

Tyrosine hydroxylase polyclonal antibody

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SA497-0100	100µl
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

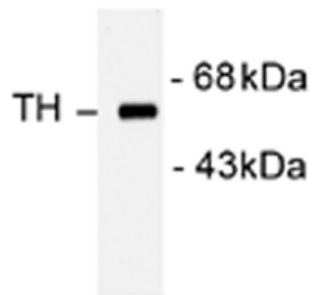


Figure 1. Western Blot analysis of rat caudate lysate using Tyrosine hydroxylase, pAb (Prod.no. BML-SA497)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name TH

Application ICC, IHC, WB

Formulation Liquid. In 10mM HEPES, pH 7.5, containing 150mM NaCl, 10% BSA and 50% glycerol.

Host Rabbit

Immunogen Denatured tyrosine hydroxylase (TH) from rat pheochromocytoma.

Purity Detail Affinity purified.

Recommendation Immunohistochemistry (1:1,000)Immunofluorescence (1:1,000)Western blot
Dilutions/Conditions (1:1,000)Optimal conditions must be determined for each application individually.

Species Reactivity Mammalian

UniProt ID P04177

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



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)