TROP2 recombinant monoclonal antibody (046)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

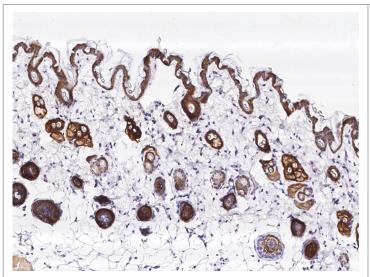
Order Online »

ENZ-ABS1025-0100

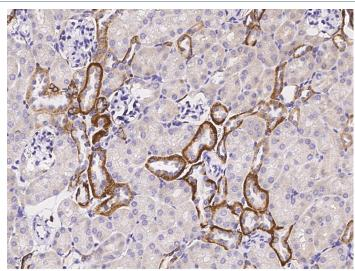
100µl

Manuals, SDS & CofA

View Online »



Immunochemical staining of mouse TACSTD2 in mouse skin with rabbit monoclonal antibody (1:20000, formalinfixed paraffin embedded sections).



Immunochemical 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.

lgG lgG

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

Phone: 800.942.0430
infousa@enzolifesciences.com

Phone: 441 61 926 8989
infoeu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com