

Titin, I-band fragment human, recombinant, *E. coli*

Cat. No.	Amount
PR-392	100 μ g

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 20 mM HEPES-KOH pH 7.0,
50 mM NaCl, 1 mM EDTA and 1 mM DTT.

Purity

>95% by SDS-PAGE

Description

Titin is the largest known protein. Next to actin and myosin, titin is the third most abundant protein in the striated muscle sarcomere, providing passive tension to muscle and acting as a template for sarcomere formation. Titin I27-I34 contains 6 Immunoglobulin-like domains from the I-band region and can be used as a force and length standard for atomic force spectroscopy.

Selected References:

- Rief *et al.* (1997) Reversible unfolding of individual titin immunoglobulin domains by AFM. *Science* **276**:1109.
Clark *et al.* (2002) Striated muscle cytoarchitecture: An intricate web of form and function. *Annual Review of Cell & Developmental Biology* **18**:637.
Gautel *et al.* (1999) Control of sarcomeric assembly: The flow of information on titin. *Reviews of Physiology Biochemistry and Pharmacology* **138**:97.
Wang *et al.* (2001) Single molecule measurements of titin elasticity. *Progress in Biophysics & Molecular Biology* **77**:1.