

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.

Isotype IgM

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](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com