

## Recombinant MAP2K2 protein

**Catalog No:** 81332, 81632

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.3 µg/µl

**Source:** Human

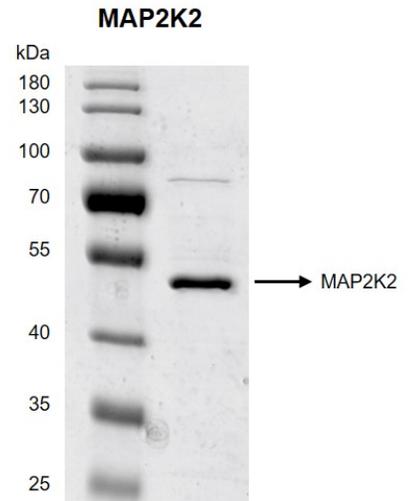
**Buffer Contents:** Recombinant MAP2K2 protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, and 0.5 mM TCEP.

**Background:** MAP2K2 (Mitogen-Activated Protein Kinase Kinase 2) is a dual specificity protein kinase that belongs to the MAP kinase kinase family. This kinase plays a critical role in mitogen growth factor signal transduction. It phosphorylates and activates MAPK1/ERK2 and MAPK2/ERK3. The activation of this kinase itself is dependent on the Ser/Thr phosphorylation by MAP kinase kinase kinases. Mutations in this gene cause cardiofaciocutaneous syndrome (CFC syndrome), a disease characterized by heart defects, cognitive disability, and distinctive facial features similar to those found in Noonan syndrome.

**Protein Details:** Recombinant MAP2K2 protein was expressed in baculovirus expression system as the full length protein (accession number NP\_109587.1) with a N-terminal FLAG Tag. The molecular weight of the protein is 45.7 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 for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.



**Recombinant MAP2K2 protein gel**  
10% SDS-PAGE with Coomassie blue staining  
MW: 45.7 kDa  
Purity: >90%