## 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 »

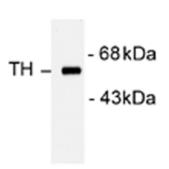


Figure 1. Western Blot analysis of rat caudate lysate using Tyrosine hydroxylase, pAb (Prod.no. BML-SA497)

## **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 Dilutions/Conditions** Immunohistochemistry (1:1,000)Immunofluorescence

(1:1,000)Western blot (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

.

Last modified: May 29, 2024

