lysosomal carboxypeptidase,
C-terminal His tag

Cat. #: 80380

Lot #: 110517-1

Conc.: 0.06 mg/ml

Formulated in: 40 mM Tris pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% Glycerol, 3 mM DTT, 200 mM Imidazole

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

References:

1. Tan, F., *et al.* 1993. *J Biol Chem* **268** (22): 16631–8.
2. Yang, H.Y., *et al.* 1968. *Nature* **218** (5148): 1224–6.
3. Wang, L., *et al.* 2006. *Am. J. Obstet. Gynecol.* **195** (1): 162–71.

Description:

Human recombinant Lysosomal Pro-X carboxypeptidase, also known as angiotensinase C (GenBank Accession #NM_005040), (31-496) with C-terminal His tag, MW=53 kDa, expressed in HEK293 cells.

Application:

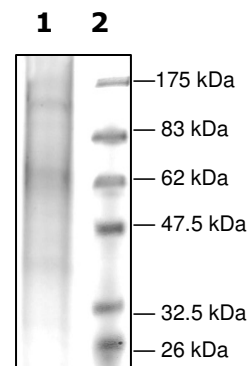
Useful for the cleavage of C-terminal amino acids linked to proline in peptides such as angiotension II, III and des-Arg9-bradykinin.

Quality Assurance

10% SDS-PAGE Coomassie staining

Lane 1:
1 μg PrCP
Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 53 kDa.
Purity: $\geq 15\%$



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



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Amino acid sequence:

NP_005031.1

SLHLPTNPTSLPAVAKNYSVLYFQQKVDHFGFNTVKTFNQRYLVADKYWKKNGGSILFYTGNEGDI
IWFCNNTGFMWDVAEELKAMLVFAEHRYYGESLPFGDNSFKDSRHLNFLTSEQALADFAELIKHLK
RTIPGAENQPVIAIGGSYGGMLAAWFRMKYPHMVVGALAASAPIWQFEDLVPCGVFMKIVTTDFRK
SGPHCSESIHRSDAINRLSNTGSGLQWLTGALHLCSP L TSQDIQHLKDWISETWVNLAMVDYPY
ASNFLQPLPAWPIKVVVCQYLKNPNVSDSLLLQNIQALNVYYNYSGQVKCLNISETATSSLGTLGW
SYQACTEVVMPFCTNGVDDMFEPHSWNLKELSDDCFQQWGVRPRPSWITMYGGKNISSHTNIV
FSNGELDPWGGGVTKDITDLVAVTISEGAHHLDLRTKNALDPMSVLLARSLEVRHMKNWIRDFY
DSAGKQH**HHHHHH**

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