RNA polymerase II monoclonal antibody (H14)

RPB1 is the catalytic and largest component of RNA polymerase II, which synthesizes mRNA precursors and many functional non-coding RNAs. It forms the polymerase active center together with RPB2, the second largest subunit. Polymerase II (Pol II) is the central component of the basal RNA polymerase II transcription machinery.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS134A-0100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name DNA-directed RNA polymerase II subunit RPB1

Application IF, IP, WB

Clone H14

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

Host Mouse

Immunogen Purified phosphorylated form of RNA polymerase II

extracted from a transformed cell line.

lsotype lgM

Purity Detail Purified by affinity chromatography

Species Reactivity Species independent

Specificity Recognizes RNA polymerase II phosphorylated at Ser⁵ in

a broad range of species (from yeast to human).

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

.

Last modified: May 29, 2024

