

Recombinant MAOB protein

Catalog No: 31503, 31907

Lot No: 25315001

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.3 µg/µl

Source: Human

Buffer Contents: Full length recombinant MAOB protein was expressed in Sf9 cells and is supplied in 25 mM Hepes pH 7.5, 300 mM NaCl, 5% Glycerol, 0.04% Triton X-100, 0.2 mM TCEP, 0.2 mg/ml 3× Flag peptide.

Background: MAOB (Monoamine Oxidase B) is one of two neighboring mitochondrial enzyme family members which catalyze the oxidative deamination of amines, such as dopamine, norepinephrine, and serotonin. The other member is MAOA. MAOB preferentially degrades benzylamine and phenylethylamine. MAOB catalyzes the oxidative deamination of biogenic and xenobiotic amines and plays an important role in the metabolism of neuroactive and vasoactive amines in the central nervous system and peripheral tissues.

Protein Details: Recombinant MAOB (accession number NP_000889.3) was expressed in Sf9 cells and contains an N-terminal FLAG-Tag with an observed molecular weight of 59.8 kDa. The recombinant protein is >85% pure by SDS-PAGE.

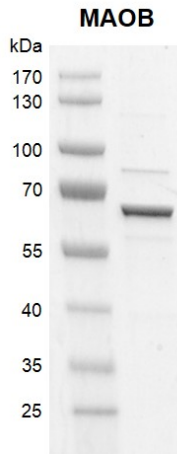
Application Notes: Recombinant MAOB protein is suitable for use in the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

Monoamine Oxidase Assay:

40 µM substrate was incubated with MAOA protein in reaction buffer for 1 hour at room temperature, then the product was incubated with Detection Reagent for 0.5 hour at temperature. The activity of MAOA was detected by Promega MAO-Glo Assay.

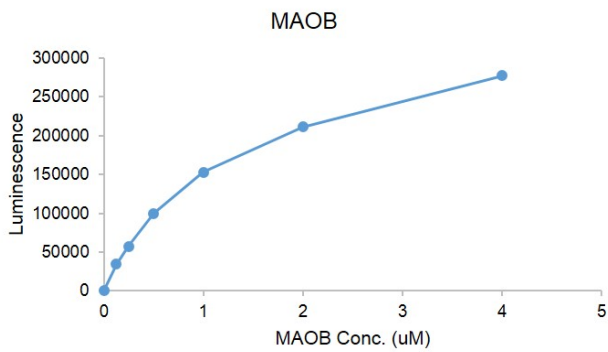
Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



Recombinant MAOB protein gel.

MAOB protein was run on an 8% SDS-PAGE gel and stained with Coomassie blue.



MAO-Glo™ Assay for MAOA activity

40 μ M substrate was incubated with MAOA protein in reaction buffer for 1 hour at room temperature, then the product was incubated with Detection Reagent for 0.5 hour at temperature. The activity of MAOA was detected by Promega MAO-Glo Assay.