Stat3 monoclonal antibody (7G8)

Stats (Signal transducers and activators of transcription) are a family of cytoplasmic latent transcription factors that are activated to regulate gene expression in response to a large number of extracellular signaling polypeptides including cytokines, interferons, and growth factors. After phosphorylation by JAK tyrosine kinases, Stats enter the nucleus to regulate transcription of many different genes.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

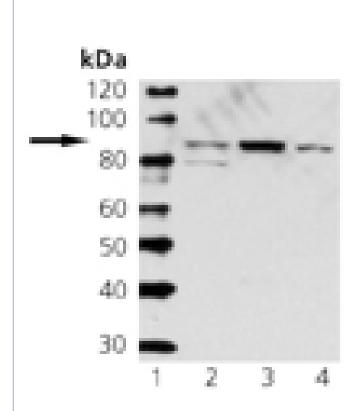
Order Online »

ADI-905-756-100

100µg

Manuals, SDS & CofA

View Online »



Western blot analysis of Stat3 Lane 1: MW, Lane 2: A431, Lane 3:HeLa, Lane 4: NIH-3T3.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Signal transducer and activator of transcription 3

Application WB

Application NotesDetects a band of ~88kDa by Western blot.

Clone 7G8

Formulation Liquid. In PBS, pH 7.2, containing BSA, 0.09% sodium

azide, and 50% glycerol.

Host Mouse

Immunogen Recombinant full length human Stat3.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Western Blot (1µg/ml, ECL)Suggested dilutions/conditions

may not be available for all applications. Optimal conditions

must be determined individually for each application.

Source Purified from mouse ascites.

Species Reactivity Human, Mouse, Rat

UniProt ID P40763

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

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com