

Recombinant MAOA protein

Catalog No: 31502, 31906

Lot No: 25315001

Expressed In: Baculovirus

Quantity: 10, 1000 µg

Concentration: 0.35 µg/µl

Source: Human

Buffer Contents: Full length recombinant MAOA 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: MAOA (Monoamine Oxidase A) 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 MAOB. Mutation of MAOA results in Brunner syndrome and maoa-related behavior disorders. MAOA has important functions in the metabolism of neuroactive and vasoactive amines in the central nervous system and peripheral tissues. MAOA preferentially oxidizes biogenic amines such as 5-hydroxytryptamine (5-HT), norepinephrine and epinephrine.

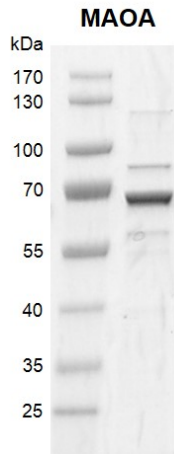
Protein Details: Recombinant MAOA (accession number NP_000231.1) was expressed in Sf9 cells and contains an N-terminal FLAG-Tag with an observed molecular weight of 60.7 kDa. The recombinant protein is >80% pure by SDS-PAGE.

Application Notes: Recombinant MAOA 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.

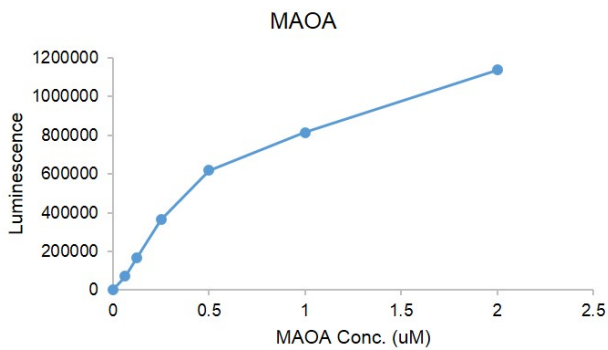
Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

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



Recombinant MAOA protein gel.

MAOA 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.