STAT6 monoclonal antibody (D-1)

STAT6 is a transcription factor activated downstream of IL-4 and IL-13 signaling. It is involved in Th2 immune responses and serves as a diagnostic marker for solitary fibrous tumors.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

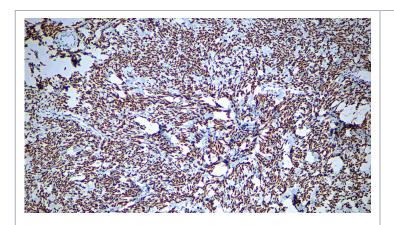
Order Online »

ENZ-ABS1084-0500

500µl

Manuals, SDS & CofA

View Online »



Formalin fixed paraffin embedded human solitary fibrous tumor stained with STAT6 antibody

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Signal transducer and activator of transcription 6, IL-4 Stat

Application IHC (PS)

Application Notes Pretreatment – EDTA Buffer pH 8.0. Incubation &

Temperature 30 min @ RT

Clone D-1

Formulation Tissue culture supernatant. Contains sodium azide as a

preservative.

Host Mouse

Immunogen Synthetic peptide corresponding to amino acids 799-823 at

the C-terminus.

Isotype IgG2b

Positive Control Solitary fibrous tumor

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections) (1:50-1:100)

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: June 6, 2025



info-

eu@enzolifesciences.com