

Calmodulin

human, recombinant, *E. coli*

Cat. No.	Amount
PR-390	100 μ g

For *in vitro* use only
Quality guaranteed for 12 months
Store at -80°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied as 0.5 mg/ml solution in 20 mM HEPES-KOH pH 7.0, 50 mM NaCl, 1 mM EDTA and 1 mM DTT.

Activity

Assayed for calcium binding and activation of MLCK.

Purity

>95% by SDS-PAGE.

Description

Calmodulin (CaM) is an intracellular receptor protein for Ca²⁺-ions. It activates the myosin light chain kinase which is the catalyst of myosin phosphorylation.

CaM participates in the activation of enzymes such as cyclic nucleotide-dependent phosphodiesterase, calcineurin, ATPase, Myosin Light Chain Kinases, and CAM kinase.

Selected References:

- Cheung, W.Y. (1980) Calmodulin plays a pivotal role in cellular regulation. *Science* **207**:19.
Vetter *et al.* (2003) Novel aspects of calmodulin target recognition and activation. *Eur. J. Biochem.* **270**:404.
Agell *et al.* (2002) Modulation of the Ras/Raf/MEK/ERK pathway by Ca²⁺ and calmodulin. *Cellular Signalling* **14**:649