

User Guide

# Anti-Troponin T Antibody, Mouse Monoclonal

Clone JLT-12, Purified from Hybridoma Cell Culture

**T6277**

## Product Description

Monoclonal Anti-Troponin T (mouse IgG1 isotype) is derived from the hybridoma produced by the fusion of mouse myeloma cells and splenocytes from an immunized mouse. Purified Troponin T from rabbit skeletal muscle was used as the immunogen.

Monoclonal Anti-Troponin T is immunospecific for Troponin T in total rabbit skeletal muscle extract as determined by an immunoblot assay. The product reacts best with alcohol-fixed, paraffin-embedded or frozen sections of rabbit skeletal muscle; however, it shows broad species cross-reactivity in immunofluorescent labeling methods.

Troponin T is a microfilament protein closely associated with actin and its microfilament accessory proteins that appear in skeletal muscle myofibrils. Troponin T has a polypeptide molecular weight of approximately 38 kDa. Troponin T and tropomyosin from skeletal muscle run close to each other in 12.5% SDS polyacrylamide gels.

Monoclonal Anti-Troponin T may be used in immunofluorescent assays to study skeletal muscle organization. It can also be used to study the membrane anchorage sites of actin.

## Product Profile

The minimum working dilution of 1:200 was determined by immunoblotting using a total rabbit skeletal muscle extract.

In order to obtain best results, it is recommended that each individual user determine the optimum working dilution for their system by titration assay.

## Reagents

Supplied as a solution in 0.01 M PBS, pH 7.4, containing 15 mM sodium azide as a preservative.

## Precautions and Disclaimer

Due to the sodium azide content, a safety data sheet (SDS) for this product has been sent to the attention of the safety officer of your institution. Consult the SDS for information regarding hazards and safe handling practices.

## Storage

For continuous use, store at 2–8 °C for up to one month. For extended storage, the solution may be frozen in working aliquots. Repeated freezing and thawing is not recommended. If slight turbidity occurs upon prolonged storage, clarify by centrifugation before use.

---

## Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

### Technical Assistance

Visit the tech service page at [SigmaAldrich.com/techservice](https://SigmaAldrich.com/techservice).

### Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at [SigmaAldrich.com/terms](https://SigmaAldrich.com/terms).

### Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/offices](https://SigmaAldrich.com/offices).

The life science business of Merck operates as MilliporeSigma in the U.S. and Canada.

Merck and Sigma-Aldrich are trademarks of Merck or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

© 2002-2026 Merck and/or its affiliates. All Rights Reserved.  
23346842 Rev 02/26

