## RNA polymerase II monoclonal antibody (8WG16)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

ENZ-ABS132A-0100

100µg

Manuals, SDS & CofA

**View Online** »

## **Handling & Storage**

**Handling** Avoid freeze/thaw cycles.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

**Application** IF, IP, WB

Clone 8WG16

**Formulation** Liquid. In PBS, pH 7.2.

**Host** Mouse

Immunogen RNA polymerase (RNA pol II) purified from wheat germ

extract.

lsotype lgG2a

Purity Detail Affinity purified

Species Reactivity Species independent

**Specificity** Recognizes RNA polymerase II in a broad range of

species (from yeast to human). Detects the C-terminal heptapeptide repeat present on the largest subunit of pol

II.

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

•

