

TROP2 recombinant monoclonal antibody (046)

This antibody is covered by our [Worry-Free Guarantee](#).

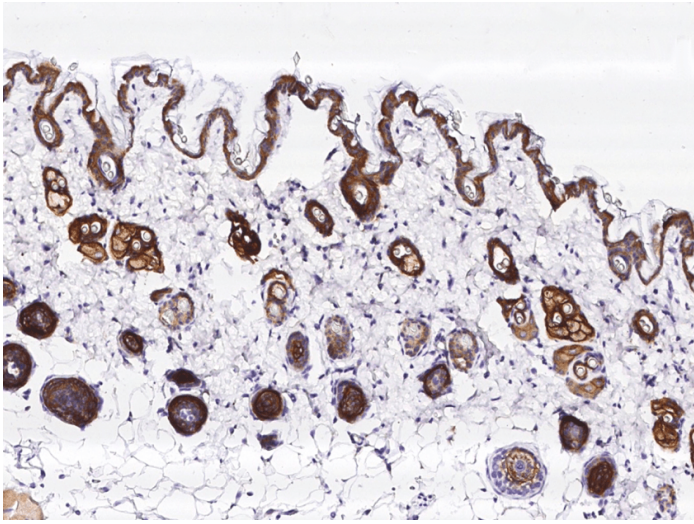
Ordering Information

[Order Online »](#)

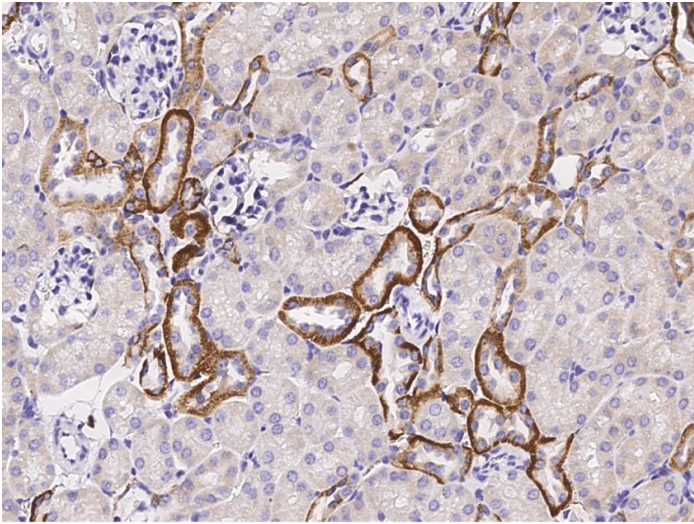
ENZ-ABS1025-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemical staining of mouse TACSTD2 in mouse skin with rabbit monoclonal antibody (1:20000, formalin-fixed paraffin embedded sections).



Immunohistochemical staining of mouse TACSTD2 in mouse kidney with rabbit monoclonal antibody (1:20000, formalin-fixed paraffin embedded sections).

Handling & Storage

Handling Protected from light. Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Tumor associated calcium signal transducer 2, GA733-1, M1S1, TACSTD2

Application IHC (PS)

Clone 046

Formulation Liquid. 0.2 µm filtered solution in PBS. Preservative free.

Host Rabbit

Immunogen Recombinant mouse TROP-2 / TACSTD2 protein.

Isotype IgG

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections) (1:10'000-1:20'000) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Mouse

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#)

.



Last modified: September 17, 2024

ENZO LIFE SCIENCES

Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office

ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands

& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France

Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany

Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland

Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com