

BACE2 (soluble) (human), (recombinant) (His- tag)

Highly active BACE2 is an aspartic protease found in both transmembrane and soluble forms. It is related to, and possesses an overlapping substrate profile with, BACE1. It is likely involved in Alzheimer's disease, and implicated in other neurological, pancreatic, and muscle pathologies.

Citations: 3

[View Online »](#)

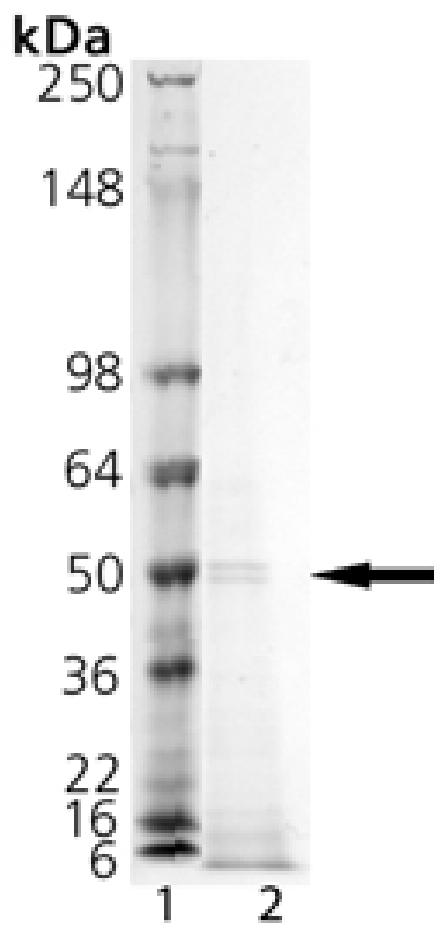
Ordering Information

[Order Online »](#)

BML-SE550-0010	10µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane 1: MW Marker, Lane 2: 1 μ g BACE2

Handling & Storage

Use/Stability	BACE2 is stable after 5 freeze-thaws at 0.3-0.4 µg/µl; freeze-thaw stability of more dilute preparations has not been tested and could lead to loss of activity.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	β-Secretase 2, β-Site APP-cleaving enzyme 2, Memapsin-1, ASP-1
Application Notes	Useful tool to study enzyme kinetics, cleave target substrates, and screen for BACE2 inhibitors or BACE1 inhibitor selectivity.
Formulation	Liquid. In 50mM TRIS-HCl, pH 7.5, containing 100mM sodium chloride and 20% glycerol.
MW	~45kDa (calculated); ~48kDa (doublet by SDS-PAGE)
Purity	≥80% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Sequence	aa Ala ²¹ -Pro ⁴⁶⁶ .
Source	Produced in insect cells. Recombinant soluble BACE2, cloned from human cDNA, secreted as zymogen from insect cells, purified using a C-terminal His-tag, and cleaved to the mature <i>active</i> form. Produced in a baculovirus expression system.
Specific Activity	≥20 U/µg enzyme. One unit will hydrolyze one pmole Mca-KPLGL-Dpa-AR substrate (Prod. No. BML-P276) (10µM) per minute at 37°C, in 50mM NaOAc, pH 4.5, 1M sodium chloride.

Last modified: May 29, 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)