

Ventricular Myosin Regulatory Light Chain LCII human, recombinant, *E. coli*

Cat. No.	Amount
PR-393	100 µg

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, 1 mM DTT.

Activity

Assayed for phosphorylation by MLCK.

Purity

>95% by SDS-PAGE.

Description

LCII (ventricular myosin regulatory light chain) is a subunit of the heterohexameric myosin protein, which modulates striated muscle contraction in response to phosphorylation by MLCK (myosin light chain kinase).

Selected References:

- Somlyo *et al.* (2000) Signal transduction by G-proteins, Rhokinase and protein phosphatase to smooth muscle and nonmuscle myosin II. *Journal of Physiology* **522**:177.
- Davis *et al.* (2001) The overall pattern of cardiac contraction depends on a spatial gradient of myosin regulatory light chain phosphorylation. *Cell* **107**:631.
- Davis *et al.* (2002) Kinetic effects of myosin regulatory light chain phosphorylation on skeletal muscle contraction. *Biophysical Journal* **83**:359.