

Recombinant COQ3 protein

Catalog No: 81457, 81557

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Source: Human

Buffer Contents: Recombinant COQ3 protein is supplied in 25 mM Tris-HCl pH 8.0, 300 mM NaCl, 10% glycerol and 0.5 mM TCEP.

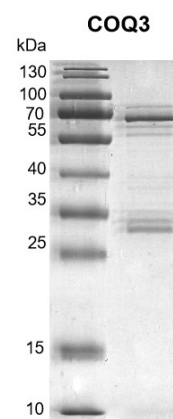
Background: COQ3(Ubiquinone biosynthesis O-methyltransferase, mitochondrial) is an O-methyltransferase that catalyzes the 2 O-methylation steps in the ubiquinone biosynthetic pathway.

Protein Details: Recombinant COQ3 protein that includes full length of human COQ3 protein (accession number NP_059117.3) was expressed was expressed in *E. coli* cells with an N-terminal GST tag. The molecular weight of the protein is 67.89 kDa.

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for a given product is shown on the lot-specific Technical Data Sheet..

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 COQ3 protein

12.5% SDS-PAGE Coomassie staining

MW: 67.89 kDa

Purity: ≥ 85%