

---

## MAOB Polyclonal Antibody, HRP Conjugated

(Catalog #A53197)

---

### Background

Catalyzes the oxidative deamination of biogenic and xenobiotic amines and has important functions in the metabolism of neuroactive and vasoactive amines in the central nervous system and peripheral tissues. MAOB preferentially degrades benzylamine and phenylethylamine.

### Description

MAOB Polyclonal Antibody, HRP Conjugated. HRP. Raised in: Rabbit.

### Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

### Specificity

Human

### Isotype

IgG

### Uniprot ID

P27338

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Amine oxidase [flavin-containing] B protein (2-489AA)

### Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Monoamine oxidase type B, MAOB

### Application

ELISA; Recommended dilution: ELISA 1:500-1:1000

---