S100B monoclonal antibody (60)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

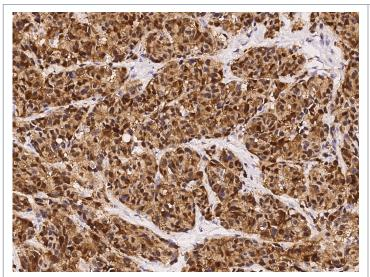
Order Online »

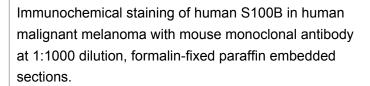
ENZ-ABS938-0100

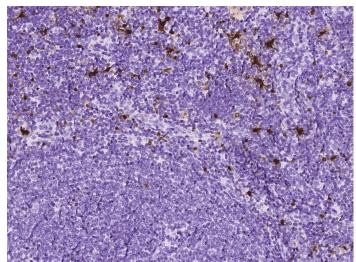
100µl

Manuals, SDS & CofA

View Online »







Immunochemical staining of human S100B in human lymphoma with mouse monoclonal antibody at 1:1000 dilution, 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

Application IHC (PS)

Clone 60

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

Host Mouse

Immunogen Recombinant human S100B protein.

lgG2a

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections) (1:200-

1:2'000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be

determined individually for each application.

Species Reactivity Human

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: September 17, 2024

