

# Neprilysin (human), (recombinant) (His- tag)

## Highly active

Neprilysin is a glycosylated zinc endopeptidase whose physiological targets include bradykinin, glucagon, GLP-1, and neuropeptides. Preferential cleavage of polypeptides between hydrophobic residues, particularly with Phe or Tyr at P1.

Neprilysin is an integral plasma membrane (or secreted soluble) protease that has a broad expression pattern, but is most abundant in the kidney. Due to its involvement in cancer, Alzheimer's and Parkinson's diseases, hypertension, diabetes, and pain, it is an important research target.

Citations: 3

[View Online »](#)

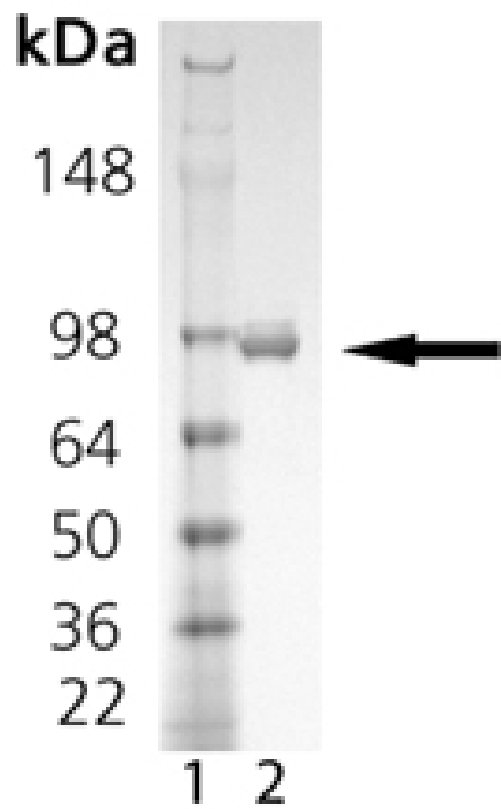
## Ordering Information

[Order Online »](#)

BML-SE532-0010	10µg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane1: MWM; Lane 2: 1.0  $\mu$ g of purified Nephilysin protein.

# Handling & Storage

<b>Use/Stability</b>	Stable on ice for at least 1h. Do not freeze in dilute form.
<b>Handling</b>	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	NEP, Neutral endopeptidase 24.11, Enkephalinase, CALLA, Common acute lymphoblastic leukemia antigen, CD10
<b>Application Notes</b>	Useful tool to study enzyme kinetics, cleave target substrates, and screen for inhibitors.
<b>Formulation</b>	Liquid. In 50mM TRIS, pH 8.0, containing 300mM sodium chloride and 20% glycerol.
<b>MW</b>	80.7 kDa (calculated) / ~85 kDa (glycosylated form on SDS-PAGE)
<b>Purity</b>	≥95% (SDS-PAGE)
<b>Purity Detail</b>	Purified as the glycosylated soluble form.
<b>Sequence</b>	NEP from human cDNA, transcript variant 1, aa 53-750 (D <sub>53</sub> – W <sup>750</sup> ; Identical to GeneBank accession NM_000902) is fused at the N-terminus to a His-tag.
<b>Source</b>	Produced in Sf9 insect cells. Produced in a baculovirus expression system.
<b>UniProt ID</b>	P08473